

**UNITED STATES
SECURITIES AND EXCHANGE COMMISSION**
Washington, D.C. 20549

FORM 10-K

(Mark One)

ANNUAL REPORT PURSUANT TO SECTION 13 OR 15(d) OF THE SECURITIES EXCHANGE ACT OF 1934

For the fiscal year ended December 31, 2025

OR

TRANSITION REPORT PURSUANT TO SECTION 13 OR 15(d) OF THE SECURITIES EXCHANGE ACT OF 1934

For the transition period from _____ to _____

Commission file number: 001-38319

QUANTERIX CORPORATION

(Exact name of registrant as specified in its charter)

Delaware

(State or other jurisdiction of
incorporation or organization)

900 Middlesex Turnpike, Billerica, MA

(Address of principal executive offices)

20-8957988

(I.R.S. Employer Identification No.)

01821

(Zip Code)

Registrant's telephone number, including area code: **(617) 301-9400**

Securities registered pursuant to Section 12(b) of the Exchange Act:

Title of each class	Trading Symbol(s)	Name of each exchange on which registered
Common Stock, \$0.001 par value per share	QTRX	The Nasdaq Global Market

Securities registered pursuant to Section 12(g) of the Exchange Act: None

Indicate by check mark if the registrant is a well-known seasoned issuer, as defined in Rule 405 of the Securities Act. Yes No

Indicate by check mark if the registrant is not required to file reports pursuant to Section 13 or Section 15(d) of the Exchange Act. Yes No

Indicate by check mark whether the registrant (1) has filed all reports required to be filed by Section 13 or 15(d) of the Securities Exchange Act of 1934 during the preceding 12 months (or for such shorter period that the registrant was required to file such reports), and (2) has been subject to such filing requirements for the past 90 days. Yes No

Indicate by check mark whether the registrant has submitted electronically every Interactive Data File required to be submitted pursuant to Rule 405 of Regulation S-T during the preceding 12 months (or for such shorter period that the registrant was required to submit such files). Yes No

Indicate by check mark whether the registrant is a large accelerated filer, an accelerated filer, a non-accelerated filer, a smaller reporting company, or an emerging growth company. See the definitions of "large accelerated filer," "accelerated filer," "smaller reporting company," and "emerging growth company" in Rule 12b-2 of the Exchange Act.

Large Accelerated filer

Non-accelerated filer

Accelerated filer

Smaller reporting company

Emerging growth company

If an emerging growth company, indicate by check mark if the registrant has elected not to use the extended transition period for complying with any new or revised financial accounting standards provided pursuant to Section 13(a) of the Exchange Act.

Indicate by check mark whether the registrant has filed a report on and attestation to its management's assessment of the effectiveness of its internal control over financial reporting under Section 404(b) of the Sarbanes-Oxley Act (15 U.S.C. 7262(b)) by the registered public accounting firm that prepared or issued its audit report.

If securities are registered pursuant to Section 12(b) of the Act, indicate by check mark whether the financial statements of the registrant included in the filing reflect the correction of an error to previously issued financial statements.

Indicate by check mark whether any of those error corrections are restatements that required a recovery analysis of incentive-based compensation received by any of the registrant's executive officers during the relevant recovery period pursuant to §240.10D-1(b).

Indicate by check mark whether the registrant is a shell company (as defined in Rule 12b-2 of the Exchange Act). Yes No

As of the last business day of the registrant's most recently completed second fiscal quarter (June 30, 2025), the aggregate market value of the voting and non-voting common equity held by non-affiliates of the registrant, based on the last reported sales price for the registrant's common stock, par value \$0.001 per share, on the Nasdaq Global Market on such date, was approximately \$239 million.

As of February 24, 2026, the registrant had 46,939,975 shares of common stock outstanding.

DOCUMENTS INCORPORATED BY REFERENCE

Portions of the registrant's proxy statement for its 2026 Annual Stockholders' Meeting, which the registrant intends to file with the Securities and Exchange Commission pursuant to Regulation 14A within 120 days after the end of the registrant's fiscal year ended December 31, 2025, are incorporated by reference into Part III of this Annual Report on Form 10-K.

QUANTERIX CORPORATION

INDEX TO FORM 10-K

Note Regarding Forward-Looking Statements

	Page
<u>PART I</u>	
Item 1. Business	1
Item 1A. Risk Factors	22
Item 1B. Unresolved Staff Comments	46
Item 1C. Cybersecurity	46
Item 2. Properties	47
Item 3. Legal Proceedings	47
Item 4. Mine Safety Disclosures	47
<u>PART II</u>	
Item 5. Market for Registrant’s Common Equity, Related Stockholder Matters and Issuer Purchases of Equity Securities	48
Item 6. Reserved	49
Item 7. Management’s Discussion and Analysis of Financial Condition and Results of Operations	50
Item 7A. Quantitative and Qualitative Disclosures about Market Risk	65
Item 8. Financial Statements and Supplementary Data	65
Item 9. Changes in and Disagreements with Accountants on Accounting and Financial Disclosure	65
Item 9A. Controls and Procedures	65
Item 9B. Other Information	68
Item 9C. Disclosure Regarding Foreign Jurisdictions that Prevent Inspections	68
<u>PART III</u>	
Item 10. Directors, Executive Officers and Corporate Governance	69
Item 11. Executive Compensation	69
Item 12. Security Ownership of Certain Beneficial Owners and Management and Related Stockholder Matters	69
Item 13. Certain Relationships and Related Transactions, and Director Independence	69
Item 14. Principal Accountant Fees and Services	69
<u>PART IV</u>	
Item 15. Exhibits and Financial Statement Schedules	70
Item 16. Form 10-K Summary	74
Signatures	75
Consolidated Financial Statements	F-1

Unless the context otherwise requires, the terms “Quanterix,” the “Company,” “we,” “it,” “us,” and “our” in this Annual Report on Form 10 K refer to Quanterix Corporation and its consolidated subsidiaries.

Note Regarding Forward-Looking Statements

This Annual Report on Form 10-K contains forward-looking statements (within the meaning of the U.S. Private Securities Litigation Reform Act of 1995) that involve risks and uncertainties. All statements other than statements of historical facts contained in this Annual Report on Form 10-K are forward-looking statements. In some cases, forward-looking statements can be identified by words such as “anticipate,” “believe,” “contemplate,” “continue,” “could,” “estimate,” “expect,” “intend,” “may,” “plan,” “potential,” “predict,” “project,” “seek,” “should,” “target,” “will,” “would,” or the negative of these words or other comparable terminology. These forward-looking statements include, but are not limited to, statements about:

- the anticipated benefits and synergies from our acquisitions of Akoya Biosciences, Inc. ("Akoya") and Emission, Inc. ("Emission");
- the commercialization and adoption of our existing products and services and the success of our new product and service offerings, such as, for example, our LucentAD Complete multiplex test in the area of Alzheimer's disease diagnostics;
- our future financial performance and estimates regarding expenses, future revenues, capital requirements, and our needs for additional financing;
- the ability of our Simoa and Spatial Biology technology to improve existing diagnostics and to enable the development of new diagnostic tests and tools;
- the potential of our Simoa and Spatial Biology technology in the field of screening/diagnostic tests linked to therapeutic drugs and adoption by healthcare professionals;
- our ability to successfully penetrate the diagnostics market;
- our ability to retain and expand our customer base and achieve sufficient market acceptance of our products;
- the anticipated timing for launch of, and features of, our next-generation instruments to upgrade our existing platforms;
- the ability of our contract manufacturers and suppliers to reliably and consistently manufacture and supply our instruments;
- our ability to maintain effective internal control over financial reporting and disclosure controls and procedures; and
- the impact of changes in U.S. government policies, including impacts of tariffs and reductions in federal research funding.

These forward-looking statements are subject to a number of risks, uncertainties, and assumptions, including those further described in the section titled “Part I, Item 1A, Risk Factors” and elsewhere in this Annual Report on Form 10-K. Moreover, we operate in a very competitive and rapidly changing environment, and new risks emerge from time to time. It is not possible for us to predict all risks, nor can we assess the impact of all factors on our business or the extent to which any factor, or combination of factors, may cause actual results to differ materially from those contained in any forward-looking statements we may make. In light of these risks, uncertainties, and assumptions, the forward-looking events and circumstances discussed in this Annual Report on Form 10-K may not occur and actual results could differ materially and adversely from those anticipated or implied in the forward-looking statements.

Readers should not rely upon forward-looking statements as predictions of future events. Although we believe that the expectations reflected in any forward-looking statements are reasonable, we cannot guarantee that the future results, levels of activity, performance, or events and circumstances reflected in forward-looking statements will be achieved or occur. We undertake no obligation to update publicly any forward-looking statements for any reason after the date of this

Annual Report on Form 10-K to conform these statements to new information, actual results, or to changes in our expectations, except as required by law.

Readers should read this Annual Report on Form 10-K and any documents referenced herein that we have filed with the Securities and Exchange Commission (“SEC”) as exhibits to this Annual Report on Form 10-K with the understanding that our actual future results, levels of activity, performance, and events and circumstances may be materially different from what we expect.

This Annual Report on Form 10-K includes statistical, industry, and market data that we obtained from industry publications and research, surveys, and studies conducted by third parties. Industry publications and third-party research, surveys, and studies generally indicate that their information has been obtained from sources believed to be reliable, although they do not guarantee the accuracy or completeness of such information. Such data involves a number of assumptions and limitations and contains projections, and estimates of the future performance of the markets in which we operate and intend to operate that are subject to a high degree of uncertainty. We caution not to give undue weight to such projections, assumptions, and estimates.

Service Marks, Trademarks, and Trade Names

“Quanterix,” “Simoa,” “Simoa HD-X,” “Simoa ONE,” “SR-X,” “SP-X,” “HD-X,” “LucentAD,” “Lucent Diagnostics,” “Akoya,” “PhenoCycler,” “PhenoImager,” “PhenoCode,” and our logo are our trademarks. All other service marks, trademarks, and trade names appearing in this Annual Report on Form 10-K are the property of their respective owners. We do not intend our use or display of other companies’ trade names, trademarks, or service marks to imply a relationship with, or endorsement or sponsorship of us, by these other companies.

PART I

ITEM 1. BUSINESS

Overview

We are a life sciences company transforming healthcare innovation by accelerating biomarker breakthroughs from discovery to diagnostics using our ultra-sensitive translational research and spatial biology instruments, consumables, and services. We continue to invest in pushing a paradigm shift in healthcare from an emphasis on later-stage treatment to a focus on earlier detection, monitoring, prognosis, and, ultimately, prevention. Our combined platforms have achieved significant commercial adoption with an installed base of over 2,500 instruments and scientific validation with citations in more than 6,200 scientific publications in areas of high unmet medical need and research interest such as neurology, oncology, immunology, and inflammation.

Our proprietary digital “Simoa” detection technology enables customers to reliably detect protein biomarkers at ultra-low concentrations in blood, serum and other fluids that, in many cases, are undetectable using conventional, analog immunoassay technologies. Multi-plexing biomarker analysis in tissue samples with our “Spatial Biology” platforms enables scientists to understand the localized interactions occurring on the cellular level. We believe our combination of technologies will enable scientists to help drive diagnostic innovation in the evolving healthcare landscape with data across the tissue to fluid continuum. Currently, the ability of our Simoa platforms to detect proteins in the femtomolar range is enabling the development of novel therapies and diagnostics and has the potential to identify early-stage disease markers before symptoms appear.

We sell our proprietary instruments and related consumables worldwide to research laboratories, contract research organizations (“CROs”), academic institutions, and bio-pharmaceutical companies. In addition, we provide contract research services and clinical laboratory testing services, including four Laboratory Developed Tests (“LDTs”), using our proprietary technology and platforms through our Accelerator Laboratory, which is certified under the Clinical Laboratory Improvement Amendments of 1988 (“CLIA”), located in Bedford, Massachusetts (the “Accelerator Laboratory”).

Recent Developments

Recent Board of Directors and Management Transitions

In the second half of 2025, four new directors joined our board of directors and four previous members of our board of directors departed, including Dr. David Walt, our founding scientist. Additionally, in January 2026 Dr. Masoud Toloue left the company, and Everett Cunningham became our Chief Executive Officer and joined our board of directors. The selection process for these new board members and our CEO reflects a deliberate enhancement of technology, commercial, spatial, and diagnostics expertise across our leadership and board of directors.

Acquisition of Akoya Biosciences, Inc.

On July 8, 2025, we acquired Akoya Biosciences, Inc. (“Akoya”), a life sciences technology company based in Marlborough, Massachusetts delivering spatial biology solutions through the power of spatial phenotyping. Spatial phenotyping refers to a rapidly evolving technology that enables scientists to detect and map the distribution of cell types and biomarkers across whole tissue samples at single-cell resolution, enabling advancements in their understanding of disease progression and patient response to therapy. Akoya commercializes proprietary instrument platforms, reagents, software, and services that offer end-to-end solutions to perform tissue analysis and spatial phenotyping from discovery through translational and clinical research and diagnostics.

Acquisition of Emission, Inc.

On January 8, 2025, we acquired Emission, Inc. (“Emission”), a life sciences manufacturing company based in Georgetown, Texas. Emission produces large-scale, highly-uniform dye-encapsulating magnetic beads designed for low and mid-plex assays and a mid-plex platform that reads its proprietary beads. The transaction secured the use of Emission’s highly controlled beads in our next generation platforms and expansion into a new multi-plex segment targeting third-party original equipment manufacturer customers.

For further information about the Akoya and Emission acquisitions, refer to Note 3 - *Acquisitions* in the Notes to Consolidated Financial Statements for the year ended December 31, 2025 included in the section titled, "Part II, Item 8. Financial Statements and Supplemental Data" of this Annual Report on Form 10-K.

FDA 510(k) Submission for a Multi-Analyte Algorithmic Blood Test for Alzheimer's Disease Detection

On January 31, 2026, we submitted a 510(k) premarket notification to the U.S. Food and Drug Administration ("FDA") for our multi-analyte algorithmic blood test for Alzheimer's disease ("AD").

This submission represents a significant milestone in the Company's mission to provide superior, non-invasive, high-performance diagnostic tools to aid in the evaluation of patients with cognitive symptoms for possible Alzheimer's disease. Our multi-analyte test previously received Breakthrough Device Designation from the FDA, a program intended to accelerate the development and review of devices that provide for more effective treatment or diagnosis of life-threatening or irreversibly debilitating diseases. Our test, if cleared by the FDA, is intended to aid in identifying whether patients with cognitive symptoms are likely to have amyloid brain plaques—a hallmark of Alzheimer's disease—providing diagnostic clarity through a non-invasive blood test.

LucentAD Complete Medicare Pricing

In November 2025, the Centers for Medicare & Medicaid Services ("CMS") approved a reimbursement rate of \$897 for our LucentAD Complete multiplex test. This milestone provides a nationally recognized reference price, an important step for coverage decisions with payers, enables broader access, and supports efforts to bring this multiplex diagnostic solution to hospitals and laboratories across the country. In addition, based on available data, we believe LucentAD Complete has the potential to obtain a Local Coverage Determination, which would support more consistent reimbursement of claims.

Cooperation Agreement with Kent Lake Partners

As previously disclosed, on August 4, 2025, we entered into a Cooperation Agreement (the "Cooperation Agreement") with Kent Lake PR LLC, the general partner of Kent Lake Partners LP (together, "Kent Lake"), pursuant to which we agreed to retain an executive search firm to identify a potential candidate to be appointed to our board of directors (the "Board"). We also agreed to cooperate with Kent Lake to select an individual from the executive search firm's list of six candidates and to appoint such person as a Class I director by December 1, 2025. On November 20, 2025, the Board appointed two new directors, including one new Class I director.

Also pursuant to the Cooperation Agreement, we sought and obtained stockholder approval at our 2025 annual meeting of stockholders of an amendment to our Amended and Restated Certificate of Incorporation that declassified the Board, and we amended our bylaws to adopt a majority voting standard for uncontested director elections, with a plurality voting standard for contested director elections.

In the Cooperation Agreement, Kent Lake agreed to abide by certain voting commitments, customary standstill obligations and mutual non-disparagement, and no litigation provisions until the date that is thirty days prior to the nomination deadline under the bylaws for the nomination of director candidates for election to the Board at the 2027 annual meeting of stockholders, unless the Cooperation Agreement is earlier terminated in accordance with its terms.

Industry Background

Proteins are versatile macromolecules that serve critical functions in nearly all biological processes. Proteins are analytes that are highly relevant physiologically, providing real-time pictures of disease, and researchers and clinicians rely extensively on protein biomarkers for use in research and as clinical and diagnostic tools. However, normal physiological levels of many proteins are not detectable in easily accessible blood samples using conventional, analog immunoassay technologies, and many of these technologies can only detect proteins once they have reached levels that reflect more advanced disease or injury. For many other low abundance proteins, these technologies cannot detect proteins even at disease- or injury-elevated levels.

Enzyme-linked immunosorbent assay ("ELISA") technology has been the most widely used method of sensitive detection of proteins for more than 50 years. In simple terms, ELISA involves using a plate coated with an antibody that binds to the target molecule. If the target is present in the sample, it attaches to the plate via the antibody. Then, a second

molecule, often an enzyme-linked antibody, is added, binding to the target. The enzyme produces a detectable signal (like fluorescence or a color change) when a specific substrate is added, indicating the presence and quantity of the target molecule in the sample. Although ELISA is widely used in medical diagnostics and research, it has very significant limitations, including limited sensitivity and narrow dynamic range (i.e., the range of concentration of proteins being detected). Efforts to increase the sensitivity of conventional ELISA have had limited success due to procedural complexity and length.

Immunohistochemistry ("IHC") is the traditional technique used for detecting and localizing antibody protein markers in tissue samples. This microscopy application allows clinicians and scientists to visualize target antigens within tissue slides helping to understand their location and distribution to help diagnose cancer or tumor types. The application generates fluorescent signals through the use of antibodies linked to a fluorophore or secondary antibody carrying a detectable enzyme or dye. Our Spatial Biology platforms have advanced this concept to a highly multiplexed capability enabling a 2D/3D functional view of tissue and biomarkers. Detecting multiple biomarkers with this technology now enables significantly more detail to be extracted from the same tissue sample for better precision and understanding in the tissue microenvironment.

Spatial proteomics and fluid-based biomarker analysis both continue to evolve and expand their use case from core research markets to clinical applications. Quanterix believes that the combination of insights gained from the use of our instrumentation will drive innovation and application expansion supporting clinical decision making into the future.

Our Technologies

Simoa Technology

Our Simoa bead-based and planar array technologies are based on traditional ELISA technology but significantly advance conventional ELISA technology and are capable of substantially greater protein detection sensitivity. We believe that our Simoa platforms are among the most sensitive commercially available multiplex protein detection platforms.

Simoa Bead-Based Technology

Simoa bead-based digital immunoassays utilize the basic principles of conventional bead-based sandwich ELISA. However, unlike ELISA, which runs the enzyme-substrate reaction on all molecules in one well, Simoa bead-based reactions are run on individual molecules in tiny microwells, 40 trillionths of a milliliter, that are 2.5 billion times smaller than traditional ELISA wells. In traditional analog ELISA measurements, the detected signal increases in intensity as the concentration of a sample increases. In Simoa bead-based digital technology measurements, however, the detected signal relies on a binary signal/no signal readout, enabling single molecule detection, and analytical sensitivity in the femtomolar range compared with nanomolar and picomolar levels of detection in conventional ELISA.

HD-X



Our HD-X is our flagship instrument that empowers biomarker research and accelerates drug development. It is a fully automated immunoassay platform with multiplexing and custom assay capability designed to deliver significant productivity and operational efficiency improvements, as well as greater user flexibility. The HD-X is based on our bead-based technology, and assays run on the HD-X are fully automated.

SR-X



Our SR-X instrument is a compact benchtop instrument with a lower price point, more flexible assay preparation, and a wider range of applications. The SR-X utilizes the same Simoa bead-based technology and assay kits as the HD-X. In contrast to the fully automated workflow of the HD-X, the assay incubation and washing steps for the SR-X are performed outside of the instrument using conventional liquid handling methods. The offline sample prep provides additional flexibility to enable researchers to apply Simoa detection in an expanded range of applications.

Simoa Planar Array Technology

Simoa planar array immunoassays utilize the basic principles of conventional microplate-based sandwich ELISA. However, unlike ELISA, which runs the enzyme-substrate reaction on all molecules coating the entire bottom surface in one well, Simoa planar array reactions are run on spatially segregated micro-spots within the bottom of microtiter plate wells that concentrate the signal to a surface area 1,000 times smaller than a traditional ELISA. The small spot size and spatial segregation of each spot enables multiplexing up to 12 different assays within a single sample well.

SP-X



Our SP-X uses Simoa planar array technology for multiplex chemiluminescent immunoassay measurement with sensitivity similar to that found in our Simoa bead-based platforms.

Simoa Assays and Consumables

Recurring revenue is derived through the sale of consumables used to run assays on our instruments and from our growing menu of Simoa digital biomarker assays. The current menu of approximately 65 analyte-specific single-plex and multi-plex assay kits for our bead-based instruments includes assays for biomarkers in the areas of neurology, infectious disease, immunology and oncology for both human and mouse samples. The current menu of assay kits for the planar array instrument includes approximately 45 biomarkers ranging from 1-10 analytes per assay in the areas of immunology, and oncology research.

In addition to the assays we have developed, the Simoa platforms allow ease and flexibility in assay design, enabling our customers to develop their own proprietary in-house assays, which we refer to as "homebrew" assays, using our homebrew assay kits. These kits include all components required for customers to run tests using their own antibodies. Our consumables portfolio for our bead-based platform also includes our proprietary Simoa disks that are unique to our bead-based platform, as well as cuvettes and disposable tips.

Spatial Technology

Our Spatial Biology solutions measure protein expression and cellular phenotypes within intact tissue while preserving spatial context and tissue architecture. By enabling the simultaneous detection of multiple biomarkers at single-cell resolution across whole tissue sections, our platforms provide quantitative insights into cellular composition, organization, and interactions within complex biological systems.

Conventional genomic and proteomic techniques, including next-generation sequencing ("NGS"), single-cell analysis, flow cytometry, and mass spectrometry, generate valuable molecular data but generally require tissue dissociation or destruction. As a result, these methods do not preserve spatial relationships between cells or localized biomarker expression within the native tissue environment. Traditional histological and non-destructive tissue analysis methods preserve tissue structure but are typically limited to measuring a small number of biomarkers per sample and often rely on qualitative interpretation by trained pathologists.

Our Spatial Biology platforms are designed to address these limitations by enabling end-to-end spatial phenotyping, allowing researchers to quantitatively analyze large numbers of protein biomarkers and cell types across intact tissue sections at single-cell resolution. The resulting data generate high-dimensional, spatially resolved maps that characterize tissue architecture, cellular heterogeneity, and biomarker co-localization, supporting the study of tissue dynamics and cell-cell interactions.

Our Spatial Biology technology platform integrates proprietary imaging hardware, scalable multiplexing enabled by barcoded antibody chemistry and classic Opal™ chemistry, and curated biomarker panels designed for standardized implementation. Together, these components support reproducible, whole-slide imaging and the generation of structured, computationally analyzable datasets.

We believe our platform capabilities support applications across the research continuum, from exploratory and discovery-stage research through translational and clinical research settings. These capabilities are intended to enable customers to advance the understanding of human biology, disease mechanisms, and therapeutic response, and to support the progression of spatial biology insights into downstream research and clinical workflows.

Spatial Biology Instruments

PhenoCycler & PhenoImager Fusion



Our PhenoCycler and PhenoImager Fusion instrument modules operate in a coordinated manner and provide the customer with a complete system for performing high-quality, high-plex tissue biomarker measurements.

Our PhenoCycler employs a barcoded antibody-based chemistry designed to enable highly multiplexed protein detection through cyclic signal readout rather than sequential antibody staining. In this workflow, all antibodies in a panel are applied to the tissue simultaneously in a single staining step. Each antibody is conjugated to a unique oligonucleotide barcode that serves as a molecular identifier for its target protein.

Following antibody binding, signal detection is achieved through iterative cycles of oligonucleotide hybridization and dehybridization. In each cycle, a defined set of fluorescently labeled reporter oligonucleotides—each complementary (antisense) to a subset of the antibody barcodes—are introduced together with a nuclear counterstain. Each cycle enables detection of multiple protein targets through reporter hybridization, while the nuclear stain provides consistent spatial reference across cycles.

After reporter hybridization, the tissue is imaged using the PhenoImager Fusion to capture spatially resolved signal from the bound reporters. The reporter oligonucleotides are then removed through controlled dehybridization, enabling subsequent cycles of reporter hybridization and imaging on the same tissue section. This process allows repeated signal acquisition without repeated antibody staining or tissue stripping.

This cyclic hybridization, imaging, and dehybridization approach enables multiplexed detection of protein biomarkers while preserving tissue integrity and spatial context. The cycling steps are automated through the PhenoCycler fluidics system and are designed to operate in conjunction with the PhenoImager Fusion reporter imaging platform, which performs whole-slide imaging following each reporter hybridization cycle.

PhenoImager HT



Our PhenoImager HT is a whole-slide imaging platform designed to support multiplexed tissue imaging using single-round staining and imaging workflows. The system enables one-shot detection of multiple protein biomarkers within a single imaging acquisition, typically supporting low- to mid-plex immunofluorescence and brightfield assays while preserving tissue morphology and spatial context.

PhenoImager HT is designed to image formalin-fixed paraffin-embedded ("FFPE") and fresh frozen tissue sections and tissue microarrays. The platform supports both fluorescence and chromogenic detection methods and is compatible with established IHC and immunofluorescence workflows commonly used in translational research and clinical laboratory environments.

Unlike cyclic imaging approaches, PhenoImager HT performs multiplexed imaging in a single imaging step, enabling streamlined workflows and reduced assay complexity. This one-shot imaging approach is intended to support applications where moderate levels of multiplexing are sufficient and where workflow simplicity, reproducibility, and compatibility with existing laboratory practices are important considerations.

Image data generated by PhenoImager HT are processed into our proprietary QTIFF file format, as well as additional standard image formats, to support downstream analysis using our software solutions and third-party image analysis platforms. The system is commonly used in translational research, biomarker validation, and LDT workflows, and has historically supported applications bridging discovery research and clinical investigation.

Spatial Biology Assays

Our Spatial Biology assays are based primarily on proprietary barcoding chemistry designed to support multiplexed protein detection in spatial biology applications. This chemistry utilizes barcoded antibodies in combination with cyclic signal detection workflows, enabling multiplexed analysis of protein expression while preserving tissue architecture and spatial context. These assays are designed for use across our Spatial Biology platforms, including PhenoCycler, PhenoImager Fusion, and PhenoImager HT, depending on the assay format and level of multiplexing required.

We offer a portfolio of ready-to-use, preconfigured antibody panels under our PhenoCode brand. For the PhenoCycler platform, these include curated discovery panels designed for high-plex spatial phenotyping applications, such as the PhenoCode IO60 panel, Mouse IO panel, Human Neuro panel, and Mouse Neuro panel. These panels are designed to support standardized implementation of multiplexed spatial assays and are intended to reduce assay development complexity for customers.

For our PhenoImager platforms, we offer PhenoCode Signature Panels, which are designed for single-round, low- to mid-plex imaging workflows. These panels are intended to support translational research, biomarker validation, and clinical research workflows where streamlined assay execution and reproducibility are important considerations.

In addition to ready-to-use panels, we sell a broad range of reagents on an à la carte basis, including barcoded antibodies, reporters, buffers, and ancillary reagents. These offerings allow customers to design and optimize custom panels and workflows tailored to specific research objectives and experimental requirements. This modular reagent approach is intended to support flexibility across discovery, translational, and applied research use cases.

We also support Opal ‘classic’ chemistry, an immunofluorescence method based on unconjugated primary antibodies, secondary antibodies conjugated to polymer HRP, and a suite of TSA-conjugated fluorophores. Opal chemistry protocols for multiplexed measurements can be automated on industry-standard autostainers. The signal amplification provides enhanced sensitivity. Opal chemistry is compatible with standalone PhenoImager Fusion and PhenoImager HT systems. This chemistry is commonly used in established immunohistochemistry and immunofluorescence workflows and provides continuity for customers with existing assay investments.

Services

Through our CLIA-certified Accelerator Laboratory, we provide research customers contract research services and we provide medical providers clinical laboratory testing services. The Accelerator Laboratory supports multiple projects and services, including:

- *Research and Clinical Trial Testing:* Utilizing either our commercially available reagent kits, or customer specific custom assays, our Accelerator Laboratory provides customers access to our Simoa and Spatial Biology platforms by running studies ranging from thousands of specimens to small *de novo* experiments with just a few samples. Partnering with clients to address critical needs on sample volume, custom assay development, and unique study designs, we support a broad range of programs from large scale clinical trial data generation to research focused cohorts supporting discovery, validation, and clinical testing of predictive biomarkers in translational research and enabling patient stratification and selection in clinical studies.
- *Lucent Diagnostics:* In July 2023, we launched Lucent Diagnostics, a suite of clinical laboratory testing services for neurology healthcare providers, initially focused on Alzheimer’s disease. To support a streamlined provider workflow, Lucent Diagnostics includes a web portal where healthcare providers can order sample collection materials, order a test, track the status of a test, and retrieve the test report. The process for testing involves a healthcare provider drawing a blood sample from an individual and shipping the sample to our Accelerator Laboratory to analyze the sample and communicate the result back to the provider. Healthcare providers can currently order the following tests through Lucent Diagnostics:
 1. *LucentAD p-Tau 181:* an LDT to quantitatively measure p-Tau 181 in plasma as an aid in diagnosis of Alzheimer’s disease.
 2. *Simoa NfL LDT:* an LDT to quantitatively measure neurofilament light chain ("NfL") in serum as an aid in the evaluation of individuals for possible neurodegenerative conditions or other causes of neuronal or central nervous system damage.
 3. *LucentAD p-Tau 217:* an LDT to quantitatively measure p-Tau 217 in plasma to aid in diagnosis of Alzheimer’s disease.
 4. *LucentAD Complete:* a multi-marker (p-Tau 217, amyloid β 42 ("A β 42"), amyloid β 40 ("A β 40"), NfL, and glial fibrillary acidic protein ("GFAP")) algorithmic LDT for high accuracy plasma detection of amyloid pathology to aid in diagnostic evaluation of patients with cognitive issues who may have Alzheimer’s disease. By using a proprietary algorithm to score five Alzheimer’s disease-related biomarkers, LucentAD Complete provides significantly better amyloid classification compared to single-marker tests alone and has lowered the intermediate zone of a single-marker test from approximately 30% to 10%, while maintaining high sensitivity, specificity, and accuracy.

Extended Warranty and Service Contracts

We also generate revenues through extended-warranty and service contracts for our installed base of instruments.

Our Strategy

Our commercial strategy is based on three core growth vectors: (i) Grow menu; (ii) Expand into adjacencies; and (iii) Translate into diagnostics. Our plans to achieve this GET strategy include:

- **Growing through rapid internal menu expansion.** In 2024 and 2025, we launched 20 and 11 new Simoa assays, respectively that were developed using improved protocols, manufacturing efficiencies, and reagent improvements to provide more consistent results and improved lot-to-lot consistency. In 2025, we also launched two Spatial Biology panels. We expect that the expansion of our assay menu to increase average consumable pull-through as additional assays become available for use on our existing installed base.
- **Continuing to grow our leadership position in the measurement of neuro-based biomarkers.** We intend to leverage the growing importance of neurological biomarkers to advance the development of therapeutics and diagnostics for neurodegenerative conditions, including Alzheimer's disease. The importance of neurological biomarkers, such as NfL, p-Tau 181, p-Tau 217 and GFAP, has increased significantly in recent years, and our ultra-sensitive Simoa platforms have allowed research of neurological disorders, previously limited primarily to cerebrospinal fluid ("CSF"), to expand significantly.
- **Expanding further into indications beyond neurology.** The ability of our Simoa technology to detect and quantify low abundance proteins with exquisite sensitivity is a distinct advantage in detecting disease earlier in other indications, in particular oncology and immunology. There are over 1020 scientific publications that reference Simoa technologies in non-neurology indications. We expect to expand our Simoa product offerings into adjacencies through internal research and development investment. The acquisition of Akoya establishes us in the spatial biology market with applications across research and clinical trial settings. The acquisition diversifies our revenues by disease type, shifting from over 90% neurology to a diversified mix of approximately 60% neurology and 40% oncology. Furthermore, we believe researchers will be able to combine the capabilities of our multiplexed biomarker detection in tissue with our ultra-sensitive biomarker detection in blood to achieve new discoveries about disease onset and progression in oncology and immunology.
- **Expanding our presence in diagnostics.** Our ultra-sensitive Simoa platforms have enabled the development of a new category of less-invasive diagnostic tests that could replace current invasive, expensive, and inconvenient diagnostic methods, including spinal tap and diagnostic imaging. We currently offer four neurological LDTs through our Accelerator Laboratory.
- **Helping to build the global infrastructure for Alzheimer's disease testing and diagnosis.** There are over 55 million people living with Alzheimer's disease, a number that is expected to double by 2050. We believe our ultrasensitive technology can enable earlier detection. We are helping to build the global infrastructure for Alzheimer's testing in two ways. First, we are enabling partners by providing Simoa technology to a number of the top prescribing hospital networks and reference labs globally. We entered into 13 of these partnerships in 2025 bringing the total number of partners at the end of 2025 to 25 and intend to continue to grow this network in 2026. Second, through Lucent Diagnostics, we are offering best-in-class diagnostic testing through our LucentAD Complete test, which was launched in November 2024.

Our Key Focus Areas

We have focused the application of our technologies on areas of high growth and high unmet need and where existing platforms have significant shortcomings that our technology addresses, including neurology, oncology, and immunology.

Neurology

The ability of Simoa technology to detect neurological biomarkers in blood at ultra-low levels, combined with our multiplexing capability, has significantly advanced neurology research, drug development, and diagnostics test development. Prior to the launch of our p-Tau 181 LDT for clinical use in July 2022, the brain was the only organ in the body for which there was not a blood-based diagnostic test. The challenge with developing blood-based tests for the brain is that the blood-brain barrier, which is formed by endothelial cells lining the cerebral microvasculature, is very tight and severely restricts the movement of proteins and other substances between these endothelial cells and into blood circulation. Accordingly, diagnosis of brain disease and injury has traditionally required either brain imaging or a spinal tap to collect

CSF. The sensitivity of Simoa technology has enabled researchers to discover that extremely small amounts of critical neural biomarkers diffuse through the blood brain barrier and are released into the blood during injury and in connection with many neurodegenerative brain diseases. However, the concentrations of many of these neural biomarkers in the blood can be so low that they are difficult to detect by conventional, analog immunoassay technologies. Furthermore, neurological pathophysiology is complex, and it has become clear that no single biomarker is sufficient to serve as comprehensive indicator of these processes. For this reason, the capability of Simoa to multiplex blood-based tests for multiple neural biomarkers into a single test has emerged as useful to facilitate biomarker profiling of neurological disorders. This capability has materially contributed to rapid research progress, notably in the Alzheimer's disease landscape.

Our Spatial Biology platforms enable high-plex, single-cell protein analysis directly within intact brain tissue, making them highly relevant to neurology research and neurodegenerative disease markets. Neurological disorders such as Alzheimer's disease, Parkinson's disease, multiple sclerosis, and amyotrophic lateral sclerosis ("ALS") are driven by complex cellular interactions—including neurons, glia, immune cells, and vascular components—that cannot be fully understood through bulk or dissociated assays. Our Spatial Biology technologies allow researchers to localize and quantify dozens of protein biomarkers simultaneously while preserving anatomical context, supporting the identification of disease-associated cell states, pathogenic microenvironments, and therapeutic targets. This capability is increasingly important as pharmaceutical and academic groups prioritize biomarker-driven drug development, patient stratification, and translational research in neurology, positioning spatial biology as a foundational tool in a large, growing, and underserved research and clinical development market.

Developments on the Alzheimer's Disease Landscape

According to the Alzheimer's Association, there are over 7 million individuals living with Alzheimer's disease and other dementias in the United States, and that figure is expected to double by 2060. The FDA's approval of Leqembi and Kisunla as disease-modifying treatments for Alzheimer's disease has underscored an urgent need for non-invasive, widely available blood tests to facilitate diagnosis in the early stages of the disease to identify patients for treatment when therapeutic intervention is most likely to provide clinical benefit. Established biomarker-based approaches to diagnostic workup for Alzheimer's include positron emission tomography ("PET") imaging and CSF biomarkers for amyloid and phosphorylated tau, both of which are invasive, expensive, and may not be widely available. Recently two blood tests for assessing the risk of brain amyloid pathology associated with Alzheimer's disease in individuals with early signs of cognitive impairment were cleared by the FDA. This is an important milestone demonstrating support by regulatory authorities for the use of high quality blood tests as part of Alzheimer's diagnostic workflows. Assessing the effectiveness of an anti-amyloid Alzheimer's therapy by testing for amyloid plaque reduction over the course of treatment with serial testing represents another potential opportunity for blood biomarker tests. The availability of a simple blood test for amyloid positivity has the potential to alleviate bottleneck and cost burdens and facilitate the accessibility of Alzheimer's therapeutic intervention for patients. We believe our Simoa technology is well-suited to meet these critical needs.

Simoa Multi-Marker Algorithmic Test

Leveraging both the multiplexing and high sensitivity capability of Simoa, we have developed a multi-marker algorithmic test that combines five biomarkers (p-Tau 217, A β 42, A β 40, NfL, and GFAP) and uses a proprietary algorithm for high accuracy plasma detection of amyloid pathology. Researchers have extensively studied these biomarkers to advance the understanding of how blood-based biomarkers reflect different aspects of Alzheimer's pathology, to both further the basic knowledge of the disease pathophysiology, and to explore the potential of these biomarkers to predict disease state, severity, and progression. This work has resulted in a large body of published evidence that has advanced the understanding of the disease and facilitated drug development through the use of these biomarkers as exploratory indicators of drug target engagement. From a diagnostic test standpoint, a central concept has emerged: since these biomarkers reflect different aspects of Alzheimer's pathology, combining their signals through logistic regression can enhance overall test accuracy. This information, together with clinical evaluations and cognitive testing, can help determine whether cognitive symptoms are the result of Alzheimer's disease or another type of dementia.

We have partnered with the Alzheimer's Drug Discovery Foundation to develop and clinically implement a multi-marker blood test to aid Alzheimer's diagnosis and differential diagnosis of patients presenting with cognitive symptoms of uncertain origin. Validation data for the clinical performance of the test for amyloid detection was presented at the Clinical Trials on Alzheimer's Disease annual meeting late 2024. The data across more than 1,000 patients showed that the test delivered best-in-class performance for amyloid detection accuracy as well as reduced uncertainty in borderline cases. We launched this multi-marker algorithmic test as an LDT in November 2024 under the brand name LucentAD Complete.

Additionally, in November 2025 CMS approved a reimbursement rate of \$897 for this test. A version of this multi-analyte algorithmic blood test received Breakthrough Device designation by the FDA, and we submitted a 510(k) premarket notification to the FDA earlier this year.

Other Neurological Conditions

Our ultra-sensitive Simoa technology has also been instrumental in advancing research into other neurological conditions, such as multiple sclerosis and traumatic brain injury ("TBI"). Evidence of the potential clinical utility of NfL as a biomarker in multiple sclerosis has progressed rapidly, and Simoa's role in that progression has been foundational. In April 2022, the FDA granted our Simoa NfL serum test Breakthrough Device designation as a prognostic aid in assessing the risk of disease activity in patients diagnosed with relapsing-remitting multiple sclerosis ("RRMS"). The test has shown promise to be used in conjunction with clinical, imaging and laboratory findings as an aid in identifying RRMS patients who are at lower or higher risk for relapse within four years. We believe this prognostic information could be clinically useful in tailoring the therapeutic approach to more effectively treat the disease.

Current methods of TBI diagnosis involve CT scans that fail to diagnose approximately 90% of mild TBI. Simoa technology has demonstrated the sensitivity to identify relevant neurological biomarkers, such as NfL, tau, GFAP, and ubiquitin C-terminal hydrolase L1 ("UCH-L1"), to more adequately aid in diagnosis of TBIs and overall brain health. Researchers in neurology have used Simoa technology to study biomarkers in the blood of athletes after concussion in many high-impact sports. Simoa can measure critical neural biomarkers in blood that correlate to repeated head trauma from both concussions and subconcussive events with poor patient outcomes, including the potential development of Chronic Traumatic Encephalopathy, which currently can only be diagnosed after death via a brain autopsy.

Additional assays for blood-based biomarkers with potential utility in Alzheimer's research and diagnostics have either been launched or are in development. These include brain-derived Tau, p-Tau 231, p-Tau 212, sTREM2, and PSD95.

Oncology and Immunology

Our Spatial Biology platforms are highly relevant to oncology research due to their ability to resolve the tumor microenvironment at single-cell resolution while preserving tissue architecture. Cancer progression, immune evasion, and therapeutic response are driven by spatially organized interactions among tumor cells, immune infiltrates, stromal cells, and vasculature—features that are not captured by bulk or non-spatial assays. Our technologies enable simultaneous measurement of dozens of protein biomarkers within intact tumor tissue, supporting biomarker discovery, mechanism-of-action studies, and identification of predictive and pharmacodynamic markers for immuno-oncology and targeted therapies. As oncology drug development increasingly relies on biomarker-guided patient selection and translational insights linking tissue biology to clinical outcomes, spatial biology has become a critical enabling technology, underpinning demand from pharmaceutical companies, translational research centers, and clinical research laboratories in one of the largest and fastest-growing life sciences markets.

Cancer immunotherapy is a promising new area that is significantly affecting cancer remission rates. One challenge of immunotherapy approaches is that the elicited immune responses are not always predictable and can vary from person to person and protocol to protocol. There exists a significant need to develop biomarker tools to monitor these drugs and their effects. Circulating protein biomarkers (serum and plasma) have the potential to be used in the field of health-oncology to stratify patients, predict response, predict recurrence, reveal mechanism of action and monitor for adverse effects. One technical challenge facing the health-oncology drug development process has been the availability of immunoassays with sufficient sensitivity to measure immunomodulatory biomarkers directly in serum and plasma. We have developed a number of tumor biomarker and immune modulation assays (cytokines and chemokines) that can be used to monitor tumor proliferation and host immune response. In particular key immune regulatory cells (T-regs, dendritic cells, macrophages) secrete very low amounts of the protein Interferon gamma ("IFN-gamma") and these levels cannot be reliably measured in serum and plasma using conventional, immunoassay technology, however they can be tracked with our Simoa IFN-gamma assay. Additionally, we have developed an ultra-sensitive assay for ubiquitin C-terminal hydrolase L1 ("IL-6"), which is one of the cytokines commonly measured for monitoring cytokine release syndrome as an adverse effect in immunotherapies. Several studies have shown that our ultrasensitive assays can be valuable tools for monitoring health-oncology drugs and protocols.

Inflammation

Inflammation underlies the response of the body to injury in a variety of diseases. Simoa assays can measure inflammatory and anti-inflammatory molecules in serum and plasma with unprecedented sensitivity. This has the potential to enable new discoveries into the role of inflammation in the biology of health and disease. Our Simoa technology measures low levels of inflammatory proteins, including cytokines and chemokines, that characterize a range of inflammatory diseases, including Crohn's disease, asthma, rheumatoid arthritis, and neuro-inflammation. We believe the sensitivity of Simoa technology can provide a clearer picture of the underlying state of the immune response and disease progression.

Combining the ultra-sensitive capabilities of our Simoa technology with our Spatial Biology platform is highly relevant to inflammation research because inflammatory diseases are driven by localized, cell-to-cell signaling events that vary by tissue compartment and disease stage. Conditions such as autoimmune disorders, fibrosis, cardiovascular disease, and inflammatory bowel disease involve complex interactions among immune cells, parenchymal cells, and stromal elements that cannot be adequately characterized using bulk or non-spatial methods. Our Spatial technologies enable high-plex, single-cell protein analysis within intact tissue, allowing researchers to map inflammatory pathways, immune cell activation states, and spatial biomarkers associated with disease progression and therapeutic response.

Research and Development

We continue to seek to improve our technologies to enable more sensitive detection and measurement of biological molecules. This effort includes examining new assay formats to increase their performance. We intend to expand our assay menu to extend the scope of applications for our platforms to biomarkers of significant interest to the scientific community. Further, we are developing new instrumentation to update our current Simoa platforms, with a design that aims to ultimately have higher sensitivity and multiplexing. For example, in late 2025, we launched an early access program to our next generation instrument "Simoa ONE" to give key evaluators hands-on experience with the technology. We continue to gather feedback on this instrument to innovate its design and determine which features are most important.

Sales and Marketing

We distribute our instruments and consumables via direct field sales and support organizations located in North America and Europe and through a combination of our own sales force and third-party distributors in additional countries, including Australia, Brazil, China, Czech Republic, Cypress, India, Hong Kong, Israel, Japan, New Zealand, Qatar, Saudi Arabia, Singapore, South Africa, South Korea, Taiwan, and UAE. In addition, we sell NFL antibodies and NFL ELISA kits produced by our Swedish subsidiary, UmanDiagnostics AB ("Uman"), directly and in conjunction with a distributor worldwide.

As of December 31, 2025, we had 151 employees in sales, sales support, and marketing, including technical field application scientists and field service personnel. This staff is primarily located in North America and Europe. We expect we will expand our sales, sales support, and marketing efforts in North America as it relates to the commercialization of our LucentAD Complete test. Additionally, we expect to continue to develop a comprehensive distribution and support network in China and other APAC region countries where significant new opportunities exist.

Manufacturing and Supply

We outsource the manufacturing of our Simoa bead-based instruments and Spatial Biology instruments to third-party manufacturers, and we manufacture our planar array instrument and all assay kits in our own facilities.

Instruments

Our HD-X instrument is supplied by STRATEC, based in Birkenfeld, Germany, and is manufactured and shipped from its Birkenfeld and Beringen, Switzerland facilities. Our SR-X instrument is supplied by Paramit, based in Morgan Hill, California, and is shipped to our global customers by Paramit. Our Spatial Biology instruments are manufactured by Columbia Tech, based in Westborough, Massachusetts. See the section titled "*Key Agreements*" for a description of our agreements with these manufacturers. Installation of, and training on, our instruments is provided by our employees where we conduct direct sales, and by distributors where sales are conducted through distributors.

We believe this manufacturing strategy is efficient and conserves capital. However, in the event it becomes necessary to utilize a different contract manufacturer for the HD-X, SR-X, or Spatial Biology instruments, we could experience additional costs, delays and difficulties in doing so, and our business could be harmed.

Our SP-X instrument is manufactured, tested, shipped and supported by us from our Billerica, Massachusetts facility. All internal components are sourced domestically except one significant component that is sourced in Germany. These components are sourced from a limited number of suppliers, including certain single-source suppliers. Although we believe that alternatives would be available, it would take time to identify and validate replacement components, which could negatively affect our ability to supply SP-X instruments on a timely basis.

Consumables

We assemble our assay kits for our Simoa bead-based platforms and Spatial Biology platforms in our Billerica, Massachusetts facility. Our bead-based assays include all components required to run an enzyme-based immunoassay, such as beads, capture and detector reagents, enzyme reagents and enzyme substrate. Reagents for our assays are sourced from a limited number of suppliers, including certain single-source suppliers. Although we believe that alternatives would be available, it would take time to identify and validate replacement reagents for our assay kits, which could negatively affect our ability to supply assay kits on a timely basis.

Simoa disks for our bead-based platforms are supplied through a single source supplier pursuant to a long-term supply agreement with STRATEC Consumables GmbH, a subsidiary of STRATEC. We believe that this agreement provides for a sufficient notification period to allow for supply continuity and the identification and tech transfer to a new supplier in the event either party wishes to terminate the relationship. Our cuvettes for our bead-based platforms are single sourced through STRATEC, and the disposable tips used in our bead-based platforms are commercially available.

We assemble our 96-well sample plate kits for our planar array platform in our Billerica, Massachusetts facility. Reagents for our planar array assays include all components required to run an enzyme-based chemiluminescent immunoassay, such as capture antibody printed plates and detector reagents, enzyme reagents and enzyme substrate. These reagents are sourced from a limited number of suppliers, including certain single-source suppliers. Although we believe that alternatives would be available, it would take time to identify and validate replacement reagents for our assay kits, which could negatively affect our ability to supply assay kits on a timely basis. Because our planar array assays have a shelf life of 12 months, we believe we are able to mitigate this risk through inventory control.

NfL antibodies and NfL ELISA Kits

The storage of Uman's proprietary NfL antibody producing hybridomas, as well as the cultivation and purification of the antibodies, is outsourced to a contract manufacturer, and bulk material of purified antibodies is delivered to Uman's site in Umeå, Sweden. Functional testing and verification of concentration are performed at Uman before the material is approved for use in production activities. The antibodies currently have a shelf life of 18 months and can be aliquoted and sold as single reagents or used for the production of Uman's NfL ELISA kits.

All components in Uman's NfL ELISA kits are manufactured in-house at Uman from starting materials sourced from suppliers that have been evaluated and approved. Uman has entered into supply agreements with critical suppliers. The kit components include buffers (sample diluents and wash solutions), ELISA 96-well plates coated with a capture antibody, detector antibodies, streptavidine conjugates, substrates (TMB) and stop reagents. The final ELISA kit products are subject to quality control procedures, which include testing of human CSF or human serum quality control samples to assure a high batch consistency. After testing and batch record review, the material is released to market. The current shelf-life of the kits is 18 months (NfL ELISA (CSF)) or 13 months (NfL Serum ELISA).

Key Agreements

Development Agreement and Supply Agreement with STRATEC

In August 2011, as amended in November 2016 and July 2025, we entered into a Strategic Development Services and Equity Participation Agreement with STRATEC, pursuant to which STRATEC undertook the development of the Simoa HD instrument. In September 2011, as amended in October 2013, we also entered into a Supply and Manufacturing Agreement with STRATEC (the "STRATEC Supply Agreement"), pursuant to which STRATEC agreed to supply HD instruments to us, and we agreed to procure those instruments exclusively from STRATEC, subject to STRATEC's ability

to supply the instruments. We are responsible for obtaining any regulatory approval necessary to sell the instruments. The instrument price stipulated in the STRATEC Supply Agreement was established based on certain specified assumptions and is subject to certain adjustments.

The STRATEC Supply Agreement is terminable by either party on 12 months' notice to the other party. The STRATEC Supply Agreement may also be terminated on the insolvency of a party or the uncured material breach of a party, or, by us, on a change of control of our company (subject to certain obligations to compensate STRATEC on such termination). On termination by us for STRATEC's insolvency or uncured material breach or termination by STRATEC for convenience, we are granted a nonexclusive royalty free license of STRATEC intellectual property to manufacture the instruments.

Paramit Manufacturing Services Agreement

In November 2016, we entered into a Manufacturing Services Agreement (the "Paramit Agreement") with Paramit to produce and test our SR-X instrument on an as-ordered basis. We also engaged Paramit to supply spare parts for the SR-X instrument. Paramit has no obligation to maintain inventory in excess of any open purchase orders or materials in excess of the amount Paramit reasonably determines will be consumed within 90 days or within the lead time of manufacturing our instrument, whichever is greater. We have an obligation to purchase any material or instruments deemed in excess pursuant to the Paramit Agreement. The price is determined according to a mutually agreed-upon pricing formula. The parties agreed to review the pricing methodology yearly or upon a material change in cost.

The Paramit Agreement is terminable by either party for convenience with written notice to the other party given at least nine months prior to the end of the then-current annual term. The agreement may also be terminated by us with three months' notice to Paramit upon the occurrence of (i) a failure of Paramit to obtain any necessary governmental licenses, registrations or approvals required to manufacture our instrument or (ii) an assignment by Paramit of its rights or obligations under the agreement without our consent. The Paramit Agreement is terminable by Paramit with 30 days' notice to us in the event of a material breach after written notice and a 60-day opportunity to cure the breach.

Columbia Tech Manufacturing Agreement

In July 2019, Akoya entered into a Manufacturing Agreement ("the CT Agreement") with Columbia Tech to manufacture its Spatial Biology instruments on an as-ordered basis. We are responsible for obtaining any regulatory approval necessary to sell the instruments. The instrument price is established by Columbia Tech and is subject to certain adjustments.

The CT Agreement is terminable by either party for convenience upon 90 days' notice to the other party. The CT Agreement may also be terminated by either party with 30 days' notice of an uncured material breach, or by us if Columbia Tech fails to deliver ordered products within 45 days of the delivery date.

Intellectual Property

Our success depends in part on our ability to obtain and maintain intellectual property protection for our products and technology.

We rely primarily on a combination of patent, copyright, trademark, and trade secret laws, as well as contractual provisions with employees and third parties, to establish and protect our intellectual property rights. Our patent strategy is to pursue broad protection for key technologies, supplemented by additional patent filings covering conceptual methods, specific aspects of current and proposed products, and forward-looking applications and technological developments, and licensing of certain patent families from third parties. We also engage in strategic analysis of our owned and licensed patent assets and pursue additional patent claims from our existing portfolio that may provide us with market and other competitive advantages. We do not rely heavily on trade secret protection but do maintain a certain amount of in-house know-how that is not disclosed publicly.

We provide products to customers, commercial, and academic collaborators pursuant to agreements with non-disclosure terms and other conditions that impose restrictions on use and disclosure. We further make use of contractual obligations that require our employees, consultants, and contractors with access to our proprietary information to execute non-disclosure, non-competition and assignment of intellectual property agreements to preserve our intellectual property

rights. We generally control access to our proprietary and confidential information through the use of internal and external controls that are subject to periodic review.

Our Simoa patent strategy protects but is not limited to: fundamental methods for detecting single molecules independent of the specific analyte to be detected, embodiments of the core technology directed to the detection of specific analytes, novel instrumentation, consumables, and manufacturing processes used in applying the invention to certain commercial products or future product opportunities and specific uses of the core technology (e.g., biomarkers and diagnostics).

Our Spatial Biology patent strategy protects but is not limited to: spatial / tissue multiplexing technologies and spatial / tissue signal amplification technologies for both protein and nucleic acids; selective labeling, RNA sequence identification and error correction visualizing tissues and compartments within cells, image, and signal processing technologies including spectral unmixing, dynamic optical correction, machine classifiers, multiple-image registration, systems and methods for whole-slide imaging, automated adjustment of imaging systems, extraction of pure spectra from sample images, focus determination in imaging systems as well as specific biomarkers or combinations thereof for use in discovery and diagnostics.

Our patent strategy is both offensive and defensive in nature, seeking to protect not only technology we currently practice but also alternative, related embodiments.

As of December 31, 2025, our active patent portfolio (owned or exclusively licensed) included approximately 154 issued patents worldwide with an additional 102 applications pending. Our owned or exclusively licensed patents and patent applications, if issued, are expected to expire between 2026 and 2043, in each case without taking into account any possible patent term adjustments or extensions and assuming payment of all appropriate maintenance, renewal, annuity, or other governmental fees.

Our core Simoa technology, directed to general methods and devices for single molecule detection, originated at Tufts University ("Tufts"), in the laboratory of Professor David Walt, who is one of the founders of Quanterix. Professor Walt and his students pioneered the single molecule array technology, including technologies that enabled the detection of single enzyme labels in arrays of microwells, thereby facilitating ultra-sensitive detection. We have exclusively licensed from Tufts the relevant patent rights related to these technologies (see the section titled "Material Licenses").

Our core Spatial Biology labeling technology PhenoCycler (formerly CODEX) originated in the laboratory of Professor Garry P. Nolan at Stanford, who is a former member of Akoya's board of directors. Patent families covering this technology are exclusively licensed from Stanford University (see the section titled "Material Licenses").

Our key Spatial Biology imaging technology PhenoImager (formerly PhenoOptics) originated at Cambridge Research and Instrumentation Inc, ("Cambridge Research"), a company that was later acquired by Caliper Life Sciences, Inc. ("Caliper Life Sciences"). Caliper Life Sciences was subsequently acquired by Perkin Elmer, Inc., subsequently known as Revvity ("PKI"). We purchased key patent assets covering this technology from PKI, Cambridge Research, and Caliper Life Sciences, and also licensed certain supplemental patents from PKI, Cambridge Research, and VisEn Medical Inc. ("VisEn Medical"). Some of the supplemental patents are exclusively licensed and others are non-exclusively licensed (see the section titled "Material Licenses").

In addition to our reliance on patent protection for our inventions, products and technologies, we also rely on trade secrets, know-how, confidentiality agreements, and continuing technological innovation to develop and maintain our competitive position. For example, some elements of our manufacturing processes, analytics techniques and processes, as well as computational-biological algorithms, and related processes and software, are based on unpatented trade secrets and know-how that are not publicly disclosed. Although we take steps to protect our proprietary information and trade secrets, including through contractual means with our employees, advisors and consultants, these agreements may be breached or may be unenforceable and we may not have adequate remedies. In addition, third parties may independently develop substantially equivalent proprietary information and techniques or otherwise gain access to our trade secrets or disclose our technology. As a result, we may not be able to meaningfully protect our intellectual property. For further discussion of the risks relating to our intellectual property, see the section titled "Risk Factors — Risks Related to our Intellectual Property."

Material Licenses

Tufts University

In June 2007, as amended in April 2013, August 2017, and September 2020, we entered into a license agreement with Tufts, pursuant to which we obtained an exclusive, worldwide license to all patent rights to the Simoa bead-based technology owned by Tufts, as well as a non-exclusive license to related know-how. The rights licensed to us are for all fields of use and are sublicensable.

Under the terms of the agreement, as amended, we paid a one-time, non-refundable upfront fee and issued Tufts shares of our stock. We are required to pay Tufts low single-digit royalties on all net sales of products and services that use the licensed technology, as well as a portion of any sublicensing revenues. We are also obligated to pay annual maintenance fees, which are fully creditable against any royalty payments made by us, and a milestone payment upon any sublicense by us. We were also required to reimburse Tufts for all patent prosecution costs.

The term of the license agreement will continue on a country-by-country basis so long as there is a valid claim of a licensed patent in such country. Tufts may terminate the agreement or convert to a non-exclusive license in the event (1) we fail to pay any undisputed amount when required and fail to cure such non-payment within 60 days after receipt of notice from Tufts, (2) we are in breach of any material provision of the agreement and fail to remedy such breach within 60 days after receipt of notice from Tufts, (3) we do not demonstrate diligent efforts to develop a product incorporating the licensed technology, (4) we are found on five separate audits to have underpaid pursuant to the terms of the agreement, (5) we cease to carry on the business related to the licensed technology either directly or indirectly, or (6) we are adjudged insolvent, make an assignment for the benefit of creditors or have a petition in bankruptcy filed for or against us that is not removed within 60 days. We may terminate the agreement at any time upon at least 60 days' written notice. Upon termination of the agreement, all rights revert to Tufts.

Stanford University

In November 2015, Akoya entered into an exclusive (equity) agreement with Stanford University ("Stanford"), pursuant to which Stanford granted Akoya an exclusive, sublicensable (subject to certain requirements), worldwide license under certain patent rights owned by Stanford relating to oligonucleotide-based biological sample labeling to make, use and sell products and services that are covered by such patent rights (the "Stanford Licensed Products") in all fields of use. The patents are related to oligonucleotide-based labeling technology, and we refer to this technology as the CODEX 1 technology.

In November 2016, the agreement was amended to include an exclusive, sublicensable (subject to certain requirements), worldwide license granted to Akoya by Stanford under additional patent rights owned by Stanford relating to oligonucleotide-based biological sample labeling to make, use, and sell products and services that are covered by such patent rights, in all fields of use (such products and services are also included in the Stanford Licensed Products). We refer to the technology disclosed in the additional patents as the CODEX 2 technology. Akoya is obligated to use commercially reasonable efforts to develop, manufacture, sell and develop markets for Stanford Licensed Products, including with respect to accomplishing specific goals with specific deadlines set forth in the agreement.

Akoya made one-time upfront payments upon the execution of the agreement and amendments. We are required to pay Stanford annual license maintenance fees in the mid-five figures. We further agreed to make certain milestone payments based on specified patent issuance and sales milestone events. We are also obligated to pay Stanford a low single-digit percentage royalty on net sales of Stanford Licensed Products and a portion of any of its sublicensing income.

Subject to Stanford's approval, we control the prosecution and maintenance of the licensed patents and, if we develop Stanford Licensed Products, we have the first right to institute a suit, or defend any declaratory judgment action, related to third-party infringement of the licensed patents.

The agreement will continue until the expiration, revocation, invalidation or abandonment of the last patent or patent application that is licensed to us, unless terminated earlier in accordance with its terms. The last licensed patent is set to expire in 2036. We may terminate the agreement at any time by providing advance written notice of at least 30 days. Stanford may terminate the agreement if we violate or fail to perform any material terms thereof or for our failure to achieve certain milestones or use commercially reasonable efforts to develop and commercialize the Stanford Licensed Products and fail to cure such violation or failure within 30 days of written notice from Stanford.

PKI/Revvity, Cambridge Research, and VisEn Medical

In September 2018, in connection with the acquisition of the Quantitative Pathology Solutions ("QPS") technology from PKI, Akoya entered into a license and royalty agreement with PKI, Cambridge Research, and VisEn Medical, (collectively, the "Licensor"), pursuant to which the Licensor granted Akoya an exclusive, nontransferable, sublicensable (subject to certain conditions), worldwide license within certain fields of use under certain patent rights and know-how owned by the Licensor to make, use, and sell products within such fields of use, as well as a similar, non-exclusive license under certain other patent rights. The licensed patents relate to methods and systems for analyzing biological samples, and in particular, slide-mounted tissue samples.

We are obligated to pay the Licensor royalties ranging from low to mid single-digit percentages on net sales of products covered by either license on a decreasing schedule that ends upon the expiration of the last valid claim of the licensed patents, at which point the agreement shall terminate and our rights and licenses thereunder shall survive on a fully-paid up, royalty-free basis. The last licensed patent is set to expire in 2036. Neither we nor the Licensor has the right to terminate the agreement prior to such expiration.

The Licensor has the first right to control the prosecution, maintenance and defense of the licensed patents. We have the first right to enforce any exclusively licensed patent with respect to third-party infringement occurring solely within its licensed field of use, and Licensor has the first right to enforce the license patents with respect to any other third-party infringement. If any exclusively licensed patent is believed to be infringed by the development, manufacture, use, offer for sale, sale or importation of a product by the third-party solely inside field of use worldwide, the Licensor has the first right to institute, prosecute and control any action or proceeding with respect to such infringement of such patent.

University of Washington

In June 2018, Akoya entered into an exclusive patent license agreement with the University of Washington (the "University"), pursuant to which the University granted Akoya an exclusive, sublicensable (subject to certain conditions), worldwide license in certain fields of use under certain patent rights owned by the University relating to technology for molecular profiling of cells and tissue specimens, to make, use and sell products that are covered by such patent rights, or the Washington Licensed Products. The licensed patents are related to the detection of biomolecules, particularly proteins and nucleic acids, in biological samples.

We are obligated to pay the University a low single-digit percentage running royalty on net sales of Washington Licensed Products, subject to certain minimum annual royalty payments and potential reductions based on a royalty-stacking allowance for certain third -party rights that are required to be obtained to make, use, sell or import Washington Licensed Products. We are also obligated to make cumulative one-time payments to the University upon the achievement of certain commercial milestones, as well as sharing a portion of any of its non-royalty sublicensing income.

We are obligated to use commercially reasonable efforts to commercialize the inventions covered by the licensed patent rights and to make and sell Washington Licensed Products as soon as practicable and maximize sales thereof, including with respect to accomplishing specific goals with specific deadlines set forth in the agreement.

The University must conduct the prosecution of the licensed patents per our instructions and at our expense, subject to certain exceptions. We have the first right to defend and enforce the licensed patents at its expense.

The agreement will expire when all licensed patent rights have terminated, unless terminated earlier in accordance with the terms thereof. The last licensed patent is set to expire in 2032. We may terminate the agreement at any time by providing advance written notice of at least 60 days. The University may terminate the agreement if we violate or fail to perform any material term thereof and fail to cure such violation or failure within 60 days of written notice from the University. In addition, the University may terminate the exclusive license agreement upon 10 days' prior written notice upon certain insolvency-related events involving us or should we challenge the validity of the licensed patents.

Competition

We compete with both established and development-stage life science companies that design, manufacture, and market instruments for proteomics discovery and clinical research applications. Companies such as Bio-Techne, MesoScale Discovery, SEER, Bio-Rad Laboratories, Alamar, Spear Bio, and others, have products for protein measurements in biofluids that compete in certain segments of the market in which we sell our products. In the spatial biology market, there

are companies, both established and early-stage, that have indicated that they are designing, manufacturing, and marketing products for, among other things, tissue analysis, single-cell analysis and spatial analysis. These companies include 10x Genomics, Vizgen, Bio-Techne, Bruker, Miltenyi Biotec, and Standard BioTools, each of which has products that compete to varying degrees with some but not all of our product solutions, as well as a number of other emerging and established companies.

Our Accelerator Laboratory competes with other research laboratories such as LabCorp, Covance, Q2 Solutions, Rules Based Medicine, Monogram Biosciences, Frontage, Thermo Fisher Scientific, and others, some of which are customers of ours. In addition, as we or our partners expand the applications for our products to include diagnostics, we expect to compete with companies such as Roche, Fujirebio, Danaher, and C2N Diagnostics. Furthermore, our technology and products are showing promise for non-invasive early disease detection, and in the future, we could experience competition from companies that develop and market imaging and other molecular detection technologies.

Many of the companies with which we compete or will compete have substantially greater resources than we have. However, we believe we are differentiated from our competitors for several reasons, including our position as a leader in a large and growing market, proprietary technologies, rigorous product development processes, scalable infrastructure, and positive customer experience. We believe our customers view our products and company favorably because of these differentiators.

The life science instrumentation and lab services industries in which we operate are highly competitive and expected to grow more competitive with the increasing knowledge gained from ongoing research and development. We believe the principal competitive factors for us include:

- sensitivity;
- cost of instruments and consumables;
- reputation among customers and key opinion leaders;
- innovation in product offerings;
- accuracy and reproducibility of results; and
- customer support infrastructure.

We believe that we are well positioned with respect to these competitive factors and expect to enhance our position through ongoing global expansion, innovative new product introductions and ongoing collaborations, and partnerships with key opinion leaders.

For further discussion of the risks we face relating to competition, see the section titled “Risk Factors — Risks Related to our Business and Industry — Our market is highly competitive, and if we cannot compete successfully with our competitors, we may be unable to increase or sustain our revenue, or achieve and sustain profitability.”

Government Regulation

The majority of our products are currently intended for research use only ("RUO") applications, although our customers may use our products to develop their own products that are subject to regulation by the FDA or CMS. Although in vitro diagnostic products intended for RUO are not currently required to obtain premarket clearance or approval by the FDA, products labeled as RUO are subject to the FDA's premarket review requirements if they are determined to be intended for use for clinical rather than non-clinical research purposes. Consequently, other than our four LDTs intended for clinical testing, our products are labeled and intended “For Research Use Only. Not for Diagnostic Procedures.”

The FDA has issued Final Guidance for Industry and Food and Drug Administration Staff on "Distribution of In Vitro Diagnostic Products Labeled for Research Use Only or Investigational Use Only" (the "RUO/IUO Guidance"). The purpose of this FDA guidance document is to provide the FDA's current thinking on when IVD products are properly labeled for RUO or for investigational use only ("IUO") and when products labeled RUO or IUO will be viewed by the FDA as intended for clinical use. The RUO/IUO Guidance explains that the FDA will review the totality of the circumstances when evaluating whether equipment and testing components are properly labeled as RUO. Merely including

a labeling statement that a product is intended for research use only will not necessarily exempt the device from the FDA's premarket notification and clearance process (510(k)), premarket approval, or other requirements, if the circumstances surrounding the distribution of the product indicate that the manufacturer intends its product to be used for clinical diagnostic use. These circumstances may include written or verbal marketing claims or links to articles regarding a product's performance in clinical applications, a manufacturer's provision of technical support for clinical validation or clinical applications, or solicitation of business from clinical laboratories, all of which could be considered evidence of intended uses that conflict with RUO labeling. We believe that our labeling and promotion of our products, including the custom assay RUO products developed by the Accelerator Laboratory, is consistent with the RUO/IUO Guidance because we have not promoted our products for clinical use in humans.

In vitro diagnostics ("IVD") tests that are commercially distributed and intended for clinical diagnostic use are regulated by the FDA as medical devices, however, the FDA has historically not regulated most laboratory tests referred to as LDTs (as discussed further below). We currently offer four LDTs through our CLIA-certified laboratory: (1) an LDT to quantitatively measure p-Tau 181 in plasma as an aid in diagnostic evaluation of Alzheimer's disease; (2) an LDT to quantitatively measure NfL in serum as an aid in the evaluation of individuals for possible neurodegenerative conditions or other causes of neuronal or central nervous system damage; (3) an LDT to quantitatively measure p-Tau 217 in plasma to aid in the diagnosis of Alzheimer's disease; and (4) a multi-marker (p-Tau 217, A β 42, A β 40, NfL, and GFAP) algorithmic LDT, under the brand name LucentAD Complete, for high accuracy plasma detection of amyloid pathology to aid in diagnostic evaluation of patients with cognitive issues who may have Alzheimer's disease.

Clinical Laboratory Improvement Amendments of 1988, Regulation of LDTs and State Regulation

We own and operate a CLIA-certified laboratory. CLIA, along with federal regulations promulgated under CLIA, apply to all clinical laboratory testing performed on humans in the United States (with the exception of research testing that does not report patient specific results). A clinical laboratory is defined by CLIA as any facility that performs examinations of specimens obtained from humans for the purpose of providing information for the diagnosis, prevention, or treatment of any disease or impairment of, or the assessment of health of, human beings. CLIA requires such laboratories to be certified by the federal government (or a CMS-approved accreditation organization) and mandates compliance with various operational, personnel, facilities administration, quality and proficiency testing requirements intended to ensure that testing services are accurate, reliable and timely. CLIA certification also is a prerequisite to be eligible to bill state and federal health care programs, as well as many private insurers, for laboratory testing services.

In addition, CLIA requires certified laboratories to enroll in an approved proficiency testing program for each of the specialties and subspecialties for which it is certified. If a laboratory fails to achieve a passing score on a proficiency test, then its CLIA certificate may be suspended, limited or revoked, or other sanctions may be imposed.

As a condition of CLIA certification, laboratories are subject to survey and inspection every other year (except laboratories with only a certificate of waiver or certificate of provider-performed microscopy procedures are not subject to biennial inspections), in addition to being subject to additional random inspections. The biennial survey is conducted by CMS, a CMS agent (typically a state agency), or a CMS-approved accreditation organization.

High complexity, CLIA-certified laboratories, such as ours, frequently develop testing procedures to provide diagnostic results to customers. These tests have been offered by high-complexity laboratories for the last few decades as LDTs, the validation and performance of which are subject to CMS oversight through its enforcement of CLIA. The FDA also has claimed that it has regulatory authority over LDTs under the agency's medical device authorities, but historically did not enforce those requirements with respect to most LDTs. However, in April 2024, the FDA ended its policy of enforcement discretion for LDTs and finalized a rule that would subject nearly all LDTs—including the LDTs that we offer—to regulation as devices under the Federal Food, Drug, and Cosmetic Act ("FDCA"). In response to FDA's final rule, several organizations filed legal challenges, asserting that the FDA rule violated the Administrative Procedure Act because LDTs do not qualify as "devices" under the FDCA. On March 31, 2025, a federal district court agreed with the challengers, holding that LDTs are laboratory services and do not meet the definition of a "device" in the FDCA. Therefore, the district court held that FDA's rule was invalid and FDA could not regulate LDTs as devices. See *ACLA v. FDA*, No. 4:24-cv-00479-SDJ (E.D. Tex. Mar. 31, 2025). The FDA did not appeal that district court decision and subsequently withdrew the LDT rule.

Notwithstanding this district court decision, and FDA's decision to not appeal, it is possible that in the future the FDA could again attempt to assert that LDTs are medical devices and must comply with the requirements of the FDCA. It is also possible that the FDA could attempt to assert that some LDTs do qualify as an "LDT." Historically, FDA defined an

LDT as a test that is intended for clinical use and is designed, manufactured, and used within a single high-complexity, CLIA-certified laboratory. If a diagnostic test commercialized by a clinical laboratory does not meet all elements of this definition, it is possible that the FDA could assert that the test is not an LDT and is subject to its jurisdiction as a medical device.

If a laboratory is out of compliance with state laws or regulations governing licensed laboratories or with CLIA, it may be subject to enforcement actions that may include suspension, limitation, or revocation of the license or CLIA certificate, assessment of financial penalties or fines, or imprisonment. Loss of a laboratory's CLIA certificate or state license may also result in the inability to receive payments from state and federal health care programs as well as private third-party payors.

When we perform clinical diagnostic testing, we may be subject to HIPAA, depending on the types of transactions we engage in, as well as additional federal and state laws that impose a variety of fraud and abuse prohibitions on healthcare providers, including clinical laboratories.

Europe/Rest of World Government Regulation

We must obtain the requisite approvals from regulatory authorities in non-U.S. countries prior to the commencement of clinical trials or marketing of our product for clinical diagnostic use in those countries. The regulations in other jurisdictions vary from those in the U.S. and may be easier or more difficult to satisfy and are subject to change. For example, in the European Union ("E.U.") new regulations recently entered into force that introduce greater regulation of medical devices and IVDs. The IVD regulation is significantly different from the IVD Directive that it replaces in that it ensures that the new requirements apply uniformly and on the same schedule across the member states, includes a risk-based classification system and increases the requirements for conformity assessment.

The CE registration for the NfL ELISA assay kit of our subsidiary Uman was approved in March 2014 under the IVD Directive. Under the IVD Directive, the assay is classified as a general IVD product and required self-certification with no involvement of a notified body/authority. The IVD Regulation introduces a new classification system for IVDs and assessment by a notified body is required for class B, C, and D products. Uman's NfL ELISA assay kit is classified as a class B product and must fully comply with (and have a CE mark issued under) the IVD Regulation by May 2027 (subject to extension of the transitional periods in the IVD Regulation). The new requirements include an ISO 13485 certification of the quality system (which Uman received in July 2018) and increased technical evidence and follow-up of performance of the specific product (e.g., clinical evidence and post-market activities). The work to evaluate and to meet the new technical requirements is on-going. When all requirements are met, a notified body will be contacted, and the certification initiated.

The NF-light Serum ELISA is currently sold only as a RUO product (not intended for diagnostic use). Work is on-going to prepare a technical file compliant with the IVD Regulation for this product as well.

Other Governmental Regulation

Privacy and Data Security Laws and Regulations

As a business with a global footprint, compliance with evolving regulations and standards in privacy and data security has resulted, and may continue to result, in increased costs, new compliance challenges, and the threat of increased regulatory enforcement activity. Our business relies on various safeguards to secure electronic transmission, storage and hosting of sensitive information, including personal information, protected health information, financial information, intellectual property, and other sensitive information related to our customers and workforce.

For example, in the U.S., the collection, maintenance, protection, use, transmission, disclosure, and disposal of certain personal information and the security of medical devices are regulated at the U.S. federal and state, international, and industry levels. U.S. federal and state laws protect the confidentiality of certain patient health information, including patient medical records, and restrict the use and disclosure of patient health information by health care providers. Privacy and Security Rules under HIPAA, as amended by HITECH, govern the use, disclosure, and security of protected health information by "Covered Entities," (which include health care providers that submit electronic claims, health plans, and health care clearinghouses) and by their "Business Associates" (which is anyone that performs a service on behalf of a Covered Entity involving the use or disclosure of protected health information and is not a member of the Covered Entity's workforce). Rules under HIPAA and HITECH include specific security standards and breach notification requirements. The U.S. Department of Health and Human Services (through the Office for Civil Rights) has direct civil enforcement

authority against Covered Entities and Business Associates with regard to both the Security and Privacy Rules. The U.S. Department of Justice has criminal enforcement authority against Covered Entities, Business Associates, and certain other entities and individuals. In addition, State Attorneys General may bring enforcement actions under HIPAA. Generally we are not a Covered Entity, however, we may operate as a Business Associate to Covered Entities under certain circumstances.

In addition, a number of states have also adopted laws and regulations that may affect our privacy and data security practices for personal information, such as comprehensive state privacy laws that govern the use, disclosure and protection of personal information, such as certain health information, Social Security numbers, and credit card account data. State consumer protection laws and consumer health privacy laws also establish privacy and security standards for use and management of personal information or consumer health data, including information related to consumers and care providers.

Outside the U.S., we are impacted by the privacy and data security requirements at the international, national, and regional level, and on an industry specific basis. Legal requirements in foreign countries relating to the collection, storage, handling and transfer of personal data and potentially intellectual property continue to evolve with increasingly strict enforcement regimes. More privacy and security laws and regulations are being adopted, and more are being enforced, with potential for significant financial penalties. In the E.U., stringent data protection and privacy rules which substantially impact the use of patient data across the healthcare industry became effective in May 2018. The General Data Protection Regulation ("GDPR") applies uniformly across the E.U. and includes, among other things, a requirement for prompt notice of data breaches to data subjects and supervisory authorities in certain circumstances and significant fines for non-compliance. The GDPR also requires companies processing personal data of individuals residing in the E.U. to comply with E.U. privacy and data protection rules. In the area of health data, the GDPR is supplemented by national laws and regulations that are less harmonized.

Because data privacy laws and regulations continue to expand, differ from jurisdiction to jurisdiction, and are subject to evolving (and at times inconsistent) governmental interpretation, compliance with these laws and regulations may require significant additional expenditures or changes in products or business that increase competition or reduce revenue. Noncompliance could result in the imposition of fines, penalties, or orders to stop noncompliant activities.

Environmental Health and Safety Laws

We are subject to federal, state, and local laws and regulations related to the protection of the environment, the health and safety of employees and the handling, transportation, storage and disposal of medical specimens, infectious and hazardous waste and radioactive materials. For example, the U.S. Occupational Safety and Health Administration ("OSHA"), has established extensive requirements relating specifically to workplace safety for healthcare employers in the U.S. This includes requirements to develop and implement multi-faceted programs to protect workers from exposure to blood-borne pathogens, including preventing or minimizing any exposure through needle stick injuries. OSHA also regulates the use of hazardous chemicals in the workplace. Likewise, we are subject to the U.S. Environmental Protection Agency and state requirements relating to the management and disposal of hazardous waste, and state requirements relating to the disposal of regulated medical waste. For purposes of transportation, some biological materials and laboratory supplies are classified as hazardous materials and are subject to regulation by one or more of the following agencies: the U.S. Department of Transportation, the U.S. Public Health Service, the U.S. Postal Service, and the International Air Transport Association. We generally use third-party vendors to dispose of regulated medical waste, hazardous waste, and radioactive materials that we may use during our research and services.

Employees and Human Capital

As of December 31, 2025, we had 450 full-time employees, of which 151 worked in sales, sales support, field service, and marketing, 24 worked in engineering and research and development, 198 worked in manufacturing and operations and 77 worked in general and administration. Of our 450 full-time employees, 382 were located in the United States and 68 were located in 14 foreign countries. None of our employees are represented by a labor union or subject to a collective bargaining agreement. Our culture emphasizes the impact our work has on the detection of neurological and other critical disorders.

Our success depends upon our ability to attract and retain highly qualified employees. Talent management is critical to our ability to execute our long-term growth strategy, and we seek to cultivate a high performing pool of talent by providing career growth, on-the-job learning opportunities and competitive total rewards.

Workforce Compensation and Pay Equity

We provide competitive compensation and benefits programs to help recruit and retain our high performing employees. We utilize third party benchmark compensation data to assist in the evaluation of market wages. Our compensation is designed to attract, retain, and motivate employees to achieve results while balancing short- and long-term company performance. All our employees are eligible for an annual bonus and/or commission plan, a matching 401(k) Plan (in the case of U.S. employees), healthcare and insurance benefits, paid time off, family leave, employee assistance programs, and behavioral health services. Additionally, some of our employees are eligible for annual equity-based grants with vesting conditions designed to award our employees' performance and encourage retention.

Company Culture

We are committed to an inclusive culture which is grounded on our values of integrity, respect, and equality. In support of our inclusive culture, we sponsor an internal group of culture champions comprised of employees and executives whose mission is to provide respectful workplace training and awareness to strengthen employee understanding and knowledge of these values. As of December 31, 2025, approximately 45% of our employees were women and approximately 33% of our employees were people of color.

We expect all personnel working at Quanterix, employees, interns, and contractors, to observe the highest levels of business ethics, integrity, and mutual respect. Our employee handbook and Corporate Code of Conduct and Ethics set forth policies that reflect our values and provide guidance for registering complaints in the event of any violation of our policies. An "open door" policy is maintained at all levels of the organization, and any form of retaliation against an employee for making a good faith complaint is strictly prohibited.

Employee Engagement and Wellness

The success of our business is dependent on the physical and mental well-being of our employees. Accordingly, we are committed to creating a safe and healthy workplace for all personnel. We provide our employees with a wide range of policies and practices to ensure an environment of physical and psychological safety and well-being.

Corporate Information

We were incorporated under the laws of the State of Delaware in April 2007 under the name "Digital Genomics, Inc.". In August 2007, we changed our name to "Quanterix Corporation". Our principal executive offices are located at 900 Middlesex Turnpike, Billerica, Massachusetts 01821, and our telephone number is (617) 301-9400.

Information Available on the Internet

Our Internet website address is www.quanterix.com. The information contained on, or that can be accessed through, our website is not a part of or incorporated by reference in this Annual Report on Form 10-K. We have included our website address in this Annual Report on Form 10-K solely as an inactive textual reference. We make available free of charge through our website our Annual Report on Form 10-K, Quarterly Reports on Form 10-Q, Current Reports on Form 8-K and amendments to those reports filed or furnished pursuant to Sections 13(a) and 15(d) of the Securities Exchange Act of 1934, as amended ("Exchange Act"). We make these reports available through the "Investors—Financial Information—SEC Filings" section of our website as soon as reasonably practicable after we electronically file such reports with, or furnish such reports to, the SEC. We also make available, free of charge on our website, the reports filed with the SEC by our executive officers, directors and 10% stockholders pursuant to Section 16 under the Exchange Act as soon as reasonably practicable after copies of those filings are provided to us by those persons. Our electronically filed reports and other information that we file with the SEC can be viewed on the SEC's website at www.sec.gov.

ITEM 1A. RISK FACTORS

The following risk factors and other information included in this Annual Report on Form 10-K should be carefully considered. The risks and uncertainties described below are not the only ones we face. Additional risks and uncertainties not presently known to us or that we presently deem less significant may also impair our business operations. Please see page ii of this Annual Report on Form 10-K for a discussion of some of the forward-looking statements that are qualified by these risk factors. If any of the following risks occur, our business, financial condition, results of operations, and future growth prospects could be materially and adversely affected.

Risk Factor Summary

- Our quarterly and annual operating results and cash flows have fluctuated in the past and might continue to fluctuate, which could cause the value of our common stock to fluctuate or decline significantly.
- We have incurred annual losses since we were formed and expect to incur losses in the future. We cannot be certain that we will achieve or sustain profitability.
- Inherent limitations associated with our internal control over financial reporting could result in material misstatements in our financial statements.
- Integrating our business with that of Akoya may be more difficult, costly or time-consuming than expected and we may fail to realize the remaining anticipated benefits of the acquisition, which may adversely affect our business results and negatively affect the value of our common stock.
- We may not achieve the remaining expected cost savings and related benefits from our cost reduction actions, and the consequences of those actions may adversely impact our business.
- Recent turnover in our board of directors and management could result in changes in our strategic plan, product focus, and investment priorities, and we might not realize the anticipated benefits from any such changes.
- If our products fail to achieve and sustain sufficient market acceptance, our revenue will be adversely affected.
- Sales of our Simoa-based assays for neurological indications have become increasingly important to our business, and there can be no assurance that we will be able to continue to generate meaningful revenue from the sale of such products.
- We may not be successful in penetrating the diagnostics market.
- The sales cycle for our instruments can be lengthy and variable, which makes it difficult for us to forecast revenue and other operating results.
- Purchase of our instruments by our customers requires a significant capital investment which can impact sales in times of constrained spending.
- Because a significant portion of our revenue comes from a few large customers, any significant decrease in sales to these customers, due to industry consolidation or otherwise, could harm our operating results.
- Our long-term results depend upon our ability to improve existing products and introduce and market new products successfully.
- We may experience delays in launching and commercializing our next-generation instruments, on our anticipated timeline, which could adversely affect our business, financial condition, and results of operations.
- Defects or other quality issues in our products could lead to unforeseen costs, product recalls, adverse regulatory actions, negative publicity, and litigation, including product liability claims, any of which could cause customers to decide not to purchase our products, harm our reputation, and negatively affect our sales, operating results and financial condition.

- We generate a substantial portion of our revenue internationally and we expect this will continue in the future; as a result, our business is subject to various risks relating to our international activities, which could adversely affect our business, operating results, and financial condition.
- Our reliance on distributors for sales of our products outside of the United States could impact our revenue.
- We rely on single contract manufacturers for certain key instruments. If any of these manufacturers should fail to perform, or not perform satisfactorily, our ability to supply these instruments would be negatively and adversely affected, which could have a material adverse effect on our business, financial condition, results of operations, and reputation.
- We rely on a limited number of suppliers or, in some cases, one supplier, for some of our materials and components used in our consumable products and certain of our instruments, and we may not be able to find replacements or immediately transition to alternative suppliers if any of these suppliers fail to perform, which could have a material adverse effect on our business, financial condition, results of operations, and reputation.
- We face significant competition in the life sciences research and diagnostic markets.
- Changes in U.S. government policies, including the impacts of tariffs and potential reductions in federal research funding, could adversely affect our business. Similarly, there is substantial uncertainty regarding how the administration's initiatives might impact the FDA, its implementations of laws, regulations, policies and guidance and its personnel, which could prevent, limit or delay development and regulatory approval of our future diagnostic products.
- If the FDA determines that our products are subject to regulation as medical devices, if Congress enacts new laws that subject our LDT to FDA regulation, if we seek to market our products for clinical diagnostic or health screening use, or if we continue to expand our product, technology and service offerings and the applications and uses of our products into new fields, we expect to become subject to additional government regulations, and the regulatory approval and maintenance process would be expensive, time-consuming and uncertain both in timing and in outcome.
- Our products may in the future be subject to product recalls that could harm our reputation, business, and financial results.
- U.S. legislative, FDA, or global regulatory reforms may make it more difficult and costly for us to obtain any required regulatory approval of our product candidates and to manufacture, market, and distribute our products after approval is obtained.
- If we do not comply with governmental regulations applicable to our CLIA-certified laboratory, we may not be able to continue our laboratory operations or continue offering our LDTs.
- We expect to rely on third parties in conducting any required future studies of diagnostic products that may be required by the FDA or other regulatory authorities, and those third parties may not perform satisfactorily.
- If diagnostic procedures that are enabled by our technology are subject to unfavorable pricing regulations or third-party coverage and reimbursement policies, our business could be harmed.
- We depend on our information technology systems, and any failure of these systems could harm our business.
- Cybersecurity breaches, loss of data and other disruptions could compromise sensitive information related to our business or prevent us from accessing critical information and expose us to liability, which could adversely affect our business and our reputation.
- Inability to adapt to and effectively incorporate potential advantages of artificial intelligence ("AI") could negatively impact our ability to compete, and inability to manage the risks of AI could expose us to liability or put us at a disadvantage.

- If we are unable to protect our intellectual property, our ability to maintain any technological or competitive advantage over our competitors and potential competitors may be reduced and our business may be harmed.

Risks Related to our Financial Condition and Financial Reporting Matters

Our quarterly and annual operating results and cash flows have fluctuated in the past, and our operating results may continue to fluctuate, which could cause the value of our common stock to fluctuate or decline significantly.

Numerous factors, many of which are outside of our control, may cause or contribute to significant fluctuations in our quarterly and annual operating results. These fluctuations may make financial planning and forecasting difficult. In addition, one or more of such factors may cause our revenue or operating expenses in one period to be disproportionately higher or lower relative to the others, and comparing our operating results on a period-to-period basis might not be meaningful. Investors should not rely on our past results as indicative of our future performance. Moreover, our stock price might be based on expectations of future performance that are unrealistic or that we might not meet and, if our revenue or operating results fall below the expectations of investors or securities analysts, the price of our common stock could decline significantly.

We have incurred annual losses since our formation, and we expect to incur losses in the future. We cannot be certain that we will achieve or sustain profitability.

Quanterix incurred net losses of \$107.2 million, \$38.5 million, and \$28.4 million for the years ended December 31, 2025, 2024, and 2023, respectively. Akoya incurred net losses of \$55.4 million, and \$63.3 million for the years ended December 31, 2024, and 2023, respectively. We cannot predict if or when we will achieve profitability or if we will be able to sustain profitability once achieved. We expect that our operating losses will continue into 2026 as we execute our growth strategy. We may incur significant losses in the future for a number of reasons, many of which are beyond our control, including any need for incremental investments to fund our strategic objectives, unanticipated delays or obstacles in the integration of Akoya's business with ours, the market acceptance of our products, competitive products, future product development and our market penetration and margins, and other risks described in this Annual Report on Form 10-K and our other reports filed with the SEC.

Inherent limitations associated with, our internal control over financial reporting could result in material misstatements in our financial statements.

In our Annual Reports on Form 10-K for the years ended December 31, 2022, 2023 and 2024, we disclosed certain material weaknesses in our internal control over financial reporting relating to the operating effectiveness of our internal controls. We have remediated these previously disclosed material weaknesses.

Our efforts to maintain effective internal control over financial reporting, are ongoing; however, there are inherent limitations in all control systems and no evaluation of controls can provide absolute assurance that all deficiencies have been detected. We cannot assure you that additional material weaknesses in our internal control over financial reporting will not arise or be identified in the future. If we are unable to maintain the effectiveness of our internal control over financial reporting or our disclosure controls and procedures, we could lose investor confidence in the accuracy and completeness of our financial reports, the market price of our common stock could decline, and we could be subject to regulatory scrutiny, civil, or criminal penalties, or litigation. Future failure to maintain effective internal control over financial reporting could also result in financial statements that do not accurately reflect our financial condition or results of operations, may result in material misstatements in our financial statements, and may also restrict our future access to the capital markets.

We have incurred significant expense and dedicated significant internal resources to address the previously disclosed material weaknesses, and our remediation efforts and activities have required attention and focus from management. There can be no assurance that we will not identify additional significant deficiencies or material weaknesses that will impair our ability to report our financial condition and results of operations accurately or on a timely basis.

Our ability to use net operating losses to offset future income may be subject to certain limitations.

As of December 31, 2025, we had forecasted federal net operating loss ("NOLs") carryforwards to offset future taxable income of approximately \$597.7 million, of which approximately \$111.1 million begin to expire in 2026. A lack of future taxable income would adversely affect our ability to utilize these NOLs. In addition, under Section 382 of the

Internal Revenue Code of 1986, as amended (the “Code”), a corporation that undergoes an “ownership change” is subject to limitations on our ability to utilize our NOLs to offset future taxable income. We may have already experienced ownership changes as defined under Section 382 of the Code. Depending on the timing of any future utilization of our NOLs, the amount that can be utilized each year may be limited as a result of such previous ownership changes. In addition, future changes in our stock ownership, including changes that may be outside of our control, could result in additional ownership changes under Section 382 of the Code. Our NOLs may also be impaired under similar provisions of state law. We have recorded a full valuation allowance related to our NOLs and other deferred tax assets due to the uncertainty of the ultimate realization of the future benefits of those assets.

Risks Related to our Business

Combining the businesses of Quanterix and Akoya may be more difficult, costly or time-consuming than expected and we may fail to realize the anticipated benefits of the acquisition, which may adversely affect our business results and negatively affect the value of our common stock.

The success of our Akoya acquisition will depend on, among other things, our ability to realize the anticipated benefits, synergies and efficiencies from combining the businesses of Quanterix and Akoya. This success will depend on, among other factors, our ability to integrate our business with the business of Akoya. If we are not able to successfully integrate Akoya’s business into ours within the anticipated time frame, or at all, the anticipated synergies, efficiencies and other benefits of the acquisition may not be realized fully, or at all, or may take longer to realize than expected.

There can be no assurances that the Quanterix and Akoya businesses can be integrated successfully. It is possible that the integration process could result in the loss of key employees, the disruption of either company’s or both companies’ ongoing businesses, inconsistencies in standards, controls, procedures and policies, unexpected integration issues, higher than expected integration costs and an overall post-completion integration process that takes longer than originally anticipated. The challenges involved in this integration, which will be complex and time-consuming, also include the following:

- combining the businesses of Quanterix and Akoya, including respective operations and corporate functions, and meeting our capital requirements in a manner that permits us to achieve efficiencies anticipated to result from the acquisition, the failure of which would result in the anticipated benefits of the acquisition not being realized in the time frame currently anticipated or at all;
- integrating, retaining and, where applicable, cross-training personnel from the two companies;
- integrating the offerings and services available to customers;
- integrating each company’s technologies and technologies licensed by them from third parties;
- identifying and eliminating redundant and underperforming functions and assets;
- harmonizing each company’s operating practices, employee development and compensation programs, internal controls and other policies, procedures and processes;
- maintaining existing relationships with each company’s customers, service providers, partners, vendors and suppliers, and leveraging relationships with such third parties;
- addressing possible differences in business backgrounds, corporate cultures and management philosophies;
- consolidating each company’s administrative and information technology infrastructure;
- coordinating geographically dispersed organizations; and
- effecting actions that may be required in connection with obtaining regulatory or other governmental approvals.

Certain key employees have left the Company following completion of the acquisition, and it is possible that additional key employees may decide not to remain with us. If additional key employees terminate their employment, or if an insufficient number of employees or sales representatives are retained to maintain effective operations, our business

activities may be adversely affected and management's attention may be diverted away from integrating our business with Akoya's business, which may cause our business to suffer. In addition, we may not be able to locate suitable replacements for any key employees that leave or offer employment to potential replacements on reasonable terms. Moreover, there could be disruptions to or distractions for the workforce and management, including disruptions in the integration of employees into the combined workforce. We may not be able to attract or retain key employees to the same extent that we and Akoya were able to attract or retain their respective employees in the past.

In addition, at times the attention of management and our resources may be focused on the integration of the two businesses and diverted from day-to-day business operations or other opportunities that may have been beneficial, which may disrupt our business. An inability to realize the full extent of the anticipated benefits of the acquisition, as well as any delays or higher than expected integration costs encountered in the integration process, could have an adverse effect on our revenues, level of expenses and operating results, which may adversely affect the value of our common stock.

We may fail to achieve the remaining expected cost savings and related benefits from our cost reduction actions, and the consequences of those actions may adversely impact our business.

In 2025, we announced certain actions to reduce our operating costs, including reductions in our workforce. Certain of these cost reduction actions were part of our integration plan to realize anticipated synergies and other benefits of the Akoya acquisition. On an annualized basis, we realized approximately \$74.0 million of synergies in 2025. As we complete the integration of Akoya, we expect to take additional steps in 2026 to further reduce our operating expenses, with a goal of reducing our annualized operating expenses by a total of \$85.0 million. We incurred expenses of approximately \$8.0 million in the aggregate associated with the 2025 reductions in our workforce, substantially all of which are cash expenditures incurred in 2025 for severance and benefits. There is no guarantee that these cost reduction actions will result in the anticipated savings or other economic benefits, and we may incur unanticipated charges or make payments that were not previously contemplated. Additionally, these actions:

- may result in the loss of institutional knowledge and expertise;
- may disrupt or restrain the scope of our business activities; and
- may make it more difficult to attract and retain qualified personnel, whose duties may be expanded to include those of employees whose positions were eliminated in the reductions in force.

If we are unable to realize the anticipated benefits from these reductions in force and other operating expense reductions, or if we experience significant adverse consequences from these reductions in force and other operating expense reductions, our business, financial condition, and results of operations may be materially adversely affected.

Recent transitions in our board of directors and management could result in changes in our strategic plan, product focus, and investment priorities, and we might not realize the anticipated benefits from any such changes.

In the second half of 2025, four new directors joined our board of directors and four previous members of our board of directors departed, including Dr. David Walt, our founding scientist. In addition, in January 2026 Dr. Masoud Taloue left the Company and Everett Cunningham became our Chief Executive Officer and joined our board of directors. These board of directors and management transitions could result in a re-assessment of, and potential shift in, our strategic plan, and could thus lead to changes in our product focus and investment priorities. While any such changes would be aimed at accelerating our growth, it is possible that such changes in strategy could result in delays in new product launches, disruptions in operations, or increased expenses and may not achieve the benefits we anticipate.

If our products fail to achieve and sustain sufficient market acceptance, our revenue will be adversely affected.

Our success depends on our ability to develop and market products that are recognized and accepted by our customers and potential customers as reliable, enabling and cost-effective. Continued market acceptance of our Simoa technology platform and products, PhenoCycler and PhenoImager platform and products, and other platforms and products we may develop in the future will depend on many factors, including our ability to convince potential customers that our technologies are attractive alternatives to other available technologies. If we are unable to continue to motivate customers to use our current technologies or other technologies we may develop in the future, adoption of our technologies may be slowed and our ability to retain and grow our customer base and increase our revenue would be adversely affected.

Sales of our Simoa-based assays for neurological indications have become increasingly important to our business, and there can be no assurance that we will be able to continue to generate meaningful revenues from the sale of such products.

Neurology has been one of our primary focus areas for commercialization of our Simoa technology and the services that we provide to our customers. Sales from neurological-related biomarkers have become an increasingly important part of our business. There can be no assurance that we will continue to derive meaningful revenues from the sale of our neurological products, from services related to neurodegenerative conditions or from sales of instruments driven by customers desiring access to our technology for work relating to neurological conditions. The adoption by our customers of competitive technologies for detecting biomarkers of neurodegenerative conditions could negatively impact our revenues and have a material adverse effect on our business.

We may not be successful in penetrating the diagnostics market.

We believe our Simoa technology has the capability to enable the development of a new category of less-invasive diagnostic tests that could replace current invasive, expensive, and inconvenient diagnostic methods. Accordingly, we have begun to expand into the diagnostics market. Transitioning from research use only to also serving the diagnostics market entails significant risks, including:

- significant investments in product development, scaling manufacturing processes, marketing and sales activities, regulatory compliance, reimbursement and billing activities, and infrastructure to support the foregoing;
- navigating complex regulatory frameworks, including but not limited to FDA regulations and equivalent agencies internationally;
- competition from products that may offer superior performance, pricing, or convenience, and prevent us from penetrating target markets effectively; and
- challenges associated with obtaining adequate reimbursement from government healthcare programs and private insurers.

Further, our progress in penetrating the diagnostics market may be slower than we intend and may require a substantially larger investment than we expect. If we are unable to manage these risks effectively, our efforts to penetrate the diagnostics market may be unsuccessful, and our business, operating results and financial condition could suffer.

The sales cycle for our instruments can be lengthy and variable, which makes it difficult for us to forecast revenue and other operating results.

The sales process for our instruments generally involves numerous interactions with multiple individuals within an organization, and often includes in-depth analysis by potential customers of our technology and products and a lengthy review process. Our customers' evaluation processes often involve a number of factors, many of which are beyond our control. As a result of these factors, the capital investment required to purchase our systems, and the budget cycles of our customers, the time from initial contact with a customer to our receipt of a purchase order can vary significantly. Given the length and uncertainty of our sales cycle, we have in the past experienced, and expect in the future to experience, fluctuations in our sales on a period-to-period basis. In addition, any failure to meet customer expectations could result in customers choosing to retain their existing systems, using existing assays not requiring capital equipment, or purchasing systems other than ours.

Purchase of our instruments by our customers requires a significant capital investment which can impact sales in times of constrained spending.

The purchase of our instruments requires a significant investment by our customers, and a reduction in capital spending by potential customers can result in lower instrument sales. During periods of constrained capital spending, potential instrument customers may instead choose to engage our Accelerator Laboratory or an outside lab, or may use another instrument platform that they already have or that is less expensive than our instruments. We believe that in 2025 rapidly changing macro-economic conditions resulting from reductions in US federal research funding, reductions in research and development spending by larger pharmaceutical customers, and the impacts of new import tariffs contributed

to a constrained capital funding environment that resulted in softness in instrument sales. We expect these macro-economic factors to continue into 2026.

Because a significant portion of our revenue comes from a few large customers, any significant decrease in sales to these customers, due to industry consolidation or otherwise, could harm our operating results.

For the year ended December 31, 2025, Quanterix's top five customers accounted for approximately 14% of its total revenue. The loss of a significant amount of business from one or more of our major customers would have a material adverse effect on our business. There can be no assurance that there will not be a loss or reduction in business from one or more of our major customers. In addition, we cannot assure that net sales from customers that have accounted for significant net sales in the past, either individually or as a group, will reach or exceed historical levels in any future period.

Our long-term results depend upon our ability to improve existing products and introduce and market new products successfully and timely.

We generally sell our products in industries that are characterized by rapid technological changes, frequent new product introductions and changing industry standards. Accordingly, our business is dependent on the continued improvement of our existing products and our development of new products utilizing our current technologies or other technology we develop or acquire in the future. As we introduce new products or refine, improve or upgrade versions of existing products, we cannot predict the level of market acceptance or the amount of market share these products will achieve, if any. We cannot guarantee that we will not experience material delays in the introduction of new products in the future. In addition, introduction of new products could result in a decrease in revenues from existing products. Consistent with our strategy of offering new products and product refinements, we have invested substantial capital on research and development, and we expect to continue to use a substantial amount of capital for product research and development. Our research and development initiatives can be costly and time-consuming, and they may fail to achieve the intended benefits. If we do not develop new products and product enhancements based on technological innovation on a timely basis, our products may become obsolete over time and our revenues, cash flow, profitability and competitive position will suffer.

We may experience delays in launching and commercializing our next-generation instruments, on our anticipated timeline, which could adversely affect our business, financial condition, and results of operations.

There are various risks that could delay or prevent the successful launch and commercialization of our next generation instruments. These risks include, but are not limited to, unforeseen technical challenges, supply chain disruptions, and delays in manufacturing. Many of these risks are beyond our control. If we experience significant delays in refreshing our product portfolio and launching new products, our ability to generate revenue and achieve market adoption may be adversely impacted. Delays or setbacks could also allow competitors to introduce alternative solutions, erode our market position, or negatively affect customer confidence in our product pipeline. Additionally, if development costs exceed our expectations, or if we are unable to successfully commercialize the platform, our financial condition and results of operations could suffer.

Defects or other quality issues in our products could lead to unforeseen costs, product recalls, adverse regulatory actions, negative publicity, and litigation, including product liability claims, any of which could cause customers to decide not to purchase our products, harm our reputation, and negatively affect our sales, operating results and financial condition.

Our products are complex and may contain undetected errors or defects, especially when first introduced or as new versions or new products are released. We have periodically experienced product delays and product quality issues, and, accordingly, we have in the past devoted, and will continue to devote, funding and resources to technology development, quality assurance and manufacturing initiatives designed to ensure or improve quality. However, there can be no assurance that we will be successful in our efforts to manufacture products at a level of quality necessary for our customers or to avoid our products containing undiscovered defects or quality issues. Additionally, reduction in personnel who service our instruments may result in service delays, instrument downtime and customer dissatisfaction. Defects, errors or quality issues in our products may discourage customers from purchasing our products and could harm our reputation. We may also be subject to warranty claims and litigation involving claims for damages or incur additional costs, in each case due to errors or defects in our products. In addition, if we do not meet industry or quality standards, if applicable, our products may be subject to recall. A material liability claim, recall or other occurrence that harms our reputation or decreases market acceptance of our products could harm our business and operating results.

Use of our products or services by us or a customer for diagnostic purposes could result in a product liability claim alleging that one of our products contained a design or manufacturing defect that resulted in the failure to adequately perform, leading to death or injury. A product liability claim could result in substantial damages and be costly and time-consuming to defend, either of which could materially harm our business or financial condition. We cannot guarantee that our product liability insurance would adequately protect our assets from the financial impact of defending a product liability claim. Any product liability claim brought against us, with or without merit, could increase our product liability insurance rates or prevent us from securing insurance coverage in the future.

Our reliance on distributors for sales of our products outside of the United States could impact our revenue.

We have established distribution agreements for our instruments and related consumable products with distributors in a number of foreign countries, including Australia, Brazil, China, the Czech Republic, India, Hong Kong, Israel, Japan, New Zealand, Qatar, Saudi Arabia, Singapore, South Africa, South Korea, Taiwan and the United Arab Emirates. We intend to continue to grow our business internationally, and to do so we must attract additional distributors and retain existing distributors to maximize the commercial opportunity for our products. There is no guarantee that we will be successful in attracting or retaining desirable sales and distribution partners or that we will be able to enter into such arrangements on favorable terms. Distributors may not commit the necessary resources to market and sell our products to the level of our expectations or may choose to favor marketing the products of our competitors. If current or future distributors do not perform adequately, or if we are unable to enter into effective arrangements with distributors in particular geographic areas, we may not realize long-term international revenue growth. In addition, if our distributors fail to comply with applicable laws and ethical standards, including anti-bribery laws, this could damage our reputation and could have a significant adverse effect on our business and our revenues.

We generate a substantial portion of our revenue internationally and we expect this will continue in the future; as a result, our business is subject to various risks relating to our international activities, which could adversely affect our business, operating results and financial condition.

For the years ended December 31, 2025, 2024, and 2023, approximately 40%, 37%, and 38%, respectively, of Quanterix's total revenue was generated from customers located outside of North America. We believe that a substantial percentage of our future revenue will continue to come from international sources as we expand our overseas operations and develop opportunities in additional areas. Engaging in international business involves a number of difficulties and risks, including:

- difficulties and costs of staffing and managing foreign operations;
- required compliance with existing and changing U.S. or foreign regulatory requirements and laws;
- a shortage of high-quality salespeople and distributors;
- pricing pressure that we may experience internationally;
- difficulties in enforcing our intellectual property rights and in defending against third-party threats and intellectual property enforcement actions against us or any of our distributors, suppliers or collaborators;
- reduced or varied protection for intellectual property rights in some countries;
- compliance with anti-bribery laws, such as the U.S. Foreign Corrupt Practices Act, data privacy requirements, such as the E.U. GDPR, labor laws and anti-competition regulations;
- export or import restrictions and supply chain disruptions;
- laws and business practices favoring local companies;
- longer payment cycles and difficulties in enforcing agreements and collecting receivables through certain foreign legal systems;
- restrictions on the activities of foreign agents, representatives and distributors;

- foreign currency exchange rate fluctuations;
- the imposition of new trade restrictions;
- potentially adverse tax consequences, the impacts of new and changing levels of tariffs, customs charges, bureaucratic requirements and other trade barriers;
- the imposition of U.S. or international sanctions against a country, company, person or entity with whom we do business that would restrict or prohibit continued business with the sanctioned country, company, person or entity;
- the impact of political and economic instability and conflict, which could lead to uncertainty and instability in global markets; and
- scrutiny of foreign tax authorities which could result in significant fines, penalties and additional taxes being imposed on us.

If we are unable to manage these risks effectively, our business, operating results and financial condition will suffer.

We rely on single contract manufacturers for several of our key instruments, and we expect to that to continue in the future. If any of these manufacturers should fail to perform, or not perform satisfactorily, our ability to supply these instruments would be negatively and adversely affected, which could have a material adverse effect on our business, financial condition, results of operations and reputation.

We currently rely on a single contract manufacturer, STRATEC Biomedical AG (“STRATEC”), an analytical and diagnostic systems manufacturer located in Germany, to manufacture and supply our Simoa HD-X instruments. We currently rely on a second single contract manufacturer, Paramit Corporation (“Paramit”), a contract manufacturer located in California, to manufacture and supply all of our SR-X instruments. Additionally, we rely on a third single contract manufacturer, Columbia Tech, to manufacture our PhenoCycler and PhenoImager instruments. Certain of our manufacturing contracts do not commit our contract manufacturers to supply quantities beyond the amounts included in our forecasts or commit them to carry inventory or make available any particular quantities. Accordingly, we may not be able to obtain adequate supplies for these products in a timely manner or on commercially reasonable terms. If any of these manufacturers are not able to supply instruments to us in the quantities and/or timeframe required, our business, financial condition, results of operations and reputation could be materially adversely affected.

In the event it becomes necessary to utilize a different contract manufacturer for an instrument, we would experience additional costs, delays and difficulties in doing so as a result of needing to identify and enter into an agreement with a new supplier as well as needing to prepare such new supplier to meet the logistical requirements associated with manufacturing our instruments. These additional costs, delays and difficulties could have a material adverse effect on our business, financial condition, results of operations and reputation. We may also experience additional costs and delays in the event we need access to or rights under any intellectual property of our contract manufacturers. In addition, certain of the components used in our instruments are sourced by these manufacturers from limited or sole suppliers. If our manufacturers were to lose such suppliers, there can be no assurance that they would be able to identify or enter into agreements with alternative suppliers on a timely basis on acceptable terms, if at all. An interruption in our ability to sell and deliver instruments to customers could occur if our manufacturers encounter delays or difficulties in securing these components, or if the quality of the components supplied do not meet specifications, or if they cannot then obtain an acceptable substitute. If any of these events occur, our business, financial condition, results of operations and reputation could be materially adversely affected.

We rely on a limited number of suppliers or, in some cases, one supplier, for some of our materials and components used in our consumable products and services and certain of our instruments, and we may not be able to find replacements or immediately transition to alternative suppliers if any of these suppliers fail to perform, which could have a material adverse effect on our business, financial condition, results of operations and reputation.

We rely on limited or sole suppliers for certain reagents and other materials and components that are used in our consumable products and services and in certain of our instruments. While we have long-term contracts with some critical suppliers, we do not have contracts with all suppliers and instead rely on periodically forecasting our needs for such materials and entering into standard purchase orders with our suppliers. In addition, our use of many of the materials used

in our consumable products is limited to research use only. As we expand into diagnostic applications for our products, we will need to secure diagnostic rights to such materials. If we were to lose suppliers or were unable to secure required rights for materials from suppliers, we may be unable to identify or enter into agreements with alternative suppliers on a timely basis or on acceptable terms, if at all. An interruption in our operations could occur if we encounter delays or difficulties in securing these materials or any required rights to these materials, if the quality of the materials supplied do not meet our requirements, or if we cannot then obtain an acceptable substitute. The time and effort required to qualify a new supplier and ensure that the new materials provide the same or better quality results could result in significant additional costs. Any such interruption could significantly affect our business, financial condition, results of operations and reputation.

We could be adversely affected by violations of the U.S. Foreign Corrupt Practices Act and other worldwide anti-bribery laws by us or our agents.

We are subject to the U.S. Foreign Corrupt Practices Act (the “FCPA”), which prohibits companies and individuals from corruptly making payments, directly or indirectly through third parties, to non-U.S. government officials for the purpose of obtaining or retaining business or securing any other improper advantage. We are also subject to the FCPA’s accounting provisions, which requires us to keep accurate books and records and to maintain a system of internal accounting controls sufficient to assure management’s control, authority and responsibility over our assets. Our reliance on independent distributors to sell our products internationally demands a high degree of vigilance in maintaining our policy against participation in corrupt activity, because there are circumstances under which we could be held responsible for their actions. Other U.S. companies in the medical device and pharmaceutical fields have faced criminal penalties under the FCPA for allowing their distributors and other third parties to deviate from appropriate practices in doing business with these individuals. We are also subject to similar anti-bribery laws in the jurisdictions in which we operate, including the United Kingdom’s Bribery Act of 2010, which also prohibits commercial bribery and makes it a crime for companies to fail to prevent bribery. These laws are complex and far-reaching in nature, and any violations of these laws, or allegations of such violations, could disrupt our operations, involve significant management distraction, involve significant costs and expenses, including legal fees, and could result in a material adverse effect on our business, prospects, financial condition, or results of operations. We could also incur severe penalties, including criminal and civil penalties, disgorgement and other remedial measures.

The life sciences research and diagnostic markets are highly competitive. If we fail to effectively compete, our business, financial condition and operating results will suffer.

We face significant competition in the life sciences research and diagnostic markets. We currently compete with both established and early-stage companies that design, manufacture and market systems and consumable supplies. Many of our current competitors have competitive advantages over us, including:

- greater name and brand recognition;
- substantially greater financial and human resources;
- broader product lines;
- larger sales forces and more established distributor networks;
- more substantial intellectual property portfolios;
- larger and more established customer bases and relationships; and
- better established, larger scale and lower cost manufacturing capabilities.

We cannot guarantee that our products will compete favorably or that we will be successful in the face of increasing competition from new products and technologies introduced by our existing competitors or new companies entering our markets. In addition, we cannot guarantee that our competitors do not have or will not develop products or technologies that currently or in the future will enable them to produce competitive products with greater capabilities or at lower costs than ours. Any failure to compete effectively could materially and adversely affect our business, financial condition and operating results.

Integrating any business, product or technology we acquire can be expensive and time-consuming and can disrupt and adversely affect our ongoing business, including product sales, and distract our management.

In addition to Akoya, we have acquired, and may in the future acquire other businesses, products or technologies as well as pursue strategic alliances, joint ventures, technology licenses or investments in complementary businesses. Our ability to successfully integrate any business, product or technology we acquire depends on a number of factors, including, but not limited to, our ability to:

- minimize the disruption and distraction of our management and other employees in connection with the integration of any acquired business, product or technology;
- avoid acquisition of unanticipated liabilities related to acquired companies;
- maintain and increase sales of our existing products;
- establish or manage the transition of the manufacture and supply of any acquired product;
- identify and add the necessary sales, marketing, manufacturing, regulatory and other related personnel, capabilities and infrastructure that are required to successfully integrate any acquired business, product or technology;
- manage the transition and migration of acquired personnel and all commercial, financial, legal, regulatory and other pertinent information relating to any acquired business, product or technology;
- comply with legal, regulatory and contractual requirements applicable to any acquired business, product or technology; and
- maintain and extend intellectual property protection for any acquired product or technology.

If we are unable to perform the above functions or otherwise effectively integrate any acquired businesses, products or technologies, our business, financial condition and operating results will suffer. Also, the anticipated benefit of any acquisition may not materialize. Future acquisitions or dispositions could result in potentially dilutive issuances of our equity securities, the incurrence of debt, contingent liabilities or amortization expenses or write-offs of goodwill, any of which could harm our financial condition. We cannot predict the number, timing or size of future joint ventures or acquisitions, or the effect that any such transactions might have on our operating results.

Risks Related to Government Regulation and Diagnostic Product Reimbursement

Changes in U.S. government policies, including reductions in federal research funding and the impacts of tariffs, are adversely affecting our business, though the full extent of the impact is uncertain. Similarly, there is substantial uncertainty regarding how the current administration's initiatives might impact the FDA, its implementations of laws, regulations, policies and guidance and its personnel, which could prevent, limit or delay development and regulatory approval of our future diagnostic products.

The U.S. government has suspended or withheld disbursement of funds under certain federal research grants (or certain components of grants) and curtailed the grant of new awards, including funding and grants from the National Institutes of Health ("NIH"). These actions are negatively impacting spending within our industry and causing uncertainty, which adversely impacted our business and our financial outlook for 2026. Certain of our customers, including academic institutions and research organizations, may depend in whole or in part on federal grants to advance their medical research activities. Any prolonged suspensions or reductions in such funding could slow innovation, delay collaborations, and limit the adoption of new technologies that contribute to our business growth.

Other recent policy actions, including the imposition of, and changes in, tariffs on imported materials and goods from certain foreign countries, may also have an adverse impact on our business. The U.S. government has announced and/or implemented a range of tariffs on imports from other countries, resulting in retaliatory tariffs by certain countries. Higher tariffs on materials, goods and components used by us or our suppliers may raise production costs and could disrupt the supply chain. Because tariffs will likely increase the costs of materials, goods and components, we expect we will need

to absorb the costs in some cases and/or increase the prices of certain of our products. This could adversely impact demand for our products and our competitive positioning.

Further, there is substantial uncertainty with regard to the regulatory environment under the current administration. For example, certain initiatives have manifested to date in the form of personnel measures that could impact the FDA's ability to hire and retain key personnel, which could result in delays or limitations on our ability to obtain guidance from the FDA on our diagnostic products in development and obtain the requisite regulatory approvals in the future. The administration could issue or promulgate executive orders, regulations, policies or guidance that adversely affect us or our customers or create a more challenging or costly environment to pursue the development of new products. Alternatively, state governments may attempt to address or react to changes at the federal level with changes to their own regulatory frameworks in a manner that is adverse to our operations.

If these or similar policy changes continue or expand, or if we or our customers become negatively impacted by future governmental orders, regulations, policies or guidance, we may face increased costs, demand for our products could be impacted and there could be a material adverse effect on us and our business. Though the risks referenced above have already adversely impacted our business to some extent, the full impact of funding actions and tariffs on us and on our business partners remains highly uncertain and volatile. We cannot predict the full extent of these impacts, but any prolonged disruption could further adversely affect our business, financial condition, and results of operations.

If the FDA determines that our products are subject to regulation as medical devices, if Congress enacts new laws that subject our LDT to FDA regulation, if we seek to market our products for clinical diagnostic or health screening use, or if we continue to expand our product, technology and service offerings and the applications and uses of our products into new fields, we expect to become subject to additional government regulations, and the regulatory approval and maintenance process would be expensive, time-consuming and uncertain both in timing and in outcome.

We focused initially on the life sciences research market. This includes offering products for use by laboratories associated with academic and governmental research institutions, as well as pharmaceutical, biotechnology and contract research companies. Accordingly, the majority of our products are labeled as "Research Use Only" ("RUO"), which indicates that they are intended for use in research, not for clinical diagnostic use. None of our products are currently regulated as medical devices. However, if our products labeled as RUO are marketed or promoted for the clinical diagnosis of diseases or medical conditions, it is possible that the FDA would assert that such products qualify as medical devices under the Federal Food, Drug and Cosmetic Act ("FDCA") and must comply with all of the requirements of that statute (as discussed further below).

Further, while we focused initially on the life sciences research market and RUO products only, our strategy includes expanding our product line to encompass products that are intended to be used for the diagnosis of disease, including LDTs and in vitro diagnostic ("IVD") devices, either alone or in collaboration with third parties. IVD products are subject to regulation by the FDA, or comparable international agencies, as medical devices. If regulated as devices, our products would be subject to a number of additional requirements including requirements for regulatory clearance or approval of such products before they can be marketed and postmarket requirements, as discussed further below.

The process of obtaining regulatory clearances to market a medical device can be costly and time consuming, and we or our collaborators may not be able to obtain these clearances or approvals on a timely basis, if at all. Although some devices are exempt from premarket clearance or approval requirements, for many devices, commercial distribution of a new medical device is lawful only after the device has received clearance under Section 510(k) of the FDCA, or is the subject of an approved Premarket Approval ("PMA"). The FDA will clear marketing of a lower risk medical device through the 510(k) process if the manufacturer demonstrates that the new product is substantially equivalent to a legally marketed predicate device, which can include pre-amendment, 510(k)-exempt, 510(k) cleared products, or PMA-approved products that have subsequently been down-classified. If the FDA determines that the device is not "substantially equivalent" to a predicate device, or if the device is novel, it is automatically classified into Class III, and the device sponsor must then fulfill the much more rigorous premarketing requirements of the PMA approval process, or seek classification of the device through the de novo classification process. The PMA process is more costly, lengthy and uncertain than the 510(k) clearance process. A PMA application must be supported by extensive data, including, but not limited to, technical, preclinical, clinical trial, manufacturing and labeling data, to demonstrate to the FDA's satisfaction the safety and efficacy of the device for our intended use.

If any of our products are subject to medical device regulation, we would be subject to a substantial number of additional requirements for medical devices, including establishment registration, device listing, quality system regulations — which cover the design, testing, production, control, quality assurance, labeling, packaging, servicing, sterilization (if required), and storage and shipping of medical devices (among other activities) — product labeling, advertising, recordkeeping, post-market surveillance, post-approval studies, adverse event reporting, and correction and removal (recall) regulations. One or more of the products we may develop using our technology may also require clinical trials in order to generate the data required for a PMA, de novo classification request or 510(k) premarket notification. Further, if we sell devices for diagnostic purposes, or if our products are reimbursed by federal or state health care programs, we may in turn be subject to additional healthcare regulation and enforcement by the applicable government agencies. Such laws and regulations include, without limitation, state and federal anti-kickback, fraud and abuse, false claims, data privacy and security and transparency and reporting requirements for payments and transfers of value to physicians and certain other healthcare professionals. Complying with these requirements may be time-consuming and expensive. We may be required to expend significant resources to ensure ongoing compliance with the FDA and other healthcare regulations. Failure to comply with these requirements may subject us to a range of enforcement actions, such as warning letters, injunctions, civil monetary penalties, criminal prosecution, recall and/or seizure of products, and revocation of marketing authorization, as well as significant adverse publicity. If we fail to obtain, or experience significant delays in obtaining, regulatory approvals for IVD products, such products may not be able to be launched or successfully commercialized in a timely manner, or at all.

LDTs are tests that are offered as services by Clinical Laboratory Improvement Amendments of 1988 (“CLIA”)-certified high complexity clinical laboratories and designed and used within a single laboratory. In July 2022, we launched an LDT to quantitatively measure p-Tau 181 in plasma as an aid in diagnostic evaluation of Alzheimer’s disease, and subsequently have launched additional LDTs to quantitatively measure other biomarkers in serum as an aid in the evaluation of individuals for possible neurodegenerative conditions or other causes of neuronal or central nervous system damage. The FDA has sought to assert, and could again attempt to assert, that LDTs are medical devices and must comply with the requirements of the FDCA. It is also possible that the FDA could attempt to assert that some LDTs do not qualify as an “LDT,” which the FDA historically defined as a test that is intended for clinical use and is designed, manufactured, and used within a single high-complexity, CLIA-certified laboratory. If a diagnostic test commercialized by a clinical laboratory does not meet all elements of this definition, it is possible that the FDA could assert that the test is not an LDT and is subject to its jurisdiction as a medical device. Further, if our Accelerator Laboratory fails to comply with state laws or regulations governing licensed laboratories or with CLIA, we may be subject to enforcement actions that may include suspension, limitation, or revocation of the license or CLIA certificate, and assessment of financial and other penalties.

Foreign jurisdictions have laws and regulations similar to those described above, which may adversely affect our ability to market our products as planned in such countries. The number and scope of these requirements are increasing. As in the United States, the cost and time required to comply with regulatory requirements may be substantial, and there is no guarantee that we will obtain the necessary authorization(s) required to make our products commercially viable. In addition, the imposition of foreign requirements may also have a material adverse effect on the commercial viability of our operations.

Our products may in the future be subject to product recalls that could harm our reputation, business and financial results.

The FDA and similar foreign governmental authorities have the authority to require the recall of commercialized products, including RUO products, in the event of material deficiencies or defects in design or manufacture. In the case of the FDA, the authority to require a mandatory recall of a medical device must be based on an FDA finding that there is a reasonable probability that the device would cause serious injury or death. The FDA may also request or exert pressure on manufacturers to initiate voluntary recalls where deficiencies or defects present risks, even in situations where the FDA cannot mandate a recall. Manufacturers may, under their own initiative, recall a product if any material deficiency in a device is found. A government-mandated or voluntary recall by us or one of our distributors could occur as a result of component failures, manufacturing errors, design or labeling defects or other deficiencies and issues. Recalls of any of our products would divert managerial and financial resources and have an adverse effect on our financial condition, results of operations, and reputation.

U.S. legislative, FDA or global regulatory reforms may make it more difficult and costly for us to obtain any required regulatory approval of our product candidates and to manufacture, market and distribute our products after approval is obtained.

From time to time, legislation is drafted and introduced in Congress that could significantly change the statutory provisions governing the regulatory approval, manufacture and marketing of regulated products or the reimbursement thereof. For example, in December 2022, Congress enacted the Food and Drug Omnibus Reform Act of 2022 (“FDORA”). FDORA reauthorized the FDA to collect device user fees and contained substantive amendments to the device provisions of the FDCA, including imposing new cybersecurity and clinical trial requirements for devices. From time to time, Congress has also considered legislation to impose a new FDA regulatory framework for all diagnostics, including IVD devices and LDTs. Congress could also amend CLIA to impose new obligations on laboratories, including obligations related to LDTs. Any new statutes, regulations or revisions or reinterpretations of existing regulations may impose additional costs or lengthen review times of future products. In addition, regulations and guidance issued by FDA and the Center for Medicare & Medicaid Services (“CMS”) are often revised or reinterpreted in ways that may significantly affect our business and our products. It is impossible to predict whether legislative changes will be enacted or regulations, guidance or interpretations changed, and what the impact of such changes, if any, may be. Any change in the laws or regulations that govern the clearance and approval processes relating to our current and future products could make it more difficult and costly to obtain clearance or approval for new products, or to produce, market and distribute existing products. Significant delays in receiving clearance or approval, or the failure to receive clearance or approval for our new products would have an adverse effect on our ability to expand our business.

In addition, in the E.U. the IVD regulation (the “IVD Regulation”) replaced the European directive for IVD medical devices (the “IVD Directive”) in 2022. The IVD Regulation includes a risk-based classification system and increases the requirements for conformity assessment. The CE registration for UmanDiagnostics AB’s (“Uman’s”) NfL enzyme-linked immunosorbent assay (“ELISA”) kit for cerebral spinal fluid was approved in March 2014 under the IVD Directive. Under the IVD Directive the assay is classified as a general IVD product, and required self-certification with no involvement of a notified body/authority. However, under the IVD Regulation, assessment by a notified body is required for class B, C and D products. Uman’s NfL ELISA kit for cerebrospinal fluid (“CSF”) is classified as a class B product and must fully comply with (and have a CE mark issued under) the IVD Regulation by May 2027. Among other requirements, the IVD Regulation requires an ISO 13485 certification of our quality system (which Uman received in July 2018), technical documentation addressing, among other things, a product’s design, manufacturing, performance, and risk, and follow-up assessments of the performance of the product (e.g. clinical evidence and post-market activities). The work to evaluate and to meet the new technical requirements is ongoing.

Our failure to continue to comply with applicable foreign regulatory requirements, including those administered by authorities of the European Economic Area (“EEA”) countries, could result in enforcement actions against us, including refusal, suspension or withdrawal of our CE Certificates of Conformity by our notified body, which could impair our ability to market products in the EEA in the future.

If we do not comply with governmental regulations applicable to our CLIA-certified laboratory, we may not be able to continue our laboratory operations or continue offering our LDTs.

CLIA is a federal law that regulates clinical laboratories that perform examination of human specimens for the purpose of providing information for the diagnosis, prevention or treatment of any disease or impairment of, or the assessment of health of, human beings. The operation of our CLIA-certified laboratory is subject to regulation by numerous federal, state and local governmental authorities in the United States. This laboratory holds a CLIA certificate of compliance for high-complexity testing and is licensed by California, Maryland, Massachusetts, New York, Pennsylvania and Rhode Island. We may seek to obtain other state licenses if required in the future. Failure to comply with federal or state regulations or changes in those regulatory requirements could result in a substantial curtailment or even prohibition of the operations of our laboratory and could have an adverse effect on our business. To maintain CLIA certification, laboratories are subject to survey and inspection every two years. Moreover, CLIA inspectors may make unannounced inspections of these laboratories. If we were to lose our CLIA certification or any required state licenses, whether as a result of a revocation, suspension or limitation, we could have a material adverse effect on our business.

We expect to rely on third parties in conducting any required future studies of diagnostic products that may be required by the FDA or other regulatory authorities, and those third parties may not perform satisfactorily.

We do not have the ability to independently conduct clinical trials or other studies that may be required to obtain FDA and other regulatory clearance or approval for future diagnostic products. Accordingly, we expect that we would rely on third parties, such as clinical investigators, CROs, consultants, and collaborators to conduct such studies if needed. For example, we are currently working with the Global Alzheimer's Platform Foundation on prospective clinical trials for our neurological assays. Our reliance on these third parties for clinical and other development activities would reduce our control over these activities. If these third parties do not successfully carry out their contractual duties or regulatory obligations or meet expected deadlines, if the third parties need to be replaced or if the quality or accuracy of the data they obtain is compromised, we may not be able to obtain regulatory clearance or approval.

If diagnostic procedures that are enabled by our technology are subject to unfavorable pricing regulations or third-party coverage and reimbursement policies, our business could be harmed.

The ability of us, our customers or our collaborators to commercialize diagnostic tests based on our technology, including LDTs that we have launched or may launch in the future, will depend in part on the extent to which coverage and reimbursement for these tests will be available from government health care programs, private health insurers and other third-party payors. In the United States, the principal decisions about reimbursement for new technologies are often made by the CMS. CMS often uses third parties, such as Medicare Administrative Contractors (MACs) and other third parties to make coverage decisions. These third parties play an important role in determining coverage and reimbursement for diagnostic products and services, including reimbursement for LDTs.

Private payors often follow CMS's reimbursement policies to a substantial degree. CMS recently approved a reimbursement rate of \$897 for the LucentAD Complete multiplex test; however, it can be difficult to predict what CMS will decide with respect to reimbursement in any particular case. However, a significant trend in the U.S. healthcare industry and elsewhere is cost containment. Government authorities and third-party payors have attempted to control costs by limiting coverage and the amount of payments for particular products and procedures. We cannot be sure that coverage will be available for any diagnostic tests based on our technology, and, if coverage is available, the level of reimbursement. Payor coverage and reimbursement decisions may impact the demand for those tests. If coverage is not available or the reimbursement amount is inadequate, any tests for which marketing authorization is received may not be able to be successfully commercialized.

For example, as a result of the Protecting Access to Medicare Act ("PAMA") of 2014, CMS payments for certain diagnostic testing services, such as the performance of LDTs and other tests, may be reduced. Although reductions resulting from PAMA were delayed in prior years, unless new legislation is enacted, cuts are scheduled to resume in 2027.

Risks Related to our Operations

We depend on our information technology systems, and any failure of these systems could harm our business.

We depend on information technology and telecommunications systems to operate our business. Our enterprise software systems affect a broad range of business processes and functional areas, including, for example, systems handling human resources, accounting, manufacturing, inventory control, financial controls and reporting, sales administration, and other infrastructure operations. We maintain preventive and detective security controls and seek to enhance such controls by, for example, augmenting the monitoring and alerting functions, network design, and automatic countermeasure operations of our technical systems. We also periodically assess the adequacy of our hardware and systems and are planning to upgrade hardware and systems where appropriate. These information technology and telecommunications systems support a variety of functions, including manufacturing operations, quality control, customer service support, finance, and other general administrative activities.

Information technology and telecommunications systems are vulnerable to damage from a variety of sources, including telecommunications, systems or network failures, malicious human acts, and natural disasters. Moreover, despite network security and back-up measures, some of our servers are potentially vulnerable to physical or electronic break-ins, computer viruses, and similar disruptive problems. Despite the precautionary measures we have taken to prevent unanticipated problems that could affect our information technology and telecommunications systems, those measures may be inadequate and failures or significant downtime of our information technology or telecommunications systems or those used by our third-party suppliers could prevent us from operating our business and managing the administrative aspects of

our business. Loss of data or a material delay in our access to our data due to a security breach or other interruption could also prevent us from operating our business. Any disruption or loss of information technology or telecommunications systems on which critical aspects of our operations depend could have an adverse effect on our business.

Cybersecurity breaches, loss of data and other disruptions could compromise sensitive information related to our business or prevent us from accessing critical information and expose us to liability, which could adversely affect our business and our reputation.

In the ordinary course of our business, we collect and store sensitive data, and intellectual property and proprietary business information owned or controlled by us or our customers. This data encompasses a wide variety of business-critical information including research and development information, operational information, commercial information, and business and financial information. We face four primary risks relative to protecting this critical information: loss of access; inappropriate disclosure; inappropriate modification; and inadequate monitoring of our controls over the first three risks.

The secure processing, storage, maintenance, and transmission of critical information is vital to our operations and business strategy, and we devote significant resources to protecting such information. Although we take measures to protect sensitive information from unauthorized access or disclosure, our information technology and infrastructure may be vulnerable to attacks by hackers or viruses, breaches, interruptions due to employee error, malfeasance, faulty password management, lapses in compliance with privacy and security mandates, or other disruptions. The risk of a security breach or disruption, particularly through cyber-attack or cyber intrusion, including by computer hackers, foreign governments, and cyber terrorists, has generally increased as the number, intensity and sophistication of attempted attacks and intrusions from around the world have increased. Our IT networks and related systems are essential to the operation of our business and our ability to perform day-to-day operations. Although we make efforts to maintain the security and integrity of these types of IT networks and related systems, and we have implemented various measures to manage the risk of a security breach or disruption, no security measure is infallible and there can be no assurance that our security efforts and measures will be effective or that attempted security breaches or disruptions will not be successful or damaging. Our information technology systems may have vulnerabilities, and we may not have the resources or technical sophistication to anticipate or prevent rapidly evolving types of cyberattacks, such as ransomware attacks. Although we have experienced cybersecurity incidents from time to time that have not had a material adverse effect on our business, financial condition, or results of operations, there can be no assurance that a cyber-attack, security breach, or other cybersecurity incident will not have a material adverse effect on us in the future. A significant cyber incident, including system failure, security breach, disruption by malware or other damage, could interrupt or delay our operations, result in a violation of applicable cybersecurity and privacy and other laws, damage our reputation, cause a loss of customers, expose sensitive customer data, or give rise to monetary fines and other penalties, which could be significant.

Third parties may attempt to fraudulently induce employees or other persons into disclosing usernames, passwords, or other sensitive information, which may in turn be used to access our information systems, commit identity theft or carry out other unauthorized or illegal activities. Any such breach could compromise our networks and the information stored there could be accessed, publicly disclosed, lost, or stolen. We engage third-party vendors and service providers to store and otherwise process some of our data, including sensitive and personal information. Our vendors and service providers may also be the targets of the risks described above, including cyberattacks, malicious software, phishing schemes, and fraud. Our ability to monitor our vendors and service providers' data security is limited, and third parties may be able to circumvent any security measures, resulting in the unauthorized access to, misuse, disclosure, loss or destruction of our data, including sensitive and personal information, and disruption of our or third-party service providers' systems. We and our third-party service providers may face difficulties in identifying, or promptly responding to, potential security breaches and other instances of unauthorized access to, or disclosure or other loss of, information. Any hacking or other attack on our or our third-party service providers' or vendors' systems, and any unauthorized access to, or disclosure or other loss of, information suffered by us or our third-party service providers or vendors, or the perception that any of these have occurred, could result in legal claims or proceedings, loss of intellectual property, liability under laws that protect the privacy of personal information, negative publicity, disruption of our operations and damage to our reputation, which could divert our management's attention from the operation of our business and materially and adversely affect our business, revenues and competitive position.

Any security breach or interruption, as well as any action by us or our employees or contractors that might be inconsistent with the rapidly evolving data privacy and security laws and regulations applicable within the United States and elsewhere where we conduct business, could result in enforcement actions by state or federal governments or foreign governments, liability or sanctions under data privacy laws that protect personally identifiable information, regulatory

penalties, other legal proceedings such as but not limited to private litigation, the incurrence of significant remediation costs, disruptions to our development programs, business operations and collaborations, diversion of management efforts and damage to our reputation. Because of the rapidly moving nature of technology and the increasing sophistication of cybersecurity threats, our measures to prevent, respond to and minimize such risks may be unsuccessful.

In addition, our insurance may be insufficient to cover our losses resulting from cyber-attacks, breaches, or other interruptions, and any incidents may result in loss of, or increased costs of, such insurance. The successful assertion of one or more large claims against us that exceed available insurance coverage, the occurrence of changes in our insurance policies, including premium increases or the imposition of large deductible or co-insurance requirements, or denials of coverage, could have a material adverse effect on our business, including our financial condition, results of operations and reputation.

We are currently subject to, and may in the future become subject to additional, U.S. federal and state and international laws and regulations imposing obligations on how we collect, store and process personal information. Our actual or perceived failure to comply with such obligations could harm our business. Ensuring compliance with such laws could also impair our efforts to maintain and expand our future customer base, and thereby decrease our revenue.

In the ordinary course of our business, we collect, store, transfer, use or process sensitive data, including personally identifiable information of employees and others, and intellectual property and proprietary business information owned or controlled by us and other parties. The secure processing, storage, maintenance, and transmission of this critical information are vital to our operations and business strategy. We are, and may increasingly become, subject to various laws and regulations, as well as contractual obligations, relating to data privacy and security in the jurisdictions in which we operate. The regulatory environment related to data privacy and security is increasingly rigorous, with new and constantly changing requirements applicable to our business, and enforcement practices are likely to remain uncertain for the foreseeable future. These laws and regulations may be interpreted and applied differently over time and from jurisdiction to jurisdiction, and it is possible that they will be interpreted and applied in ways that may have a material adverse effect on our business, financial condition, results of operations and prospects.

In the United States, various federal and state regulators, including governmental agencies like the Federal Trade Commission, have adopted, or are considering adopting, laws and regulations concerning personal information and data security. Certain state laws may be more stringent or broader in scope, or offer greater individual rights, with respect to personal information than federal, international, or other state laws, and such laws may differ from each other, all of which may complicate compliance efforts. For example, the California Consumer Privacy Act (the “CCPA”), which increases privacy rights for California residents and imposes obligations on companies that process their personal information, came into effect on January 1, 2020. Among other things, the CCPA requires covered companies to provide disclosures to California consumers regarding the processing of their personal data, as well as data protection and privacy rights, including the ability to opt-out of certain sales or sharing of personal information. The CCPA provides for civil penalties for violations, as well as a private right of action for certain data breaches that result in the loss of personal information. This private right of action may increase the likelihood of, and risks associated with, data breach litigation. In November 2020, California also passed the California Privacy Rights Act (the “CPRA”), which became effective on January 1, 2023 and significantly expands the CCPA, including by introducing additional obligations such as data minimization and storage limitations and granting additional rights to consumers. More recently, other states, including Connecticut, Colorado, Utah and Virginia have passed comprehensive state data privacy laws, and states like Washington and Nevada have enacted consumer health privacy laws. Most of these laws are enforced by state attorneys general, but there is the potential for private actions by plaintiffs in some circumstances under certain laws, including under Washington’s consumer health data privacy law. In addition, laws in all 50 U.S. states require businesses to provide notice to consumers whose personal information has been disclosed as a result of a data breach. State laws are changing rapidly and there is discussion in the U.S. Congress of a new comprehensive federal data privacy law to which we would become subject if it is enacted. These and future laws and regulations may increase our compliance costs and potential liability.

Furthermore, regulations promulgated pursuant to the Health Insurance Portability and Accountability Act of 1996 (“HIPAA”) establish privacy and security standards that limit the use and disclosure of individually identifiable health information (known as “protected health information”) and require the implementation of administrative, physical, and technological safeguards to protect the privacy of protected health information and ensure the confidentiality, integrity and availability of electronic protected health information. Determining whether information constitutes protected health information has been handled in compliance with applicable privacy standards and our contractual obligations can require complex factual and statistical analyses and may be subject to changing interpretation. Although we take measures to protect sensitive data from unauthorized access, use or disclosure, our information technology and infrastructure may be

vulnerable to attacks by hackers or viruses or breached due to employee error, malfeasance or other malicious or inadvertent disruptions. Any such breach or interruption could compromise our networks and the information stored there could be accessed by unauthorized parties, manipulated, publicly disclosed, lost, or stolen. Any such access, breach or other loss of information could result in legal claims or proceedings, and liability under federal or state laws that protect the privacy of personal information, such as, if applicable, the HIPAA, the Health Information Technology for Economic and Clinical Health Act of 2009 (“HITECH”), and regulatory penalties. Where such laws are applicable, notice of breaches must be made to affected individuals, the Secretary of the Department of Health and Human Services, and for extensive breaches, notice may need to be made to the media. Such a notice could harm our reputation and our ability to compete.

Outside of the United States, many countries have privacy and data security laws and regulations concerning the collection and use of personal data, including but not limited to the GDPR and China’s Personal Information Protection Law. The GDPR, which governs the collection and use of personal data in the E.U. and is wide-ranging in scope, imposes several requirements relating to the consent of the individuals to whom the personal data relates, the information provided to the individuals, the security and confidentiality of the personal data, data breach notification and the use of third-party processors in connection with the processing of the personal data. The GDPR also imposes strict rules on the transfer of personal data out of the E.U. to the United States, enhances enforcement authority and imposes large penalties for noncompliance, including the potential for fines of up to €20 million or 4% of the annual global revenues of the infringer, whichever is greater. While we have taken steps to comply with the GDPR, including reviewing our security procedures and entering into data processing agreements with relevant contractors, we cannot guarantee that our compliance efforts will be fully successful.

Risks Related to Intellectual Property

Inability to adapt to and effectively incorporate potential advantages of artificial intelligence (“AI”) could negatively impact our ability to compete, and inability to manage the risks of AI could expose us to liability or put us at a disadvantage.

Artificial intelligence could disrupt certain aspects of our business and the markets in which we operate and change use of technology in ways that are not yet known. Currently, AI tools are not integral to our product design and development, data analysis, or product manufacturing. If we are not able to adapt and effectively incorporate potential advantages of AI in our business, it may negatively impact our ability to compete.

AI technologies are subject to a variety of laws, including intellectual property, privacy, data protection and cybersecurity, consumer protection, competition, and equal opportunity laws, and are expected to be subject to increased regulation and new laws or new applications of existing laws. Such laws and regulations may present a variety of compliance risks. The use of AI may also result in litigation, ethical concerns, and other legal and business risks. If we are not able to effectively manage the risks of AI, we may suffer harm to our results of operations and reputation.

If we are unable to protect our intellectual property, our ability to maintain any technological or competitive advantage over our competitors and potential competitors may be reduced, and our business may be harmed.

We rely on patent protection as well as trademark, copyright, trade secret and other intellectual property rights protection and contractual restrictions to protect our proprietary technologies, all of which provide limited protection and may not adequately protect our rights or permit us to gain or keep any competitive advantage. If we fail to protect our intellectual property, third parties may be able to compete more effectively against us, we may lose our technological or competitive advantage, or we may incur substantial litigation costs in our attempts to recover or restrict use of our intellectual property.

Our currently pending or future patent applications may not result in granted patents, and we cannot predict how long it will take for such patents to be granted. It is possible that, for any of our patents that have granted or that may grant in the future, others will design around our patented technologies. Further, other parties may challenge any patents granted to us and courts or regulatory agencies could hold our patents to be invalid or unenforceable. We may not be successful in defending challenges made against our patents and patent applications. Any successful third-party challenge to our patents could result in the unenforceability or invalidity of such patents, or to such patents being interpreted narrowly or otherwise in a manner adverse to our interests. Our ability to establish or maintain a technological or competitive advantage over our competitors may be diminished because of these uncertainties. For these and other reasons, our intellectual property may not provide us with any competitive advantage. To the extent our intellectual property offers inadequate protection, or is found to be invalid or unenforceable, we would be exposed to a greater risk of direct competition. If our intellectual

property does not provide adequate coverage over our products and protection against our competitors' products, our competitive position could be adversely affected, as could our business.

In addition to pursuing patents on our technology, we also rely upon trademarks, trade secrets, copyrights and unfair competition laws, as well as license agreements and other contractual provisions, to protect our intellectual property and other proprietary rights. Despite these measures, any of our intellectual property rights could be challenged, invalidated, circumvented or misappropriated. In addition, we take steps to protect our intellectual property and proprietary technology by entering into confidentiality agreements and intellectual property assignment agreements with our employees, consultants, corporate partners and, when needed, our advisors. Such agreements may not be enforceable or may not provide meaningful protection for our trade secrets or other proprietary information in the event of unauthorized use or disclosure or other breaches of the agreements, and we may not be able to prevent such unauthorized disclosure. Moreover, if a party having an agreement with us has an overlapping or conflicting obligation to a third party, our rights in and to certain intellectual property could be undermined. Monitoring unauthorized disclosure is difficult, and we do not know whether the steps we have taken to prevent such disclosure are, or will be, adequate. If we were to enforce a claim that a third party had illegally obtained and was using our trade secrets, it would be expensive and time-consuming, the outcome would be unpredictable, and any remedy may be inadequate. In addition, courts outside of the United States may be less willing to protect trade secrets.

Some of our owned and in-licensed intellectual property has been discovered through government-funded programs and thus is subject to federal regulations such as "march-in" rights, certain reporting requirements, and a preference for U.S. industry. Compliance with such regulations may limit our exclusive rights, subject us to expenditure of resources with respect to reporting requirements, and limit our ability to contract with non-U.S. manufacturers.

Some of the intellectual property rights we own and have in-licensed has been generated through the use of U.S. government funding and are therefore subject to certain federal regulations. For example, some of the issued U.S. patents we own and all of the intellectual property rights licensed to us under our license agreement with Tufts have been generated using U.S. government funds. As a result, the U.S. government has certain rights to intellectual property embodied in our current or future products pursuant to the Bayh-Dole Act of 1980. These U.S. government rights in certain inventions developed under a government-funded program include a non-exclusive, non-transferable, irrevocable worldwide license to use inventions for any governmental purpose. In addition, the U.S. government has the right to require us to grant exclusive, partially exclusive, or non-exclusive licenses to any of these inventions to a third party if the government determines that: (i) adequate steps have not been taken to commercialize the invention; (ii) government action is necessary to meet public health or safety needs; or (iii) government action is necessary to meet requirements for public use under federal regulations (also referred to as "march-in rights"). The U.S. government also has the right to take title to these inventions if we fail, or the applicable licensor fails, to disclose the invention to the government, elect title, and file an application to register the intellectual property within specified time limits. In addition, the U.S. government may acquire title to these inventions in any country in which a patent application is not filed within specified time limits. Intellectual property generated under a government funded program is also subject to certain reporting requirements, compliance with which may require us, or the applicable licensor, to expend substantial resources. In addition, the U.S. government requires that any products embodying the subject invention or produced through the use of the subject invention be manufactured substantially in the United States. The manufacturing preference requirement can be waived if the owner of the intellectual property can show that reasonable but unsuccessful efforts have been made to grant licenses on similar terms to potential licensees that would be likely to manufacture substantially in the United States or that under the circumstances domestic manufacture is not commercially feasible. This preference for U.S. manufacturing may limit our ability to license the applicable patent rights on an exclusive basis under certain circumstances.

We depend heavily on intellectual property licensed from third parties, including our license agreements with Tufts University for our Simoa bead-based technology, Stanford University for our PhenoCycler product, and the University of Washington and Revvity (formerly Perkin Elmer, Inc.) for our PhenoImager product, and our licensors may not always act in our best interest. If such owners do not properly or successfully obtain, maintain or enforce the patents underlying such licenses, or if they retain or license to others any competing rights, our competitive position and business prospects may be adversely affected.

We are heavily dependent on patents, know-how and proprietary technology licensed from others. For example, we are a party to an agreement with Tufts pursuant to which we in-license patents for our Simoa bead-based technology. We are also a party to license agreements with Stanford, pursuant to which we in-license key patents and patent applications for our proprietary PhenoCycler product, as well as possible future products and other technology used in our

PhenoCycler product, and with the University of Washington and Revvity pursuant to which we have in-licensed important patents that protect key aspects of our current and future PhenoImager technologies.

Our success will depend in part on the ability of our licensors to obtain, maintain, protect and enforce patent protection for our licensed intellectual property, in particular, those patents to which we have secured exclusive rights. Our licensors may not successfully prosecute the patent applications licensed to us. If we or our licensors fail to adequately protect our licensed intellectual property, our ability to commercialize our current or potential products and technology could suffer. In addition, we may not have the right to control the maintenance, prosecution, preparation, filing, enforcement, defense and litigation of patents and patent applications that we license from other third parties. For example, in our agreement with Revvity, we do not maintain control over the prosecution and maintenance of the licensed patents. We thus cannot be certain that activities such as the maintenance and prosecution by our licensors have been or will be conducted consistent with our best interests or in compliance with applicable laws and regulations, or will result in valid and enforceable patents and other intellectual property rights. It is possible that our licensors' infringement proceedings or defense activities may be less vigorous than had we conducted them ourselves or may not be conducted in accordance with our best interests. If our licensors fail to maintain such patents or patent applications, determine not to pursue litigation against other companies that are infringing these patents, pursue litigation less aggressively than we would, or lose rights to those patents or patent applications, the rights we have licensed may be reduced or eliminated, and our right to develop and commercialize any current or future product or potential products that are the subject of such licensed rights and our right to exclude third parties from commercializing competing products could be adversely affected. Any of the foregoing could have a material adverse effect on our business, financial condition, results of operations and prospects.

Additionally, certain of our licenses may not provide us with exclusive rights to use the licensed intellectual property and technology, or may not provide us with exclusive rights to use such intellectual property and technology in all relevant fields of use and in all territories in which we may wish to develop or commercialize our technology and products in the future. In addition, the intellectual property portfolio licensed to us by our licensors, including certain intellectual property licensed by Stanford, at least in some respects, may be used by such licensors. Thus, patents licensed to us could be put at risk of being invalidated or interpreted narrowly in litigation filed by or against our licensors or another licensee or in administrative proceedings brought by or against our licensors or another licensee in response to such litigation or for other reasons. As a result, we may not be able to prevent competitors or other third parties from developing and commercializing competitive products, including in territories covered by our licenses.

Our rights to develop and commercialize our products and technologies are subject, in part, to the terms and conditions of licenses granted to us by others.

We have in-licensed certain intellectual property rights from third parties, including Tufts, with respect to our Simoa bead-based technology, Stanford and the University of Washington, with respect to our PhenoCycler platform, and Revvity, Cambridge Research and VisEn Medical Inc. with respect to our PhenoImager platform, and we may license intellectual property rights from others in the future. See the section titled "Part I. Item 1. Business — Licenses" for more information regarding such agreements. If, for any reason, our license agreements are terminated or we otherwise lose the rights associated with such licenses, it could adversely affect our business. Our current and any future license agreements may impose various development, commercialization, funding, diligence, sublicensing, insurance, patent prosecution and enforcement or other obligations on us, as well as milestone, royalty, annual maintenance and other payment obligations. If we breach any material obligations, or use the intellectual property licensed to us in an unauthorized manner, or if, in spite of our efforts, a collaborator or licensor concludes that we have materially breached our obligations under such agreement, we may be required to pay damages and the licensor may have the right to terminate the license, which could result in us being unable to develop, manufacture and commercialize products that are covered by the licensed technology or having to negotiate new or reinstated licenses on less favorable terms, or enable a competitor or other third-party to gain access to the licensed technology.

Licensing of intellectual property is of high importance to our business and involves complex legal, business and scientific issues. Disputes may arise between us and our licensors regarding intellectual property subject to a license agreement, including:

- the scope of rights granted under the license agreement and other interpretation-related issues;
- our compliance with reporting and financial obligations under our license agreements;

- whether and the extent to which our products and technologies infringe on, misappropriate or otherwise violate intellectual property of the licensor that is not subject to the license agreement;
- our right to sublicense the applicable intellectual or proprietary rights to third parties;
- our diligence obligations with respect to the use of the licensed technology in relation to our development and commercialization of our products and technologies, and what activities satisfy those diligence obligations;
- our right to transfer or assign the license;
- the inventorship and/or ownership of patents, inventions, know-how and other intellectual property and proprietary rights resulting from activities performed by our licensors, us and our partners; and
- the priority of invention of patented technology.

These agreements may be complex, and certain provisions in such agreements may be susceptible to multiple interpretations. The resolution of any contract interpretation disagreement that may arise could narrow what we believe to be the scope of our rights to the relevant intellectual property or technology, or increase what we believe to be our financial or other obligations under the relevant agreement. Moreover, if disputes over intellectual property that we have licensed prevent or impair our ability to maintain our licensing arrangements on acceptable terms, we may not be able to successfully develop and commercialize the affected product or potential products. In addition, certain of our agreements may limit or delay our ability to consummate certain transactions, may impact the value of those transactions, or may limit our ability to pursue certain activities. Any of the foregoing could have a material adverse effect on our business, financial condition, results of operations and prospects.

Additionally, we may find it necessary or prudent to acquire or obtain licenses from third-party intellectual property holders. However, we may be unable to acquire or secure such licenses to any intellectual property rights from third parties that we identify as necessary for our products or any future products we may develop. The acquisition or licensing of third-party intellectual property rights is a competitive area, and our competitors may pursue strategies to acquire or license third-party intellectual property rights that we may consider attractive or necessary. Our competitors may have a competitive advantage over us due to their size, capital resources and greater development and commercialization capabilities. In addition, companies that perceive us to be a competitor may be unwilling to assign or license rights to us. We also may be unable to acquire or license third-party intellectual property rights on terms that would allow us to make an appropriate return on our investment or at all. Any of the foregoing could have a material adverse effect on our business, financial condition, results of operations and prospects.

If we or any of our partners are sued for infringing intellectual property rights of third parties, the resulting litigation would be costly and time-consuming, and an unfavorable outcome in that litigation could have a material adverse effect on our business.

Our success depends on our ability to develop, manufacture, market and sell our products and perform our services without infringing upon the proprietary rights of third parties. As part of a business strategy to impede our successful commercialization and entry into new markets, competitors have claimed, and may claim in the future, that our products and/or services infringe their intellectual property rights and have suggested, and may suggest in the future, that we enter into license agreements. We believe any such claims made to date are without merit. However, even if such claims are without merit, they could divert the attention of our management and technical personnel and we could incur substantial costs in defending against or settling such claims. Any adverse ruling by a court or administrative body, or perception of an adverse ruling, may have a material adverse impact on our ability to conduct our business and our finances. Moreover, third parties making claims against us may be able to obtain injunctive relief against us, which could block our ability to offer one or more products or services and could result in a substantial award of damages against us. In addition, since we sometimes indemnify customers, collaborators or licensees, we may have additional liability in connection with any infringement or alleged infringement of third party intellectual property.

Because patent applications can take many years to issue, there may be pending applications, some of which are unknown to us, that may result in issued patents upon which our products or proprietary technologies may infringe. Moreover, we may fail to identify issued patents of relevance or incorrectly conclude that an issued patent is invalid or not infringed by our technology or any of our products. There is a substantial amount of litigation involving patent and other

intellectual property rights in our industry. If a third party claims that we or any of our licensors, customers or collaboration partners infringe upon a third party's intellectual property rights, we may have to:

- seek to obtain licenses that may not be available on commercially reasonable terms, if at all;
- abandon any infringing product or redesign our products or processes to avoid infringement;
- pay substantial damages, including, in exceptional cases, treble damages and attorneys' fees;
- pay substantial royalties or fees or grant cross-licenses to our technology;
- or defend litigation or administrative proceedings that may be costly whether we win or lose, and which could result in a substantial diversion of our financial and management resources.

We may be involved in lawsuits to protect or enforce our patents or the patents of our licensors, which could be expensive, time-consuming and unsuccessful.

Competitors may infringe our patents or the patents that we license. In the event of infringement or unauthorized use, we may file one or more infringement lawsuits. Patent litigation can be very costly and time-consuming, and the outcome is uncertain. In addition, if we or any of our partners were to initiate legal proceedings against a third party to enforce a patent covering one of our products or services, the defendant in such litigation could counterclaim that our patent is invalid and/or unenforceable. In patent litigation, defendant counterclaims alleging invalidity and/or unenforceability are commonplace. The outcome following legal assertions of invalidity and unenforceability is unpredictable. If a defendant were to prevail on a legal assertion of invalidity and/or unenforceability, we would lose at least part, and perhaps all, of the challenged patent. Such a loss of patent protection could have a material adverse impact on our business.

We may not be able to protect our intellectual property rights throughout the world, which could have a material adverse effect on our business.

Filing, prosecuting and defending patents on current and future products in all countries throughout the world would be prohibitively expensive, and our intellectual property rights in some countries outside of the United States can be less extensive than those in the United States. In addition, the laws of some foreign countries do not protect intellectual property rights to the same extent that federal and state laws do in the United States. Consequently, regardless of whether we are able to prevent third parties from practicing our inventions in the United States, we may not be able to prevent third parties from practicing our inventions in all countries outside of the United States, or from selling or importing products made by using our inventions in and into the United States or other jurisdictions. Competitors may use our technologies in jurisdictions where we have not pursued and obtained patent protection to develop their own products, and further, may export otherwise infringing products to territories where we have patent protection, but enforcement is not as strong as it is in the United States. These products may compete with our products and our patents or other intellectual property rights may not be effective or sufficient to prevent them from competing. Even if we pursue and obtain issued patents in particular jurisdictions, our patent claims or other intellectual property rights may not be effective or sufficient to prevent third parties from competing. Patent protection must ultimately be sought on a country-by-country basis, which is an expensive and time-consuming process with uncertain outcomes. Accordingly, we may choose not to seek patent protection in certain countries, and we will not have the benefit of patent protection in such countries.

Many companies have encountered significant problems in protecting and defending intellectual property rights in foreign jurisdictions. The legal systems of certain countries, such as China and certain developing countries, do not favor the enforcement of patents and other intellectual property protection, particularly those relating to biotechnology, which could make it difficult for us to stop the infringement of our patents or marketing of competing products in violation of our proprietary rights generally. Proceedings to enforce our patent rights in foreign jurisdictions could result in substantial costs and divert our efforts and attention from other aspects of our business, put our patents at risk of being invalidated or interpreted narrowly and our patent applications at risk of not issuing, and provoke third parties to assert claims against us. We may not prevail in any lawsuits that we initiate and the damages or other remedies awarded, if any, may not be commercially meaningful. Accordingly, our efforts to enforce our intellectual property rights around the world may be inadequate to obtain a significant commercial advantage from the intellectual property that we develop or license and may adversely impact our business.

In addition, we and our partners also face the risk that our products are imported or reimported into markets with relatively higher prices from markets with relatively lower prices, which would result in a decrease of sales and any payments we receive from the affected market. Recent developments in U.S. patent law have made it more difficult to stop these and related practices based on theories of patent infringement.

We use third-party software that may be difficult to replace or may cause errors or failures of our products that could lead to lost customers or harm to our reputation. Additionally, our use of “open source” software could adversely affect our ability to offer our products and technologies and subject us to possible litigation.

We use software licensed from third parties in our products. In the future, this software may not be available to us on commercially reasonable terms, or at all. Any loss of the right to use any of this software could result in delays in the production of our products until equivalent technology is either developed by us, or, if available, is identified, obtained and integrated, which could harm our business. In addition, any errors or defects in third-party software or other third-party software failures could result in errors, defects or cause our products to fail, which could harm our business and be costly to correct. Many of these providers attempt to impose limitations on their liability for such errors, defects or failures, and if enforceable, we may have additional liability to our customers or third-party providers that could harm our reputation and increase our operating costs.

Additionally, we use open source software in connection with our products and technologies. Companies that incorporate open source software into their technologies have, from time to time, faced claims challenging the use of open source software and/or compliance with open source license terms. As a result, we could be subject to suits by parties claiming ownership of what we believe to be open source software or claiming non-compliance with open source licensing terms. Some open source software licenses require users who distribute software containing open source software to publicly disclose all or part of the source code to such software and/or make available any derivative works of the open source code, which could include valuable proprietary code of the user, on unfavorable terms or at no cost. While we monitor our use of open source software and try to ensure that none is used in a manner that would require us to disclose our internally developed source code or that would otherwise breach the terms of an open source agreement, such use could inadvertently occur, in part because open source license terms are often ambiguous. In addition to risks related to license requirements, use of certain open source software can lead to greater risks than use of third-party commercial software, as open source licensors generally do not provide warranties or controls on the origin of software which, thus, may contain security vulnerabilities or infringing or broken code. Any requirement to publicly disclose our internally developed source code or pay damages for breach of contract could have a material adverse effect on our business.

Risks Related to our Common Stock and Being a Public Company

The market price of our common stock has fluctuated significantly and may continue to fluctuate significantly.

The market price of shares of our common stock has been and could continue to be subject to wide fluctuations in response to many factors listed in this section, and others beyond our control, including:

- actual or anticipated fluctuations in our financial condition and operating results;
- announcements by us, our partners or our competitors of new products, significant contracts, restructuring plans, strategic partnerships, joint ventures, collaborations, acquisitions, commercial relationships or capital commitments;
- competition from existing products or new products that may emerge;
- failure to meet or exceed financial estimates and projections of the investment community or that we may provide to the public;
- issuance of new or updated research or reports by securities analysts or recommendations with respect to our stock;
- positive or adverse regulatory announcements;
- disputes or other developments related to proprietary rights, including patents, litigation matters and our ability to obtain patent protection for our technologies; commencement of, or our involvement in, litigation;

- fluctuations in the valuation of companies perceived by investors to be comparable to us;
- conditions in our markets;
- manufacturing disputes or delays, product defects or material product quality control issues;
- any future sales of our common stock or other securities;
- any change to the composition of our board or key personnel;
- general economic conditions and slow or negative growth of our markets;
- a material cybersecurity incident;
- share price and volume fluctuations attributable to inconsistent trading volume levels of our shares;
- announcement or expectation of additional debt or equity financing efforts; and
- other factors described in this Risk Factors section of this Annual Report on Form 10-K or in our other reports filed with the SEC.

These and other market and industry factors may cause the market price and demand for our common stock to fluctuate substantially, regardless of our actual operating performance, which may limit or prevent investors from readily selling their shares of our common stock and may otherwise negatively affect the liquidity of our common stock. In addition, the stock market in general, and life science companies in particular, have experienced extreme price and volume fluctuations that have often been unrelated or disproportionate to the operating performance of these companies. In the past, when the market price of a stock has been volatile, holders of that stock have on occasion instituted securities class action litigation against the company that issued the stock. If any of our stockholders were to bring a lawsuit against us, the defense and disposition of the lawsuit could be costly and divert the time and attention of our management and harm our operating results.

We have never paid dividends on our capital stock, and we do not anticipate paying any dividends in the foreseeable future. Consequently, any gains from an investment in our common stock will likely depend on whether the price of our common stock increases.

We have not paid dividends on any of our classes of capital stock to date and we currently intend to retain our future earnings, if any, to fund the development and growth of our business. As a result, capital appreciation, if any, of our common stock will be the sole source of gain for the stockholders in the foreseeable future. Consequently, in the foreseeable future, stockholders will likely only experience a gain from an investment in our common stock if the price of our common stock increases.

Anti-takeover provisions contained in our restated certificate of incorporation and restated by-laws, as well as provisions of Delaware law, could impair a takeover attempt.

Our restated certificate of incorporation, restated by-laws and Delaware law contain provisions which could have the effect of rendering more difficult, delaying or preventing an acquisition deemed undesirable by our board. Our corporate governance documents provisions include: authorizing our board to issue up to 5,000,000 shares of preferred stock without stockholder approval upon the terms and conditions and with the rights, privileges and preferences as our board may determine; specifying that special meetings of our stockholders can be called only by our board and that our stockholders may not act by written consent; establishing an advance notice procedure for stockholder proposals to be brought before an annual meeting of our stockholders, including proposed nominations of persons for election to our board; providing that directors may be removed only for cause; providing that our board may create new directorships and that vacancies on the board may be filled only by a majority of directors then in office, even though less than a quorum; and providing that our board may amend our bylaws without approval of our stockholders. These provisions, alone or together, could delay or prevent hostile takeovers and changes in control or changes in our management. As a Delaware corporation, we are also subject to provisions of Delaware law, including Section 203 of the DGCL, which prevents certain stockholders holding 15% or more of our outstanding common stock from engaging in certain business combinations

without approval of the holders of substantially all of our outstanding common stock. Any provision of our certificate of incorporation, our bylaws or the DGCL that has the effect of delaying or deterring a change in control could limit the opportunity for our stockholders to receive a premium for their shares of our common stock, and could also affect the price that some investors are willing to pay for our common stock.

ITEM 1B. UNRESOLVED STAFF COMMENTS

Not applicable.

ITEM 1C. CYBERSECURITY

We maintain a cybersecurity risk management program designed to identify, assess, manage, mitigate and respond to cybersecurity threats. This program incorporates policies, processes, and activities over domains such as access control, facility and data protection, IT systems and data transmission security, threat intelligence and incident response, third-party risk management, disaster recovery, and vulnerability management. We have a comprehensive cybersecurity policy in place that includes robust measures to protect our data and systems from cyber threats, ensuring the integrity and confidentiality of our information.

We have implemented a risk-based approach to identify and assess the cybersecurity threats that could affect our business and information systems. We base our program on multiple security frameworks including the National Institute of Standards and Technology ("NIST"), HIPAA, and privacy laws such as the E.U.'s GDPR. We also require that third-party service providers with access to personal or proprietary information implement and maintain cybersecurity practices consistent with applicable legal standards.

Cybersecurity is complex and subject to constantly evolving threats. Accordingly, we engage a range of external experts, including cybersecurity consultants and auditors in evaluating and testing our risk management systems. These partnerships enable us to leverage specialized knowledge and insights. Our collaboration with these third-party experts includes regular audits, threat assessments, and consultation on security enhancements.

The head of our IT department is tasked with ensuring that the highest levels of management and our board of directors are informed about the cybersecurity posture and potential risks facing the Company. Our IT team regularly briefs our CEO about cybersecurity risk management. The IT staff regularly informs the Head of IT about the latest developments in cybersecurity, including potential threats and risk management techniques. In the event of a cybersecurity incident, the Head of IT is informed promptly following its detection, and our response is governed by a detailed incident response plan that includes prompt actions to mitigate the impact of the incident and longer-term strategies for remediation and prevention of future incidents. The Head of IT has also chartered an Information Security Steering Committee made up of cross-functional executive leaders that meets quarterly on topics such as the current cybersecurity landscape and emerging threats, status of ongoing cybersecurity initiatives and strategies, incident reports and learnings from any cybersecurity events, and compliance with regulatory requirements and industry standards.

Cybersecurity risk management is integrated into our broader risk management framework. Our audit committee of our board of directors, which has responsibility for oversight of risk management, also has responsibility for oversight of our program, policies and procedures related to information security and data protection. On a regular basis, the Audit Committee of our Board of Directors receives reports on cybersecurity risks as well as mitigation strategies and the status of initiatives to strengthen our information security systems. We also mandate a rigorous cybersecurity training program for all employees, which includes regular updates and assessments to ensure they are equipped to recognize and respond to potential cyber threats.

For a discussion of our risks related to cybersecurity, see the section titled, "Part 1. Item 1A. Risk Factors – *Risks Related to our Operations – Cybersecurity breaches, loss of data and other disruptions could compromise sensitive information related to our business or prevent us from accessing critical information and expose us to liability, which could adversely affect our business and our reputation.*" included in this Annual Report on Form 10-K.

ITEM 2. PROPERTIES

Our corporate headquarters is currently located in Billerica, Massachusetts, which consists of an approximately 91,600 square foot facility we lease for office and manufacturing purposes. We also lease office and laboratory space domestically in Bedford and Marlborough, Massachusetts and Georgetown, Texas, and internationally in the Netherlands, Sweden, and China. We do not currently occupy one of our Bedford facilities or two of the facilities acquired from Akoya (in Menlo Park, California and Marlborough, Massachusetts) and currently sublease a portion of these facilities.

We believe our facilities are adequate and suitable for our current operations and that, should it be needed, additional or alternative space is available to accommodate our operations.

ITEM 3. LEGAL PROCEEDINGS

In the ordinary course of business, we are from time to time involved in lawsuits, claims, investigations, proceedings, and threats of litigation consisting of intellectual property, contractual, employment, and other matters. While the outcome of any such actions or proceedings cannot be predicted with certainty, as of the date of this Annual Report on Form 10-K, we were not party to any legal proceedings, the outcome of which would be expected to have a material adverse effect on our financial condition or results of operations. Regardless of any outcome, litigation can have a material adverse effect on us due to defense and settlement costs, diversion of management resources, and other factors.

ITEM 4. MINE SAFETY DISCLOSURES

Not applicable.

PART II

ITEM 5. MARKET FOR REGISTRANT'S COMMON EQUITY, RELATED STOCKHOLDER MATTERS, AND ISSUER PURCHASES OF EQUITY SECURITIES

Market Information

Our common stock is traded on The Nasdaq Global Market under the symbol "QTRX."

Holders of Record

As of February 24, 2026, there were approximately 68 holders of record of our common stock.

Dividends

The declaration of dividends is at the discretion of our Board of Directors and depend on actual cash flow from operations, our financial condition, capital requirements, and any other factors our Board of Directors may consider relevant. Since our inception, we have not declared or paid any dividends.

Securities Authorized for Issuance under Equity Compensation Plans

Refer to the section titled "Part III, Item 12. Security Ownership of Certain Beneficial Owners and Management and Related Stockholder Matters" of this Annual Report on Form 10-K for information required by Item 201(d) of Regulation S-K.

Unregistered Sales of Securities

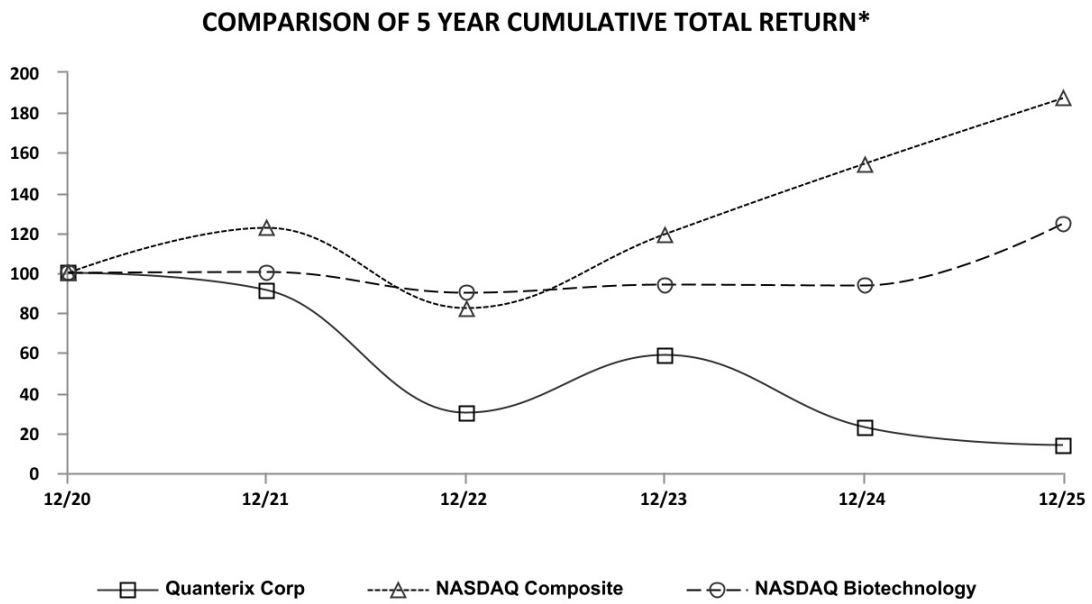
There were no unregistered sales of equity securities during the year ended December 31, 2025.

Issuer Purchases of Equity Securities

Not applicable.

Stock Performance Graph

The following graph compares the cumulative total shareholder returns over the past five years for our common stock, the NASDAQ Composite Index, and the NASDAQ Biotechnology Index, assuming \$100 invested on December 31, 2020, and reinvestment of dividends, if paid:



*\$100 invested on 12/31/20 in stock or index, including reinvestment of dividends.
Fiscal year ending December 31.

ITEM 6. RESERVED

Not applicable.

ITEM 7. MANAGEMENT’S DISCUSSION AND ANALYSIS OF FINANCIAL CONDITION AND RESULTS OF OPERATIONS

The following discussion and analysis provides information management believes to be relevant to understanding the financial condition and results of operations of Quanterix Corporation for the years ended December 31, 2025 and 2024. For a full understanding of our financial condition and results of operations, this discussion and analysis should be read in conjunction with our Consolidated Financial Statements and accompanying notes included in the section titled “Part II. Item 8. Financial Statements and Supplementary Data” of this Annual Report on Form 10-K. Certain columns and rows may not add due to the use of rounded numbers. Percentages presented are calculated from the underlying unrounded numbers. In addition to historical information, the following discussion and analysis contains forward-looking statements that involve risks, uncertainties, and assumptions. Our actual results, performance, or experience may differ materially from those discussed below due to various important factors, risks, and uncertainties, including, but not limited to, those set forth under the sections titled “Part I, Item 1A. Risk Factors” and “Note Regarding Forward-Looking Statements” included in this Annual Report on Form 10-K. Unless the context otherwise requires, the terms “Quanterix,” the “Company,” “we,” “it,” “us,” “and “our” in this Annual Report on Form 10-K refer to Quanterix Corporation and its consolidated subsidiaries.

The discussion and analysis of our financial condition as of, and results of operations for the year ended December 31, 2024 as compared to the year ended December 31, 2023 is included in the section titled “Part II, Item 7. Management’s Discussion and Analysis of Financial Condition and Results of Operations” in our Annual Report on [Form 10-K for the year ended December 31, 2024](#), as filed with the U.S Securities and Exchange Commission (the “SEC”) on March 17, 2025.

Overview

We are a life sciences company transforming healthcare innovation by accelerating biomarker breakthroughs from discovery to diagnostics using our ultra-sensitive translational research and spatial biology instruments, consumables, and services. We continue to invest in pushing a paradigm shift in healthcare from an emphasis on later-stage treatment to a focus on earlier detection, monitoring, prognosis, and, ultimately, prevention. Our combined platforms have achieved significant commercial adoption with an installed base of over 2,500 instruments and scientific validation with citations in more than 6,200 scientific publications in areas of high unmet medical need and research interest such as neurology, oncology, immunology, and inflammation.

Our proprietary digital “Simoa” detection technology enables customers to reliably detect protein biomarkers at ultra-low concentrations in blood, serum and other fluids that, in many cases, are undetectable using conventional, analog immunoassay technologies. Multi-plexing biomarker analysis in tissue samples with our “Spatial Biology” platforms enables scientists to understand the localized interactions occurring on the cellular level. We believe our combination of technologies will enable scientists to help drive diagnostic innovation in the evolving healthcare landscape with data across the tissue to fluid continuum. Currently, the ability of our Simoa platforms to detect proteins in the femtomolar range is enabling the development of novel therapies and diagnostics and has the potential to identify early-stage disease markers before symptoms appear.

Our instruments are designed to be used either with assays fully developed by us, including all antibodies and supplies required to run the assays, or with “homebrew” assay kits where we supply some of the components required for testing, and the customer supplies the remaining required elements. Accordingly, our installed instruments generate a recurring revenue stream. As the installed base of our instruments increases, we expect total consumables revenue to increase.

We also provide contract research services and clinical laboratory testing services, including four Laboratory Developed Tests (“LDT”), using our proprietary Simoa and Spatial Biology technology through our Accelerator Laboratory which is certified under the Clinical Laboratory Improvement Amendments of 1988 (“CLIA”) (the “Accelerator Laboratory”). To date, we have completed over 2,600 projects for more than 500 customers from all over the world using our platforms.

We have an extensive base of worldwide customers including research laboratories, contract research organizations ("CROs"), academic institutions, and bio-pharmaceutical companies. We sell our instruments, consumables, and services through a direct field sales and support organizations in North America and Europe, and through our own sales force and distributors in additional countries, including Australia, Brazil, China, Czech Republic, India, Hong Kong, Israel, Japan, New Zealand, Qatar, Saudi Arabia, Singapore, South Africa, South Korea, Taiwan, and the United Arab Emirates.

Our total revenues were \$138.9 million and \$137.4 million for the years ended December 31, 2025 and 2024, respectively. Since our inception, we have incurred annual net losses, including net losses of \$107.2 million and \$38.5 million for the years ended December 31, 2025 and 2024, respectively.

We expect to incur operating losses into 2026 as we incur costs related to:

- expanding our research and development efforts to improve our existing, or to develop and launch, new assays and instruments. These expenses could be particularly significant if any of our products become subject to additional or more burdensome regulation by the U.S Food and Drug Administration (the "FDA");
- investing in Lucent Diagnostics, additional laboratory developed tests ("LDTs"), and other diagnostics initiatives including entry into translational pharma and clinical diagnostic markets;
- seeking Premarket Approval ("PMA"), de novo classification, or 510(k) clearance from the FDA for our existing or new products, including new assays and instruments, if or when we decide to market products for use in the prevention, diagnosis, or treatment of a disease or other condition;
- strategically acquiring and integrating companies or technologies that may be complementary to our business;
- making required earnout payments under the Emission, Inc. ("Emission") acquisition agreement, which are contingent upon the achievement of certain performance milestones;
- entering into collaboration arrangements, or in-license other products and technologies; and
- adding or enhancing operational, financial, and management information systems.

As further described in the section titled "Recent Business Developments - Restructuring Costs", in 2025 we implemented actions to realize many of the synergies of the acquisition of Akoya Biosciences, Inc. ("Akoya"). As a result of these actions and our continued integration activities, we expect to be cash flow breakeven in 2026, although our ability to achieve this goal is dependent on our success in meeting both revenue and expense objectives.

Recent Business Developments

Acquisitions

Akoya Biosciences, Inc.

On July 8, 2025, we acquired Akoya, a life sciences technology company based in Marlborough, Massachusetts delivering spatial biology solutions through the power of spatial phenotyping. Spatial phenotyping refers to a rapidly evolving technology that enables academic and biopharma scientists to detect and map the distribution of cell types and biomarkers across whole tissue samples at single-cell resolution, enabling advancements in their understanding of disease progression and patient response to therapy. Akoya commercializes proprietary instrument platforms, reagents, software, and services that offer end-to-end solutions to perform tissue analysis and spatial phenotyping from discovery through translational and clinical research and diagnostics.

Pursuant to the acquisition agreement, the consideration transferred was \$151.0 million, which consisted of Quanterix common stock with a fair value of \$49.9 million, \$18.9 million of cash to Akoya shareholders, and \$82.1 million of cash to settle Akoya's debt financing arrangement with Midcap Financial Trust (the "Midcap Trust Term Loan"). The debt payment included a \$75.1 million payoff of the principal outstanding and \$7.0 million in early termination, legal, and prepayment fees.

Emission, Inc.

On January 8, 2025, we acquired Emission, a life sciences manufacturing company based in Georgetown, Texas. Emission produces large-scale, highly-uniform dye-encapsulating magnetic beads designed for low and mid-plex assays and a mid-plex platform that reads its proprietary beads. The transaction is part of our plans to secure the use of Emission's highly controlled beads in our next generation platforms and expansion into a new multi-plex segment targeting third-party original equipment manufacturer customers. As part of the acquisition of Emission, we made an upfront payment of \$9.0 million, with up to an additional \$1.0 million payable at the end of the holdback period, and a \$10.0 million payment in the fourth quarter of 2025 upon completion of certain technical milestones. Additionally, the selling shareholders of Emission (collectively, the "Emission Shareholders") may receive up to an additional \$50.0 million in earnout payments through December 31, 2029, contingent upon the achievement of certain performance milestones.

In connection with the closing of the acquisition, the parties entered into a call option agreement (the "Option Agreement"), in which the Emission Shareholders have the right to repurchase all of the outstanding capital stock of Emission for \$10.0 million after five years if Emission's revenues do not exceed \$5.0 million in any one year during such five-year period. If the Emission Shareholders exercise the right to repurchase Emission under the Option Agreement and consummate the repurchase, we will retain a perpetual, fully-paid, irrevocable license to all Emission intellectual property required to continue to manufacture and commercialize our products.

For further information about the Akoya and Emission acquisitions, refer to Note 3 - *Acquisitions* in the Notes to Consolidated Financial Statements included in this Annual Report on Form 10-K.

FDA 510(k) Submission for a Multi-Analyte Algorithmic Blood Test for Alzheimer's Disease Detection

On January 31, 2026, we submitted a 510(k) premarket notification to the U.S. Food and Drug Administration ("FDA") for a multi-analyte algorithmic blood test for Alzheimer's disease.

This submission represents a significant milestone in the Company's mission to provide superior, non-invasive, high-performance diagnostic tools to aid in the evaluation of patients with cognitive symptoms for possible Alzheimer's disease. The multi-analyte test previously received Breakthrough Device Designation from the FDA, a program intended to accelerate the development and review of devices that provide for more effective treatment or diagnosis of life-threatening or irreversibly debilitating diseases. The test is intended to aid in identifying whether patients with cognitive symptoms are likely to have amyloid brain plaques—a hallmark of Alzheimer's—providing diagnostic clarity through a non-invasive blood test.

LucentAD Complete Medicare Pricing

In November 2025, the Centers for Medicare & Medicaid Services ("CMS") approved a reimbursement rate of \$897 for our LucentAD Complete multiplex test. This milestone provides a nationally recognized reference price, an important step for coverage decision with private issuers, enables broader access and supports efforts to bring this multiplex diagnostic solution to hospitals and laboratories across the country. In addition, based on available data, we believe LucentAD Complete has the potential to obtain a Local Coverage Determination, which would support more consistent reimbursement of claims.

Goodwill Impairment

During 2025, we assessed several events and circumstances, including a larger than expected decline in the Company's revenue and bookings primarily due to the rapidly changing macro-economic conditions resulting from reductions in U.S. federal research funding, reductions in research and development spending by larger pharmaceutical customers, and new import tariffs. Based on an impairment test of our goodwill as of June 30, 2025, which estimated the implied fair value of our then single reporting unit under a market valuation approach (using inputs such as our quoted stock price), we determined our entire goodwill balance was impaired. As a result, we recorded a \$6.4 million impairment charge in the second quarter of 2025.

Restructuring Costs

In 2025, we announced actions to reduce operating costs, preserve cash, and realize anticipated synergies and other benefits of the Akoya acquisition, including reductions in our workforce. The actions we implemented in 2025 reduced our annualized operating expenses by approximately \$74.0 million. As we continue to integrate Akoya, we expect to take additional steps into 2026 to further reduce our operating expenses, with a goal of reducing our annualized operating expenses by a total of approximately \$85.0 million.

We incurred expenses of \$8.0 million related to the reductions in force in 2025, substantially all of which relate to cash expenditures for severance and related benefits. These reductions in force were completed in 2025.

Components of Results of Operations

Revenues

During 2025, we experienced a larger than expected decline in our revenue and bookings, primarily due to rapidly changing macro-economic conditions. These macro-economic conditions, including reductions in U.S. federal research funding and new tariffs, impacted demand from academic customers and caused decreased demand from pharmaceutical customers as research and development spending and clinical trials slowed down. These conditions caused decreased revenue for our products and Accelerator Laboratory services.

We continue to see strong opportunities with our customers and expect that revenue and bookings in 2026 will be consistent with 2025.

Product Revenue

Our product revenues are generated from sales of (1) instruments, including the related installation, and (2) consumables and related revenues. Our products are sold directly to customers and are also sold through distributors in EMEA and Asia Pacific regions.

Instrument revenues consist of sales of our instruments. We currently sell our products for research use only ("RUO") applications directly to customers or through distributors. Customers' purchase processes for certain of our instruments can be long and as a result, instrument revenue can vary from period-to-period and can be concentrated to a small number of customers in any given period. Instruments sold directly to customers include an initial year service-type warranty, which is recorded in services and other revenue on the Consolidated Statements of Operations. Instruments sold to distributors include a license to import and resell the instruments and an initial year assurance-type warranty. Costs related to assurance-type warranties are recorded in cost of product revenue on the Consolidated Statements of Operations. Instrument sales may also be bundled with assays and other consumables, training, installation, and/or an extended service warranty.

Consumable and other revenues consist of sales of assays fully developed by us, including all antibodies and supplies required to run the assays, or with "homebrew" assay kits where we supply some of the components required for testing, and the customer supplies the remaining required elements. Consumable and other revenues also consist of replacement parts, reagents, and antibodies.

Service and Other Revenue

Service revenues generally consist of fixed fee contract research services through our Accelerator Laboratory, initial service-type warranties, extended service warranty contracts, repair services, and other services such as training.

Collaboration and License Revenue

Collaboration and license revenues consist of licensing our technology, intellectual property, and know-how associated with our instruments to third parties and for related services. License arrangements consist of sales or usage-based fees and/or future royalties.

Cost of Goods Sold and Services

Cost of Product Revenue

Cost of product revenue consists of manufacturing and assembly costs for instruments, related reagents, other consumables, contract manufacturer costs, personnel costs, overhead, and other direct costs related to product sales. Raw material part costs include inbound shipping and handling costs associated with purchased goods. Cost of product revenue also includes amortization expense for acquired intangibles that relate to the sale of products as well as royalty fees due to third parties from revenue generated by collaboration or license deals.

Cost of Service and Other Revenue

Cost of services and other revenue consists of direct costs associated with operating our Accelerator Laboratory on behalf of customers, including raw materials, personnel costs, royalties, allocated overhead and other related costs. Additional costs include costs related to warranty services and other costs of servicing equipment at customer sites.

Research and Development Expense

Research and development expense consists of personnel costs, research supplies, third-party development costs for new products, materials for prototypes, quality assurance, and allocated overhead costs that include facility and other related costs. Our research and development efforts have focused primarily on supporting development and commercialization of new and existing products and improved product quality. We believe that our continued investment in research and development is essential to our long-term competitive position. We have made substantial investments in research and development since our inception and we expect to continue to drive innovation through investments in future product development.

We believe that our continued investment in research and development is essential to our long-term competitive position. We expect that the realization of anticipated synergies from the acquisition of Akoya should enable us to maintain research and development expense at a more consistent level period to period in the future.

Selling, General and Administrative Expense

Selling, general and administrative expense consists of personnel costs for our sales and marketing, finance, legal, human resources, and general management teams, shipping and handling for product sales, acquisition related costs, other general and administrative costs, as well as professional services costs, such as marketing, advertising, legal and accounting services, and allocated overhead costs that include facility and other related costs.

We expect to increase the size of our selling, general and administrative functions to support the growth in our business. However, we intend to manage the rate of increase in selling, general and administrative expenses in the future so that it remains below any future rate of increase in revenues.

The classification of shipping and handling costs for product sales varies from company to company, with some companies recording these as selling, general and administrative expenses and others recording such expenses within costs of goods sold for products. To the extent our classification of these shipping and handling costs differs from the classification used by other companies, our gross margins may not be comparable with those reported by such other companies.

Other Lease Costs

Other lease costs consist of the amortization of operating lease right-of-use assets and other facility operating expenses from leased facilities we are not using as a result of restructurings.

Impairment and Restructuring Costs

Impairment and restructuring costs primarily consists of charges recorded as a result of 2025 restructuring actions and the corresponding impairment of a portion of our goodwill and long-lived assets (including operating lease right-of-use assets, property and equipment), which were determined to have carrying values exceeding their fair values.

Comparison of Results of Operations for Years Ended December 31, 2025 and 2024:

The following table sets forth select Consolidated Statements of Operations data, and such data as a percentage of total revenues (in thousands, except percentages):

	Year Ended December 31,				Increase (Decrease)	
	2025	% of revenue	2024	% of revenue	Amount	%
Revenues:						
Product revenue	\$ 92,941	67 %	\$ 79,740	58 %	\$ 13,201	17 %
Service and other revenue	44,212	32 %	51,244	37 %	(7,032)	(14)%
Collaboration and license revenue	1,501	1 %	4,452	3 %	(2,951)	(66)%
Grant revenue	243	— %	1,985	1 %	(1,742)	(88)%
Total revenues	138,897	100 %	137,421	100 %	1,476	1 %
Costs of goods sold and services:						
Cost of product revenue	50,981	37 %	33,304	24 %	17,677	53 %
Cost of service and other revenue	22,957	16 %	21,013	16 %	1,944	9 %
Total costs of goods sold and services	73,938	53 %	54,317	40 %	19,621	36 %
Gross profit	64,959	47 %	83,104	60 %	(18,145)	(22)%
Operating expenses:						
Research and development	35,922	26 %	31,082	23 %	4,840	16 %
Selling, general and administrative	138,008	99 %	101,618	74 %	36,390	36 %
Other lease costs	844	1 %	3,020	2 %	(2,176)	(72)%
Impairment and restructuring	15,727	11%	—	—%	15,727	100 %
Total operating expenses	190,501	137 %	135,720	99 %	54,781	40 %
Loss from operations	(125,542)	(90)%	(52,616)	(39)%	(72,926)	139 %
Interest income	8,567	6 %	14,655	11 %	(6,088)	(42)%
Change in fair value of contingent liabilities	4,547	3 %	—	— %	4,547	100 %
Other income (expense), net	157	— %	(136)	— %	293	(215)%
Loss before income taxes	(112,271)	(81)%	(38,097)	(28)%	(74,174)	195 %
Income tax benefit (expense)	5,121	4 %	(434)	— %	5,555	(1280)%
Net loss	\$ (107,150)	(77)%	\$ (38,531)	(28)%	\$ (68,619)	178 %

Revenues

Total revenues increased \$1.5 million, or 1%, to \$138.9 million for the year ended December 31, 2025, compared to \$137.4 million for the year ended December 31, 2024.

For the year ended December 31, 2025, product revenue of \$92.9 million consisted of instrument sales of \$17.9 million and sales of consumables and other products of \$75.1 million. Product revenue increased \$13.2 million, or 17%, compared to \$79.7 million for the year ended December 31, 2024. The increase in product revenue was primarily due to the acquisition of Akoya, which added \$21.9 million of product revenue. Product revenue for the legacy Quanterix business decreased \$8.7 million, or 11%, primarily due to decreased demand for consumables as a result of the changing macro-economic conditions. Legacy Quanterix instrument sales in 2025 were consistent with 2024.

Service and other revenue decreased \$7.0 million, or 14%, to \$44.2 million for the year ended December 31, 2025, compared to \$51.2 million for the year ended December 31, 2024. This decrease was primarily due to a decrease of \$18.8 million, or 37%, in the legacy Quanterix business due to the macro-economic conditions impacting demand for large projects from pharmaceutical customers and the completion of an agreement with Eli Lilly and Company which generated \$4.5 million of revenue in 2024. The decrease was partially offset by the acquisition of Akoya, which added \$11.8 million of service and other revenue. Akoya's service and other revenue includes \$2.7 million of non-cash revenue from the off-

market component of an acquired contract and represents a portion of the purchase accounting fair value adjustment that would have been recognized by a market participant.

Collaboration and license revenue decreased \$3.0 million, or 66%, to \$1.5 million for the year ended December 31, 2025, compared to \$4.5 million for the year ended December 31, 2024. The decrease was primarily due to LDT and other diagnostic related license revenues in 2024 that did not repeat in 2025.

Grant revenue decreased \$1.7 million, or 88%, to \$0.2 million for the year ended December 31, 2025, compared to \$2.0 million for the year ended December 31, 2024. The decrease was primarily due to completion of milestones in 2024 under certain grants.

Cost of Goods Sold and Services

Total cost of goods sold and services increased \$19.6 million, or 36%, to \$73.9 million for the year ended December 31, 2025 compared to \$54.3 million for the year ended December 31, 2024.

Cost of product revenue increased \$17.7 million, or 53%, to \$51.0 million for the year ended December 31, 2025, compared to \$33.3 million for the year ended December 31, 2024. The increase was primarily due to the acquisition of Akoya, which added \$14.1 million to cost of product revenue. Cost of product revenue for the legacy Quanterix business increased \$3.9 million primarily due to lower cost absorption as a result of lower output and increased costs due to the introduction of new assays. These increases were partially offset by decreases in headcount and related compensation and benefit costs from the restructuring implemented in 2025.

Cost of service and other revenue increased \$1.9 million, or 9%, to \$23.0 million for the year ended December 31, 2025, compared to \$21.0 million for the year ended December 31, 2024. This increase was primarily due to the acquisition of Akoya, which added \$7.6 million to cost of service and other revenue. Cost of service revenue for the legacy Quanterix business decreased \$5.6 million primarily due to decreases in headcount and related compensation and benefit costs from the restructuring implemented in 2025 and decreased demand for Accelerator Laboratory services.

Research and Development

Research and development expense increased \$4.8 million, or 16%, to \$35.9 million for the year ended December 31, 2025, compared to \$31.1 million for the year ended December 31, 2024. This increase was primarily due the acquisition of Akoya, which added \$3.2 million of research and development expenses. Research and development expense for the legacy Quanterix business increased \$1.4 million primarily due to a \$5.1 million charge associated with the contingent compensation payable pursuant to the terms of the Emission acquisition. This increase was partially offset by a \$2.9 million decrease in costs of outside services, research lab supplies, and equipment related to product development and \$0.8 million due to decreases in headcount and related compensation and benefit costs from the restructuring implemented in 2025.

Selling, General and Administrative

Selling, general and administrative expense increased \$36.4 million, or 36% to \$138.0 million for the year ended December 31, 2025, compared to \$101.6 million for the year ended December 31, 2024. Included within selling, general and administrative expense are \$5.6 million and \$8.1 million of shipping and handling costs for product sales for the years ended December 31, 2025 and 2024, respectively.

The increase was partially due to the acquisition of Akoya, which added \$14.8 million of selling, general and administrative expenses. Selling, general and administrative expense for the legacy Quanterix business increased \$21.9 million primarily due to (1) \$16.4 million of non-recurring acquisition and integration costs related to the acquisition of Akoya, (2) a \$4.9 million charge associated with the contingent compensation payable pursuant to the terms of the Emission acquisition, (3) a \$1.6 million increase in software and information technology expenses, and (4) a \$1.3 million increase related to a leased facility we began using in the fourth quarter of 2024. These increases were partially offset by a \$2.5 million decrease in shipping and handling costs primarily due to lower sales and a \$1.5 million decrease in audit fees including the remediation of the prior year material weaknesses in our internal control over financial reporting.

Other Lease Costs

Other lease costs decreased \$2.2 million, or 72%, to \$0.8 million for the year ended December 31, 2025, compared to \$3.0 million for the year ended December 31, 2024. In the fourth quarter of 2024, we began using one of the leased facilities that we did not occupy as a result of a restructuring in 2022. Accordingly, as of the fourth quarter of 2024, the amortization of the operating lease right-of-use asset and related leased facility operating expenses at this facility are no longer recorded in other lease costs.

Impairment and Restructuring

We recorded impairment and restructuring costs of \$15.7 million for the year ended December 31, 2025 relating to a goodwill impairment charge, severance and related benefit expenses from the 2025 restructuring actions, and impairment of leased facilities acquired in the Akoya acquisition that we are not using. No such costs were recorded in the year ended December 31, 2024.

Income Tax Benefit (Expense)

Income tax benefit was \$5.1 million for the year ended December 31, 2025 compared to income tax expense of \$0.4 million for the year ended December 31, 2024. The change was primarily due to the release of a portion of our valuation allowance on deferred tax assets due to temporary tax differences related to the acquisitions of Emission and Akoya.

Liquidity and Capital Resources

Our principal sources of liquidity are cash, cash equivalents, marketable securities, and funds generated from sales of our products and services. As of December 31, 2025, we had \$29.8 million of cash and cash equivalents and \$88.4 million of marketable securities. Historically, we have also financed our operations through equity offerings and borrowings from credit facilities. Our liquidity requirements have consisted, and we expect that they will continue to consist, of sales and marketing expenses, research and development expenses, working capital, general corporate expenses, and contingent payments related to our acquisition activity.

We believe our cash, cash equivalents, and marketable securities, along with funds generated from sales of our products and services, will be sufficient to meet our anticipated operating cash requirements for at least 12 months from the date of this Annual Report on Form 10-K.

As a result of the acquisition of Akoya, along with actions already taken to reduce operating costs, preserve cash, and realize anticipated synergies and other benefits of acquisition, we expect to be cash flow breakeven in 2026, although our ability to achieve this goal is dependent on our success in meeting both revenue and expense objectives.

Our future capital requirements will depend on many factors, including, but not limited to, our pace of growth, expansion, or introduction of new instruments, assays, and services, including Lucent Diagnostics and advancing access to our diagnostic tests, market acceptance of our products and services, regulatory requirements, regulatory approval of our products or services, and the effects of competition, technological developments, and broader market and economic trends.

We regularly assess other potential acquisitions and may need capital to pursue acquisitions of complementary businesses, services, and technologies. To the extent our existing cash, cash equivalents, and marketable securities are insufficient to fund future activities or requirements to continue operating our business, we may need to raise additional capital. If the conditions for raising capital are favorable, we may seek to finance future cash needs through public or private equity, debt offerings, or other financings.

If needed, we cannot guarantee that we will be able to obtain additional funds on acceptable terms, or at all. If we raise additional funds by issuing equity or equity-linked securities, our stockholders may experience dilution. Future debt financing, if available, may involve covenants restricting our operations or our ability to incur additional debt. Any debt or equity financing that we raise may contain terms that are not favorable to us or our stockholders. If we raise additional funds through collaboration and licensing arrangements with third parties, it may be necessary to relinquish some rights to our technologies or our products, or grant licenses on terms that are not favorable to us. If we do not have or are not able to obtain sufficient funds, if needed, we may have to delay development or commercialization of our products and services. We also may have to reduce marketing, customer support, or other resources devoted to our products, or cease operations.

Cash Flows

The following table summarizes our cash flows (in thousands):

	Year Ended December 31,	
	2025	2024
Net cash used in operating activities	\$ (77,236)	\$ (35,164)
Net cash provided by (used in) investing activities	50,231	(82,265)
Net cash provided by (used in) financing activities	(708)	456
Net decrease in cash, cash equivalents, and restricted cash	\$ (27,713)	\$ (116,973)

Operating Activities

We derive cash flows from operations primarily from the sale of our products and services. Our cash flows from operating activities are also significantly influenced by our use of cash for operating expenses to develop new products and services, invest in process and product improvements, and increase our sales and marketing efforts. We have historically experienced negative cash flows from operating activities as we have developed our technology, expanded our business, and built our infrastructure, however we expect to be cash flow breakeven in 2026, although our ability to achieve this goal is dependent on our success in meeting revenue and expense objectives.

Net cash used in operating activities was \$77.2 million and \$35.2 million for the years ended December 31, 2025 and 2024, respectively. The \$42.1 million increase in net cash used in operating activities was primarily driven by an overall increase in our net loss, partially due to the inclusion of Akoya's operating results, adjusted for non-cash items. These non-cash items primarily related to stock-based compensation expense, depreciation and amortization, and impairment charges. The increase was partially offset by changes in working capital items, primarily a decrease in accounts receivable from improved collections and a decrease in inventory from lower raw materials purchases. Cash used in operations also included payments for professional fees supporting due diligence, legal, and accounting activities related to the acquisition of Akoya.

Investing Activities

Our primary investing activities consist of purchases of marketable securities. Additionally, we use funds to acquire companies and to make capital expenditures for the purchase of equipment to support our infrastructure.

Net cash provided by investing activities was \$50.2 million during the year ended December 31, 2025, which reflected \$93.2 million used for the acquisitions of Akoya and Emission, \$69.8 million of purchases of marketable securities, \$2.6 million of purchases of property and equipment, offset by \$215.8 million of proceeds from the maturities of marketable securities.

Net cash used in investing activities was \$82.3 million during the year ended December 31, 2024, which reflected \$295.6 million of purchases of marketable securities, \$3.4 million of purchases of property and equipment and \$216.7 million of proceeds from the maturities of marketable securities.

Financing Activities

Net cash used in financing activities was \$0.7 million and net cash provided by financing activities was \$0.5 million during the year ended December 31, 2025 and 2024, respectively, primarily from sales of our common stock under our employee stock purchase plan and from the exercise of options under our equity incentive plan.

Future Cash Obligations

In addition to the future cash obligations described below, we have other payables and liabilities that may be legally enforceable but are not considered contractual commitments. Refer to Note 16 - *Commitments and Contingencies* in the Notes to Consolidated Financial Statements included in this Annual Report on Form 10-K for a summary of our purchase commitments and other obligations as of December 31, 2025.

Emission Acquisition

The acquisition of Emission includes an arrangement where up to \$50.0 million of cash could be payable to the selling shareholders based on the amount and timing of certain performance targets over a five-year period ending December 31, 2029 (the "Emission Earnout"). The Emission Earnout is recorded at fair value each reporting period and at December 31, 2025 was fair valued at \$2.0 million.

PKI/Revvity License Agreement

Through the acquisition of Akoya, the Company assumed a license and contingent payment obligation ("PKI License") from Akoya's 2018 purchase of the QPS division of PerkinElmer, Inc., subsequently known as Revvity, Inc. Under the PKI License, we are required to pay single digit royalties based on net sales of certain products and services. Amounts due under the PKI License are payable annually through March 2033 and there is no limit on the amount of consideration that could be owed. Akoya's payments under the PKI License have historically been in the range of approximately \$1.0 million to \$2.0 million per year. In accordance with ASC 805 - *Business Combinations*, we measure and recognize the PKI License as a liability at fair value. At December 31, 2025, the fair value was \$3.7 million.

Operating Leases

We lease office, laboratory, and manufacturing space for our employees and operations, as well as office equipment, under non-cancellable operating lease agreements (refer to Note 15 - *Leases* in the Notes to Consolidated Financial Statements). The remaining duration of non-cancellable operating leases ranges from four months to seven years. Remaining lease payments within one year, within two to three years, within four to five years, and greater than five years from December 31, 2025 are \$10.2 million, \$17.3 million, \$15.6 million, and \$0.7 million, respectively.

Critical Accounting Policies and Estimates

Our Consolidated Financial Statements and the related notes included elsewhere in this Annual Report on Form 10-K are prepared in accordance with accounting principles generally accepted in the United States ("U.S. GAAP"). The preparation of these Consolidated Financial Statements requires us to make estimates and assumptions that affect the reported amounts of assets, liabilities, revenue, expenses, and related disclosures. We base our estimates on historical experience, worldwide economic conditions, both general and specific to the life sciences industry, and on various other assumptions we believe to be reasonable under the circumstances. We evaluate our estimates and assumptions on an ongoing basis, and changes in accounting estimates may occur from period to period. Accordingly, actual results could differ significantly from the estimates. To the extent that there are material differences between these estimates and actual results, our future financial statement presentation, financial condition, results of operations, and cash flows will be affected.

Our significant accounting policies are described in Note 2 - *Significant Accounting Policies* in the Notes to Consolidated Financial Statements included in this Annual Report on Form 10-K. We believe that the assumptions and estimates in the following critical accounting policies involve a greater degree of judgment and complexity and accordingly are the most critical to understanding and evaluating the potential impact to our Consolidated Financial Statements.

Revenue from Contracts with Customers

We generate revenue from the sale of products, services, and licenses, as further described in the section titled "Components of Results of Operations".

For contracts with customers, we recognize revenue when a customer obtains control of promised products or services, for an amount that reflects the consideration expected to be received in exchange for those products or services. We follow the five-step framework prescribed by Financial Accounting Standards Board ("FASB") Accounting Standards Codification ("ASC") Topic 606 - *Revenue from Contracts with Customers* ("ASC 606") to determine revenue recognition: (i) identify the contract(s) with a customer; (ii) identify the performance obligation(s) in the contract; (iii) determine the transaction price, including variable consideration, if any; (iv) allocate the transaction price to the performance obligations in the contract; and (v) recognize revenue when (or as) the entity satisfies a performance obligation. Revenues are presented net of any sales, value added, or similar taxes collected from customers and remitted to the government.

We determine the transaction price based on the amount of consideration we expect to be entitled to, which is generally equal to our contract amounts. In some cases, our contracts contain variable consideration which primarily relates to (1) sales and usage-based royalties related to the license of intellectual property in collaboration and license contracts and (2) contracts with minimum purchase commitments. For sales and usage-based royalties, ASC 606 provides an exception to estimating variable consideration. Under this exception, we recognize revenues from sales or usage-based royalty revenue at the later of when the sales or usage occurs or the satisfaction (or partial satisfaction) of the performance obligation to which the royalty has been allocated. All other variable amounts are constrained to the minimum guaranteed contract amount so that a reversal of cumulative revenue does not occur in future periods. Once there is no longer uncertainty over a variable amount, any incremental fees we are entitled to are allocated to the related performance obligations.

Our contracts may include either a single promise (referred to as a performance obligation) to transfer a product or service, or a combination of multiple promises to transfer products or services. We evaluate the existence of multiple promises within our contracts by using judgment to determine if (1) the customer can benefit from each contractual promise on its own or together with readily available resources and (2) the transfer of each contractual promise is separately identifiable from other promises in a contract. When both criteria are met, each promise is accounted for as a separate performance obligation.

Sales of instruments directly to customers include installation and an initial year service-type warranty (which guarantees that our instruments are free from material defects in workmanship and materials, excluding normal wear and tear, and maintenance services). We have determined that the instrument and installation are a combined performance obligation in cases where a customer has a contractual acceptance right that applies through the installation of the instrument. The included service-type warranty is considered a separate performance obligation since a customer could benefit from it independently with readily available resources and is capable of being sold on its own.

Sales of instruments to distributors include a license to import and resell the instruments and an initial year of assurance-type warranty (which guarantees that the products conform to our published specifications). We have determined that the instrument and distributor license are a combined performance obligation since the distributor only benefits from the combination of the instrument and ability to resell it. The assurance-type warranty does not create a separate performance obligation under ASC 606. Under ASC Topic 460 – *Guarantees*, we establish an accrual for estimated assurance-type warranty expense, which is recorded in cost of product revenue on the Consolidated Statements of Operations.

Instrument sales may also be bundled with assays and other consumables, training, and/or an extended service warranty, each of which is considered a separate performance obligation.

Contracts that include rights to additional products or services that are exercisable at a customer's discretion are generally considered options. We assess if these options provide a material right to the customer and if so, the material right is considered a performance obligation. The identification of material rights requires judgment to determine if the value of the option to purchase additional products and services in relation to options that may be provided to, and prices paid by, customers in the normal course of business. Material rights are recognized when they are exercised by a customer or upon expiration of the right.

For contracts that contain multiple performance obligations, the transaction price is allocated among the performance obligations on a relative basis according to their standalone selling prices ("SSP"). Determining the SSP for performance obligations requires judgment. We determine SSP based on factors including prices charged to customers in observable transactions, internal pricing objectives and list prices, pricing of similar products, expected costs to manufacture our products, and estimated margins. We have more than one range of standalone selling price for certain products and services based on the geographic location of the customer and sales channel.

The majority of our products and services are recognized at the point in time we transfer control to the customer.

For product revenues, direct instrument sales that require installation prior to contractual acceptance, the combined performance obligation is recognized upon completion of the instrument's installation. For direct instrument sales that do not contain contractual acceptance and instrument sales to distributors, revenue is recognized based on the agreed upon shipping terms (either upon shipment or delivery) as that is when transfer of control passes to the customer.

Service revenues generated from contract research services in our Accelerator Laboratory are generally recognized over time, using an output method based on the number of completed test results, since the work performed does not have an alternative use and we maintain a contractual right to payment for service performed (including a reasonable profit margin). For performance obligations that are not recognized over time, they are recognized at the point in time when we complete and deliver the research results. Service revenues generated from warranties and service contracts are recognized ratably over the service period as the customer simultaneously receives and benefits from the services.

Collaboration and license revenues are recognized at the point in time the license performance obligation is delivered as the customer has the right to use the intellectual property when it is received. Royalty revenues that are sales or usage-based are recognized at the later of when the sales or usage occurs or the satisfaction (or partial satisfaction) of the performance obligation to which the royalty has been allocated.

Acquired Goodwill, Intangible Assets, and Contingent Liabilities

When acquiring a business, we determine the estimated fair values of the assets acquired and liabilities assumed as of the acquisition date, which may include a significant amount of intangible assets such as customer relationships, developed technology, in-process research and development, trademarks and trade names, know-how, and non-compete agreements, as well as goodwill, contingent liabilities, and off-market contract assets and liabilities.

The determination of the fair values of these assets and liabilities involves significant judgment in selecting inputs used in a valuation methodology, including, but not limited to, projected revenues and expenses, future changes in technology, estimated selling prices, replacement costs or margins, customer attrition rates, covenants not to compete, obsolescence of developed technologies, the likelihood and timing of achieving milestones or performance targets, discount rates, and assumptions about the period of time a brand will continue to be used. We typically engage third-party valuation experts to assist us with the fair value analyses. Our estimates of fair value are based upon assumptions and inputs we believe to be reasonable, but which are inherently uncertain and unpredictable and, as a result, actual results may differ from estimates. A change in the inputs used could have a material impact on the estimated fair values.

Intangible assets, which consist of customer relationships, developed technology, in-process research and development, know-how, trademarks and trade names, and non-compete agreements and are recorded at their fair values as described above. Definite lived intangibles are amortized over each asset's useful life on a basis which best matches the periods in which the economic benefits are expected to be realized. Determining an intangible asset's useful life requires significant judgment and is based on evaluating a number of factors, including, but not limited to, the expected use of the asset, estimated period the asset will generate cash flows, historical client retention rates, consumer awareness, trademark and trade name history, and any contractual provisions that could limit or extend an asset's useful life. Actual useful lives may differ from estimated useful lives. Indefinite-lived intangibles, which consist of in-process research and development, are not amortized until the underlying project is completed. Upon completion, the in-process research and development asset is accounted for as a definite lived intangible asset.

Acquired customer contracts may contain off-market terms. We record an additional contract asset or liability for such favorable or unfavorable terms at their estimated acquisition date fair values. Determining the fair value involves significant judgment to assess the economic returns that could be realized in a market transaction and can include the following inputs in a valuation methodology: projected revenue, expected profit margin, and discount rates.

Business combinations may also include contingent liabilities to be paid based on the occurrence of future events, such as the completion of a technical milestone or upon meeting certain performance targets. Contingent liabilities treated as purchase price are recorded at fair value, as described above, at the acquisition date. We remeasure the fair value of outstanding contingent liabilities at each reporting period and changes are recognized in change in fair value of contingent liabilities on the Consolidated Statements of Operations.

Impairment of Goodwill and Indefinite-Lived Intangible Assets

Goodwill is required to be assessed for impairment at least annually or whenever events or circumstances indicate that there may be an impairment. Absent an event that indicates a specific impairment may exist, we have selected October 1st as the date to perform our annual goodwill impairment test.

An impairment assessment requires evaluating a potential impairment at the reporting unit level using either a qualitative assessment, to determine if it is more likely than not that the fair value of any reporting unit is less than its carrying amount, or a quantitative analysis, to determine and compare the fair value of each reporting unit to its carrying value, or a combination of both. Reporting units are determined based on the components of our operating segments that constitute a business for which financial information is available and for which operating results are regularly reviewed by segment management. Judgment is required in determining the use of a qualitative or quantitative assessment, as well as in determining a reporting unit's estimated fair value, as it requires us to make estimates of market conditions and operational performance, including items such as projected financial results, discount rates, control premiums, or valuation multiples for key financial metrics.

We completed our annual goodwill impairment test as of October 1, 2025 utilizing a qualitative assessment to determine if it was more likely than not that the fair values of each of our reporting units was less than their respective carrying values and concluded that no impairments existed.

At least annually or whenever events or circumstances change, we assess whether an indefinite-lived intangible assets has been abandoned, in which case it would be written off, or if its estimated fair value is below its carrying value, in which case it could be impaired.

Impairment of Long-Lived Assets

We continually evaluate whether events or circumstances have occurred that indicate the carrying value of any of our definite lived intangible assets or other long-lived assets may not be recoverable, or the estimated remaining useful life or estimated remaining lease term may require revision. Additionally, we continually assess the determination of our asset groups, which primarily focuses on changes in our operating structure, the way in which we expect to deploy our assets, or how we intend to recover the cost of our assets.

To assess the recoverability of a long-lived asset or asset group, we compare the estimated undiscounted future cash flows for the estimated remaining useful life, or estimated lease term, of the asset (or the primary asset in the asset group) to its carrying value. If the undiscounted cash flows are less than the carrying value, we estimate the asset's fair value using the future discounted cash flows associated with the use of the asset. To the extent that the discounted cash flows are less than the carrying value, the asset(s) are impaired and written down to their estimated fair value.

Significant assumptions that form the basis of the forecasted results utilized to calculate undiscounted cash flows include but are not limited to, estimates about future revenues, expenses, asset disposal value, expected uses of the asset, historical customer retention rates, technology roadmaps, customer awareness, trademark and trade name history, contractual provisions that could limit or extend an asset's useful life, market data, discount rates, and potential sublease opportunities including rent and rent escalation rates, time to sublease, and free rent periods. To the extent that the future cash flows are less than the carrying value, a long-lived asset or asset group is impaired and written down to its estimated fair value.

Future events could cause us to conclude that impairment indicators exist and that goodwill, intangible assets, or other long-lived assets are impaired. Any resulting impairment loss could have a material adverse impact on our results of operations.

Non-GAAP Financial Measures

To supplement our financial statements presented on a U.S. GAAP basis, we present the following non-GAAP financial measures: adjusted gross profit, adjusted gross margin, adjusted total operating expenses, and adjusted loss from operations. These non-GAAP measures are calculated by including shipping and handling costs for product sales within cost of product revenue instead of within selling, general and administrative expenses and excluding amortization of certain acquired intangible assets, acquisition and integration related costs, and certain other items which include other charges or benefits resulting from transactions or events that are highly variable, significant in size, and that we do not believe are indicative of ongoing or future business operations. Adjusted gross margin is calculated as adjusted gross profit divided by total revenues.

We believe that the presentation of these non-GAAP financial measures provides supplemental information useful to investors in understanding our underlying operating results and trends. We use these non-GAAP financial measures to evaluate our operating performance in a manner that allows for meaningful period-to-period comparison and analysis of

trends in our business and our competitors. We believe that presentation of these non-GAAP financial measures provides useful information to investors in assessing our operating performance within our industry and to allow comparability with the presentation of other companies in our industry.

The non-GAAP financial measures presented here should be considered in conjunction with, and not as a substitute for, the financial information presented in accordance with U.S. GAAP.

Set forth below is a reconciliation of adjusted gross profit, adjusted gross margin, adjusted total operating expenses, and adjusted loss from operations from their most directly comparable GAAP financial measures (in thousands, except percentages, unaudited):

	Three Months Ended December 31,		Year Ended December 31,	
	2025	2024	2025	2024
Gross profit	\$ 20,038	\$ 22,169	\$ 64,959	\$ 83,104
Shipping and handling costs	(1,400)	(1,885)	(5,581)	(8,113)
Purchase accounting impact on inventory and property and equipment (1)	356	—	391	—
Amortization of acquired intangible assets (2)	2,953	—	5,946	—
Adjusted gross profit (non-GAAP)	\$ 21,947	\$ 20,284	\$ 65,715	\$ 74,991
Total revenues	\$ 43,855	\$ 35,161	\$ 138,897	\$ 137,421
Gross margin (gross profit as % of total revenues)	45.7%	63.0%	46.8%	60.5%
Adjusted gross margin (non-GAAP) (adjusted gross profit as % of total revenues)	50.0%	57.7%	47.3%	54.6%
Total operating expenses	\$ 44,789	\$ 36,938	\$ 190,501	\$ 135,720
Shipping and handling costs	(1,400)	(1,885)	(5,581)	(8,113)
Purchase accounting impact on property and equipment (1)	(416)	—	(628)	—
Amortization of acquired intangible assets (2)	(80)	—	(153)	—
Acquisition and integration related costs (3)	(1,384)	(1,100)	(16,416)	(1,100)
Earnout recorded as compensation expense (4)	(1,871)	—	(10,000)	—
Impairments and employee separation costs (5)	(2,687)	—	(17,531)	—
Adjusted total operating expenses (non-GAAP)	\$ 36,951	\$ 33,953	\$ 140,192	\$ 126,507
Loss from operations	\$ (24,751)	\$ (14,769)	\$ (125,542)	\$ (52,616)
Purchase accounting impact on property and equipment (1)	772	—	1,019	—
Amortization of acquired intangible assets (2)	3,033	—	6,099	—
Acquisition and integration related costs (3)	1,384	1,100	16,416	1,100
Earnout recorded as compensation expense (4)	1,871	—	10,000	—
Impairments and employee separation costs (5)	2,687	—	17,531	—
Adjusted loss from operations (non-GAAP)	\$ (15,004)	\$ (13,669)	\$ (74,477)	\$ (51,516)

(1) Represents the amortization of the purchase price fair value increase of acquired inventory and property and equipment.

(2) Consists only of the amortization of intangible assets acquired in 2025.

(3) Represents acquisition and integration costs directly related to the Company's business combinations. Acquisition costs include professional and consulting fees supporting due diligence, legal, and accounting activities to execute a transaction. Integration costs include third party and internal direct costs to integrate acquired companies, employees, and their customers.

(4) Consists of the earnout recognized as compensation expense related to the Emission acquisition.

(5) Impairment charges for goodwill and acquired leased facilities not in use, as well as one-time severance and benefit costs.

Recent Accounting Pronouncements

Refer to Note 2 - *Significant Accounting Policies* in the Notes to Consolidated Financial Statements included in this Annual Report on Form 10-K for a full description of recent accounting pronouncements, including the expected dates of adoption and effects on our Consolidated Financial Statements.

ITEM 7A. QUANTITATIVE AND QUALITATIVE DISCLOSURES ABOUT MARKET RISK

We are exposed to a variety of market risks, including fluctuations in foreign currency exchange rates and interest rates affecting the return on our cash, cash equivalents, and marketable securities.

Foreign Currency Exchange Risk

As we expand internationally, our results of operations and cash flows will become increasingly subject to foreign exchange rate fluctuations. For the years ended December 31, 2025 and 2024, approximately 40% and 37%, respectively, of our total revenue was generated from customers located outside of the United States. Our expenses are generally denominated in the currencies in which our operations are located, which is primarily in the United States, with a portion of expenses incurred in Canada, China, Europe, and Japan. Our results of operations and cash flows are, therefore, subject to fluctuations due to changes in foreign exchange rates in Canadian dollars, Euros, British pounds, Swedish krona, Japanese yen, Chinese yuan, and other foreign currencies. Fluctuations in exchange rates could harm our business in the future. As of December 31, 2025, the effect of a hypothetical 10% adverse change in exchange rates on foreign denominated cash and payables would not have been material and a similar adverse change on foreign denominated receivables would decrease potential cash inflows by \$0.6 million.

To date, we have not entered into any foreign currency hedging contracts although we may do so in the future.

Interest Rate Risk

We had cash and cash equivalents of \$29.8 million and marketable securities of \$88.4 million as of December 31, 2025. All cash, cash equivalents, and marketable securities are held at large commercial banks. Marketable securities consisted entirely of highly rated debt securities including commercial paper, U.S. Treasuries, corporate notes and bonds, U.S. Government agency bonds, and similar types of debt securities. Due to the short-term nature and investment grade quality of these investments, we do not believe we have material exposure to changes in interest rates. Additionally, if needed, we have the ability to hold our marketable securities until maturity (without giving effect to any future acquisitions or mergers) and we do not hold or issue financial instruments for trading purposes. Therefore, we do not expect our operating results or cash flows to be affected materially by a sudden change in market interest rates.

Declines in interest rates, however, would reduce future investment income. If overall interest rates had decreased by a hypothetical 10% during the year ended December 31, 2025, our interest income would have decreased by approximately \$0.9 million.

ITEM 8. FINANCIAL STATEMENTS AND SUPPLEMENTARY DATA

The financial statements required to be filed pursuant to Item 8 are included in this Annual Report on Form 10-K beginning on page F-1.

ITEM 9. CHANGES IN AND DISAGREEMENTS WITH ACCOUNTANTS ON ACCOUNTING AND FINANCIAL DISCLOSURE

Not applicable.

ITEM 9A. CONTROLS AND PROCEDURES

Evaluation of Disclosure Controls and Procedures

We have established disclosure controls and procedures (as such term is defined in Rules 13a-15(e) and 15d-15(e) under the Exchange Act) that are designed to provide reasonable assurance that information required to be disclosed in the reports we file or submit under the Exchange Act is recorded, processed, summarized, and reported within the time periods specified in the rules and forms of the SEC and to ensure that such information is accumulated and communicated to management, including our Chief Executive Officer (principal executive officer) and Chief Financial Officer (principal financial officer), to allow timely decisions regarding required disclosures.

Under the supervision and with the participation of our management, including our Chief Executive Officer and Chief Financial Officer, we conducted an evaluation of the effectiveness of our disclosure controls and procedures as of

December 31, 2025. Based on this evaluation, our Chief Executive Officer and Chief Financial Officer concluded that our disclosure controls and procedures were effective as of December 31, 2025.

Management's Report on Internal Control over Financial Reporting

Our management is responsible for establishing and maintaining adequate internal control over financial reporting, as this term is defined in Rules 13a-15(f) and 15d-15(f) under the Exchange Act. Our internal control over financial reporting is a process designed to provide reasonable assurance regarding the reliability of financial reporting and the preparation of financial statements for external purposes in accordance with generally accepted accounting principles. All internal control systems, no matter how well designed, have inherent limitations. Therefore, even those systems determined to be effective can provide only reasonable assurance with respect to financial statement preparation and presentation.

Under the supervision and with the participation of our management, including our Chief Executive Officer (principal executive officer) and Chief Financial Officer (principal financial officer), we conducted an evaluation of the effectiveness of our internal control over financial reporting as of December 31, 2025, based on the framework set forth in Internal Control-Integrated Framework issued by the Committee of Sponsoring Organizations of the Treadway Commission (the "2013 Framework"). Based on our evaluation under the 2013 Framework, management concluded that our internal control over financial reporting was effective as of December 31, 2025.

The foregoing evaluation excludes our acquisition of Akoya Biosciences, Inc. ("Akoya") on July 8, 2025 (refer to Note 3 - *Acquisitions* in the Notes to Consolidated Financial Statements). The acquisition of Akoya was accounted for as a business combination and Akoya's operating results have been included in our Consolidated Financial Statements since the acquisition date. In accordance with guidance from the staff of the SEC, companies are permitted to exclude acquisitions from their evaluation of internal control over financial reporting for the first year after an acquisition. Akoya represented approximately 43% of the Company's total assets, excluding the preliminary value of goodwill, and 24% of the Company's total revenues as of and for the year ended December 31, 2025.

The Company's independent registered public accounting firm, KPMG LLP, has also issued an audit report on the Company's internal control over financial reporting, which is included elsewhere in this Annual Report on Form 10-K.

Remediation of Previously Identified Material Weaknesses

In preparing our financial statements in connection with our Annual Report on Form 10-K for the year ended December 31, 2024 and continuing through the nine months ended September 30, 2025, we previously identified and disclosed material weaknesses in our internal controls associated with the valuation of inventory, including excess and obsolescence reserves and the capitalization of labor and overhead costs and the accounting for Accelerator Laboratory revenue, a component of our service and other revenue.

Following the identification of the material weaknesses, and with the oversight of our audit committee, we commenced remediation efforts which continued during fiscal 2025 to address the material weaknesses and enhance our control environment, including our internal control over financial reporting. Our remediation efforts included:

- hiring a Vice President, SOX Transformation and personnel for our internal SOX Transformation team. This team:
 - oversaw the remediation of our material weaknesses and improvements across our internal controls;
 - evaluated and designed effective and scalable internal controls, and strengthened the documentation of our existing controls;
 - established new internal controls evaluating the accounting for inventory and enhanced inventory valuation review procedures;
 - enhanced and expanded our existing revenue recognition control procedures and attributes to sufficiently document our assessment of, and reviews over, information used to record Accelerator Laboratory revenue; and
 - provided trainings on a regular basis related to internal control over financial reporting for all control owners.
- implementing new software solutions to automate key manual inventory valuation processes and outputs;

- designing and implementing new controls;
- bringing additional internal systems into the scope of our internal control environment to reduce reliance on manual processes and controls;
- continued execution of controls that we worked to improve during fiscal year 2024 that did not have a sufficient period of time to demonstrate operating effectiveness as of December 31, 2024, including the analysis of labor and overhead cost capitalization and related controls implemented in the fourth quarter of 2024; and
- evaluating, enhancing, and adding personnel in the finance organization with a focus on the requisite experience in the areas of accounting, SEC financial reporting, and internal control compliance.

Based on these remediation actions, as well as testing the operating effectiveness of the applicable financial reporting controls over a sustained period of financial reporting cycles, we have concluded that all previously reported material weaknesses have been effectively remediated as of December 31, 2025.

Changes in Internal Control over Financial Reporting

Other than the changes outlined related to the remediation of the material weaknesses described above, there have been no changes in our internal control over financial reporting during the fiscal quarter ended December 31, 2025 that have materially affected, or are reasonably likely to materially affect, our internal control over financial reporting.

ITEM 9B. OTHER INFORMATION

Securities Trading Plans of Directors and Executive Officers

During the three months ended December 31, 2025, none of our directors or officers adopted or terminated any contract, instruction or written plan for the purchase or sale of our securities that was intended to satisfy the affirmative defense conditions of Rule 10b5-1(c) or any "non-Rule 10b5-1 trading arrangement" (as defined in Item 408(c) of Regulation S-K).

ITEM 9C. DISCLOSURE REGARDING FOREIGN JURISDICTIONS THAT PREVENT INSPECTIONS

Not applicable.

PART III

ITEM 10. DIRECTORS, EXECUTIVE OFFICERS AND CORPORATE GOVERNANCE

The information required by this Item will be set forth in our Proxy Statement for the 2026 Annual Meeting of Stockholders to be filed with the SEC within 120 days of December 31, 2025 and is incorporated into this Annual Report on Form 10-K by reference.

ITEM 11. EXECUTIVE COMPENSATION

The information required by this Item will be set forth in our Proxy Statement for the 2026 Annual Meeting of Stockholders to be filed with the SEC within 120 days of December 31, 2025 and is incorporated into this Annual Report on Form 10-K by reference.

ITEM 12. SECURITY OWNERSHIP OF CERTAIN BENEFICIAL OWNERS AND MANAGEMENT AND RELATED STOCKHOLDER MATTERS

The information required by this Item will be set forth in our Proxy Statement for the 2026 Annual Meeting of Stockholders to be filed with the SEC within 120 days of December 31, 2025 and is incorporated into this Annual Report on Form 10-K by reference.

ITEM 13. CERTAIN RELATIONSHIPS AND RELATED TRANSACTIONS, AND DIRECTOR INDEPENDENCE

The information required by this Item will be set forth in our Proxy Statement for the 2026 Annual Meeting of Stockholders to be filed with the SEC within 120 days of December 31, 2025 and is incorporated into this Annual Report on Form 10-K by reference.

ITEM 14. PRINCIPAL ACCOUNTANT FEES AND SERVICES

Our independent registered public accounting firm is KPMG LLP, New York, NY, (PCAOB ID 185).

The other information required by this Item will be set forth in our Proxy Statement for the 2026 Annual Meeting of Stockholders to be filed with the SEC within 120 days of December 31, 2025 and is incorporated into this Annual Report on Form 10-K by reference.

PART IV

ITEM 15. EXHIBITS AND FINANCIAL STATEMENT SCHEDULES

(1) *Financial Statements*

The Consolidated Financial Statements are included beginning on page F-1 attached hereto and are filed as part of this Annual Report on Form 10-K.

(2) *Financial Statement Schedules*

Financial statement schedules have been omitted since they either are not required, not applicable, or the information is otherwise included.

(3) *Exhibits*

The following is a list of exhibits filed as part of this Annual Report on Form 10-K:

Exhibit Number	Exhibit Description	Filed Herewith	Incorporated by Reference herein from Form or Schedule	Filing Date	SEC File/Reg. Number
2.1@	Share Purchase Agreement, dated December 16, 2024, by and among Quanterix Corporation, Emission Inc., the Shareholders of Emission Inc. and the Shareholder Representative		8-K	12/17/2024	001-38319
2.2#	Amended and Restated Agreement and Plan of Merger, dated March 28, 2025, by and among the Registrant, Wellfleet Merger Sub, Inc., and Akoya Biosciences, Inc.		8-K	4/29/2025	001-38319
3.1	Amended and Restated Certificate of Incorporation		8-K	10/2/2025	001-38319
3.2	Restated Bylaws		8-K	10/2/2025	001-38319
4.1	Description of Securities		10-K	3/13/2020	001-38319
4.2	Form of Common Stock Certificate		S-1	11/9/2017	333-221475
10.1.1+	2007 Stock Option and Grant Plan, as amended		S-1	11/9/2017	333-221475
10.1.2+	Form of Incentive Stock Option Agreement under the 2007 Stock Option and Grant Plan, as amended		S-1	11/9/2017	333-221475
10.1.3+	Form of Non-qualified Stock Option Agreement under the 2007 Stock Option and Grant Plan, as amended		S-1	11/9/2017	333-221475
10.1.4+	Form of Restricted Stock Agreement under the 2007 Stock Option and Grant Plan, as amended		S-1	11/9/2017	333-221475
10.2.1+	2017 Employee, Director and Consultant Equity Incentive Plan		S-1/A	11/27/2017	333-221475

[Table of Contents](#)

Exhibit Number	Exhibit Description	Filed Herewith	Incorporated by Reference herein from Form or Schedule	Filing Date	SEC File/ Reg. Number
10.2.2+	Form of Stock Option Agreement under the 2017 Employee, Director and Consultant Equity Incentive Plan		S-1/A	11/27/2017	333-221475
10.2.3+	Form of Restricted Stock Agreement under the 2017 Employee, Director and Consultant Equity Incentive Plan		S-1/A	11/27/2017	333-221475
10.2.4+	Form of Restricted Stock Unit Agreement under the 2017 Employee, Director and Consultant Equity Incentive Plan		10-K	2/29/2024	001-30319
10.3.1+	Akoya Biosciences, Inc. 2021 Equity Incentive Plan and form of stock option agreement thereunder		S-1/A	4/12/2021	333-254760
10.3.2+	Akoya Biosciences, Inc. Form of Indemnification Agreement		8-K	12/13/2024	001-40344
10.4.1+	Amended and Restated 2025 Inducement Plan		S-8	1/15/2026	333-292362
10.4.2+	Form of Stock Option Agreement under the 2025 Inducement Plan		S-8	12/22/2025	333-292362
10.4.3+	Form of Restricted Stock Unit Agreement under the 2025 Inducement Plan		S-8	12/22/2025	333-292362
10.5.1+	Employment Agreement, dated August 3, 2023, between the Registrant and Vandana Sriram		8-K	8/9/2023	001-38319
10.5.2+	Amendment, effective as of April 11, 2024, to the Employment Agreement, dated August 3, 2023, between the Registrant and Vandana Sriram		8-K	4/12/2024	001-38319
10.7.1+	Amended and Restated Employment Agreement, dated April 25, 2022, between the Registrant and Dr. Masoud Toloue		8-K	4/29/2022	001-38319
10.7.2+	Amendment, dated April 9, 2024, to the Amended and Restated Employment Agreement, dated April 25, 2022, between the Registrant and Dr. Masoud Toloue		8-K	4/12/2024	001-38319
10.7.3+	Separation Agreement by and between the Company and Masoud Toloue		8-K	1/8/2026	001-38319
10.8+	Employment Agreement by and between the Company and William Donnelly	X			
10.9+	Employment Agreement by and between the Company and Everett Cunningham		8-K	1/8/2026	001-38319

[Table of Contents](#)

Exhibit Number	Exhibit Description	Filed Herewith	Incorporated by Reference herein from Form or Schedule	Filing Date	SEC File/ Reg. Number
10.10.1*	Exclusive License Agreement, dated June 18, 2007, between the Registrant and Tufts University, as amended on April 29, 2013		S-1	11/9/2017	333-221475
10.10.2*	Second Amendment, dated August 22, 2017, to the Exclusive License Agreement between the Registrant and Tufts University		S-1	11/9/2017	333-221475
10.10.3@	Third Amendment, dated September 25, 2020, to the Exclusive License Agreement between the Registrant and Tufts University		10-Q	11/6/2020	001-38319
10.11.1*	Supply and Manufacturing Agreement, dated September 14, 2011, between the Registrant and STRATEC Biomedical AG		S-1	11/9/2017	333-221475
10.11.2	First Amendment to Supply and Manufacturing Agreement, dated October 17, 2013, between the Registrant and STRATEC Biomedical AG		S-1	11/9/2017	333-221475
10.12.1*	STRATEC Development Services and Equity Participation Agreement, dated August 15, 2011, between the Registrant and STRATEC Biomedical Systems AG		S-1	11/9/2017	333-221475
10.12.2*	First Amendment to STRATEC Development Services and Equity Participation Agreement and Second Amendment to Supply and Manufacturing Agreement, dated November 18, 2016, between the Registrant and STRATEC Biomedical AG		S-1	11/9/2017	333-221475
10.13@	Third Amendment to Supply and Manufacturing Agreement, entered into on June 17, 2025 and effective as of January 1, 2025 between the Registrant and STRATEC SE		10-Q	8/7/2025	001-38319
10.14*	Manufacturing Services Agreement, dated November 23, 2016, between the Registrant and Paramit Corporation		S-1	11/9/2017	333-221475
10.15*	Exclusivity (Equity) Agreement, dated November 17, 2015, by and between Akoya Biosciences, Inc. and The Board of Trustees of the Leland Stanford Junior University		S-1	3/26/2021	333-254760
10.16*	Amendment No. 1 to the License Agreement, dated November 18, 2016, by and between Akoya Biosciences, Inc. and the Board of Trustees of the Leland Stanford Junior University		S-1	3/26/2021	333-254760
10.17*	Manufacturing Agreement, dated July 17, 2019 between Akoya Biosciences, Inc. and Columbia Electrical Contractors, Inc.	X			

Exhibit Number	Exhibit Description	Filed Herewith	Incorporated by Reference herein from Form or Schedule	Filing Date	SEC File/ Reg. Number
10.18*	License and Royalty Agreement, dated September 28, 2018, by and among Akoya Biosciences, Inc., PerkinElmer Health Services, Inc., Cambridge Research and Instrumentation, Inc., and VisEn Medical Inc.		S-1	3/26/2021	333-254760
10.19*	Exclusive Patent License Agreement, dated June 26, 2018, by and between Akoya Biosciences, Inc. and the University of Washington		S-1	3/26/2021	333-254760
10.20	Cooperation Agreement, by and between Quanterix Corporation, Kent Lake PR LLC and Kent Lake Partners LP, dated August 4, 2025		8-K	8/4/2024	001-38319
10.21+	Form of Indemnification Agreement		S-1/A	11/27/2017	333-221475
10.22	Lease Agreement by and between SSI 900 Middlesex MA LP and the Registrant, dated October 2, 2018		8-K	10/5/2018	001-38319
10.23	Lease Agreement by and between the Registrant and XChange Owner LLC, dated January 28, 2022		8-K	1/31/2022	001-38319
10.24+	Amended and Restated Non-Employee Director Compensation Policy	X			
10.25.1#	Voting and Support Agreement, dated April 28, 2025, by and among the Registrant and certain stockholders of Akoya Biosciences, Inc. named therein.		8-K	4/29/2025	001-38319
10.25.2*	Consent and Waiver under the Voting and Support Agreement, dated April 28, 2025, by and among the Registrant and certain stockholders of Akoya Biosciences, Inc. named therein		8-K	4/29/2025	001-38319
10.26.1	Securities Purchase Agreement between the Registrant and Akoya Biosciences, Inc. dated April 2, 2025		8-K	4/4/2025	001-38319
10.26.2	Amendment No. 1, dated April 28, 2025, to the Securities Purchase Agreement between the Registrant and Akoya Biosciences, Inc. dated April 2, 2025		8-K	4/29/2025	001-38319
19.1	Quanterix Insider Trading Policy	X			
21.1	Subsidiaries of Registrant	X			
23.1	Consent of KPMG LLP	X			
23.2	Consent of Ernst & Young LLP	X			

Exhibit Number	Exhibit Description	Filed Herewith	Incorporated by Reference herein from Form or Schedule	Filing Date	SEC File/Reg. Number
31.1	Certification of the Principal Executive Officer pursuant to Section 302 of the Sarbanes-Oxley Act of 2002	X			
31.2	Certification of the Principal Financial Officer pursuant to Section 302 of the Sarbanes-Oxley Act of 2002	X			
32.1	Certifications of the Chief Executive Officer and Chief Financial Officer pursuant to Section 906 of the Sarbanes-Oxley Act of 2002	X			
97.1	Policy relating to recovery of erroneously awarded compensation		10-K	2/29/2024	001-38319
101.INS	Inline XBRL Instance Document	X			
101.SCH	Inline XBRL Taxonomy Extension Schema Document	X			
101.CAL	Inline XBRL Taxonomy Extension Calculation Linkbase Document	X			
101.DEF	Inline XBRL Taxonomy Extension Definition	X			
101.LAB	Inline XBRL Taxonomy Extension Label Linkbase Document	X			
101.PRE	Inline XBRL Taxonomy Presentation Linkbase Document	X			
104	Cover Page Interactive Data File (formatted as Inline XBRL and contained in Exhibit 101)				

+ Management contract or compensatory plan or arrangement.

* Confidential treatment has been granted for portions of this Exhibit. Redacted portions have been filed separately with the SEC.

@ Portions of this document (indicated by "[***]") have been omitted because such information is not material and is the type of information that the Registrant treats as private or confidential.

Certain exhibits and schedules have been omitted pursuant to Item 601(b)(2) of Regulation S-K. The Registrant hereby undertakes to furnish supplemental copies of any of the omitted exhibits and schedules upon request by the SEC; provided, however, that the Registrant may request confidential treatment pursuant to Rule 24b-2 of the Securities Exchange Act of 1934, as amended, for any exhibits or schedules so furnished.

ITEM 16. FORM 10-K SUMMARY

Not applicable.

QUANTERIX CORPORATION
INDEX TO CONSOLIDATED FINANCIAL STATEMENTS

	<u>Page</u>
Report of Independent Registered Public Accounting Firm (PCAOB ID: 185)	F-2
Report of Independent Registered Public Accounting Firm (PCAOB ID: 42)	F-5
Consolidated Balance Sheets	F-6
Consolidated Statements of Operations	F-7
Consolidated Statements of Comprehensive Loss	F-8
Consolidated Statements of Stockholders' Equity	F-9
Consolidated Statements of Cash Flows	F-10
Notes to Consolidated Financial Statements	F-11

REPORT OF INDEPENDENT REGISTERED PUBLIC ACCOUNTING FIRM

To the Stockholders and the Board of Directors
Quanterix Corporation:

Opinions on the Consolidated Financial Statements and Internal Control Over Financial Reporting

We have audited the accompanying consolidated balance sheet of Quanterix Corporation and subsidiaries (the Company) as of December 31, 2025, the related consolidated statements of operations, comprehensive loss, stockholders' equity, and cash flows for the year ended December 31, 2025, and the related notes collectively, the consolidated financial statements. We also have audited the Company's internal control over financial reporting as of December 31, 2025, based on criteria established in Internal Control – Integrated Framework (2013) issued by the Committee of Sponsoring Organizations of the Treadway Commission.

In our opinion, the consolidated financial statements referred to above present fairly, in all material respects, the financial position of the Company as of December 31, 2025 and the results of its operations and its cash flows for the year ended December 31, 2025, in conformity with U.S. generally accepted accounting principles. Also in our opinion, the Company maintained, in all material respects, effective internal control over financial reporting as of December 31, 2025 based on criteria established in Internal Control – Integrated Framework (2013) issued by the Committee of Sponsoring Organizations of the Treadway Commission.

The Company acquired Akoya Biosciences, Inc. during 2025, and management excluded from its assessment of the effectiveness of the Company's internal control over financial reporting as of December 31, 2025, Akoya Biosciences, Inc.'s internal control over financial reporting associated with 43% of total assets, excluding the preliminary value of goodwill and 24% of the total revenues included in the consolidated financial statements of the Company as of and for the year ended December 31, 2025. Our audit of internal control over financial reporting of the Company also excluded an evaluation of the internal control over financial reporting of Akoya Biosciences, Inc.

Basis for Opinions

The Company's management is responsible for these consolidated financial statements, for maintaining effective internal control over financial reporting, and for its assessment of the effectiveness of internal control over financial reporting, included in the accompanying Management's Report on Internal Control over Financial Reporting. Our responsibility is to express an opinion on the Company's consolidated financial statements and an opinion on the Company's internal control over financial reporting based on our audits. We are a public accounting firm registered with the Public Company Accounting Oversight Board (United States) (PCAOB) and are required to be independent with respect to the Company in accordance with the U.S. federal securities laws and the applicable rules and regulations of the Securities and Exchange Commission and the PCAOB.

We conducted our audits in accordance with the standards of the PCAOB. Those standards require that we plan and perform the audits to obtain reasonable assurance about whether the consolidated financial statements are free of material misstatement, whether due to error or fraud, and whether effective internal control over financial reporting was maintained in all material respects.

Our audit of the consolidated financial statements included performing procedures to assess the risks of material misstatement of the consolidated financial statements, whether due to error or fraud, and performing procedures that respond to those risks. Such procedures included examining, on a test basis, evidence regarding the amounts and disclosures in the consolidated financial statements. Our audit also included evaluating the accounting principles used and significant estimates made by management, as well as evaluating the overall presentation of the consolidated financial statements. Our audit of internal control over financial reporting included obtaining an understanding of internal control over financial reporting, assessing the risk that a material weakness exists, and testing and evaluating the design and operating effectiveness of internal control based on the assessed risk. Our audits also included performing such other procedures as we considered necessary in the circumstances. We believe that our audits provide a reasonable basis for our opinions.

Definition and Limitations of Internal Control Over Financial Reporting

A company's internal control over financial reporting is a process designed to provide reasonable assurance regarding the reliability of financial reporting and the preparation of financial statements for external purposes in accordance with generally accepted accounting principles. A company's internal control over financial reporting includes those policies and procedures that (1) pertain to the maintenance of records that, in reasonable detail, accurately and fairly reflect the transactions and dispositions of the assets of the company; (2) provide reasonable assurance that transactions are recorded as necessary to permit preparation of financial statements in accordance with generally accepted accounting principles, and that receipts and expenditures of the company are being made only in accordance with authorizations of management and directors of the company; and (3) provide reasonable assurance regarding prevention or timely detection of unauthorized acquisition, use, or disposition of the company's assets that could have a material effect on the financial statements.

Because of its inherent limitations, internal control over financial reporting may not prevent or detect misstatements. Also, projections of any evaluation of effectiveness to future periods are subject to the risk that controls may become inadequate because of changes in conditions, or that the degree of compliance with the policies or procedures may deteriorate.

Critical Audit Matters

The critical audit matters communicated below are matters arising from the current period audit of the consolidated financial statements that were communicated or required to be communicated to the audit committee and that: (1) relate to accounts or disclosures that are material to the consolidated financial statements and (2) involved our especially challenging, subjective, or complex judgments. The communication of critical audit matters does not alter in any way our opinion on the consolidated financial statements, taken as a whole, and we are not, by communicating the critical audit matters below, providing separate opinions on the critical audit matters or on the accounts or disclosures to which they relate.

Recoverability of an asset group

As discussed in Notes 2 and 4 to the consolidated financial statements, the Company reviews its long-lived assets to evaluate whether events or circumstances have occurred that indicate that the carrying amount of such assets may not be fully recoverable or that the estimated remaining useful lives (or estimated remaining lease term) may warrant revision. To assess the recoverability of a long-lived asset or asset group, the Company compares the estimated undiscounted future cash flows for the estimated remaining useful life, or estimated lease term, of the asset (or the primary asset in the asset group) to its carrying value. As of December 31, 2025, long-lived assets included property and equipment, net, definite lived intangible assets, net, and operating lease right-of-use assets of \$23.7 million, \$112.5 million, and \$16.6 million, respectively.

We identified the assessment of the recoverability of an asset group as a critical audit matter. Subjective and challenging auditor judgment, including the use of specialized skills and knowledge, was required to assess the accounting analysis and measurement of the asset disposal value used in the undiscounted cash flows associated with the asset group. Changes in the asset disposal value could have had a significant effect on the Company's assessment of the recoverability of the asset group.

The following are the primary procedures we performed to address this critical audit matter. We evaluated the design and tested the operating effectiveness of an internal control over the accounting analysis and measurement of the disposal value of the asset group. We assessed the reasonableness of the Company's accounting analysis of the asset disposal value used in the undiscounted cash flows associated with the asset group by evaluating the relevant accounting literature. We involved valuation professionals with specialized skills and knowledge, who assisted in evaluating the measurement of the asset group's disposal value used in the undiscounted future cash flows by comparing it to a range of disposal values that were independently developed using publicly available market data for comparable companies.

Acquisition-date fair value of the Akoya developed technology intangible asset

As discussed in Notes 3 and 4 to the consolidated financial statements, the Company acquired Akoya Biosciences, Inc. (Akoya) in July 2025, for total consideration of \$151.0 million. The transaction was accounted for using the acquisition method of accounting with acquired assets and liabilities, including identifiable intangible assets, recorded at their estimated fair values as of the acquisition date. Identifiable intangible assets included developed technology for which the Company estimated the acquisition-date fair value using a multi-period excess earnings valuation methodology. The acquisition-date fair value of the developed technology intangible asset recorded by the Company was \$99.6 million.

We identified the assessment of the acquisition-date fair value of the developed technology intangible asset acquired in the Akoya acquisition as a critical audit matter. A high degree of subjective auditor judgment was required to evaluate certain assumptions used to estimate the acquisition-date fair value of the developed technology intangible asset. Specifically, there was subjectivity in evaluating the projected revenues in certain future periods, and the evaluation of the discount rate required specialized skills and knowledge. Changes to these assumptions could have had a significant effect on the Company's determination of the acquisition-date fair value of the developed technology intangible asset.

The following are the primary procedures we performed to address this critical audit matter. We evaluated the design and tested the operating effectiveness of certain internal controls related to the Company's process to estimate the acquisition-date fair value of the developed technology intangible asset. This included controls related to the determination of projected revenues in certain future periods and discount rate assumptions used to develop the estimate. We performed sensitivity analyses over projected revenues in certain future periods to assess the impact of changes in those assumptions on the Company's determination of the acquisition-date fair value of the developed technology intangible asset. We evaluated the reasonableness of projected revenues in certain future periods by comparing them to Akoya's historical results, industry related third-party data, and revenue trends of comparable companies. In addition, we involved valuation professionals with specialized skills and knowledge, who assisted in:

- evaluating the discount rate used by the Company by comparing it against a discount rate range that was independently developed using publicly available market data for comparable companies
- assessing the acquisition-date fair value of the developed technology intangible asset using the Company's cash flow forecasts and our independently developed discount rate range and comparing the result to the Company's acquisition-date fair value.

/s/ KPMG LLP

We have served as the Company's auditor since 2025.

Boston, Massachusetts

March 2, 2026

REPORT OF INDEPENDENT REGISTERED PUBLIC ACCOUNTING FIRM

To the Stockholders and the Board of Directors of Quanterix Corporation

Opinion on the Financial Statements

We have audited the accompanying consolidated balance sheet of Quanterix Corporation (the Company) as of December 31, 2024, the related consolidated statements of operations, comprehensive loss, stockholders' equity and cash flows for each of the two years in the period ended December 31, 2024, and the related notes (collectively referred to as the "consolidated financial statements"). In our opinion, the consolidated financial statements present fairly, in all material respects, the financial position of the Company at December 31, 2024, and the results of its operations and its cash flows for each of the two years in the period ended December 31, 2024, in conformity with U.S. generally accepted accounting principles.

Basis for Opinion

These financial statements are the responsibility of the Company's management. Our responsibility is to express an opinion on the Company's financial statements based on our audits. We are a public accounting firm registered with the PCAOB and are required to be independent with respect to the Company in accordance with the U.S. federal securities laws and the applicable rules and regulations of the Securities and Exchange Commission and the PCAOB.

We conducted our audits in accordance with the standards of the PCAOB. Those standards require that we plan and perform the audit to obtain reasonable assurance about whether the financial statements are free of material misstatement, whether due to error or fraud. Our audits included performing procedures to assess the risks of material misstatement of the financial statements, whether due to error or fraud, and performing procedures that respond to those risks. Such procedures included examining, on a test basis, evidence regarding the amounts and disclosures in the financial statements. Our audits also included evaluating the accounting principles used and significant estimates made by management, as well as evaluating the overall presentation of the financial statements. We believe that our audits provide a reasonable basis for our opinion.

/s/ Ernst & Young LLP

We served as the Company's auditor from 2008 to 2025.

Boston, Massachusetts
March 17, 2025

QUANTERIX CORPORATION
CONSOLIDATED BALANCE SHEETS
(amounts in thousands, except per share data)

	December 31, 2025	December 31, 2024
ASSETS		
Current assets:		
Cash and cash equivalents	\$ 29,839	\$ 56,709
Marketable securities	88,393	232,413
Accounts receivable, net of allowance for expected credit losses	29,972	32,141
Inventory	54,763	32,775
Prepaid expenses and other current assets	9,290	9,556
Total current assets	212,257	363,594
Restricted cash	3,341	2,610
Property and equipment, net	23,672	17,150
Intangible assets, net	131,787	4,031
Goodwill	26,376	—
Operating lease right-of-use assets	16,664	16,339
Other non-current assets	4,669	2,809
Total assets	\$ 418,766	\$ 406,533
LIABILITIES AND STOCKHOLDERS' EQUITY		
Current liabilities:		
Accounts payable	\$ 13,568	\$ 6,953
Accrued compensation and benefits	14,979	12,620
Accrued expenses and other current liabilities	17,571	8,851
Deferred revenue	20,728	8,827
Operating lease liabilities	7,916	4,756
Total current liabilities	74,762	42,007
Deferred revenue, net of current portion	5,830	1,073
Operating lease liabilities, net of current portion	29,323	32,615
Non-current portion of contingent liabilities	5,024	—
Other non-current liabilities	8,097	800
Total liabilities	123,036	76,495
Commitments and contingencies (Note 16)		
Stockholders' equity:		
Common stock, \$0.001 par value per share:		
Authorized: 120,000 shares; issued and outstanding: 46,744 and 38,573 shares at December 31, 2025 and 2024, respectively	47	39
Additional paid-in capital	873,637	803,160
Accumulated other comprehensive loss	(723)	(3,080)
Accumulated deficit	(577,231)	(470,081)
Total stockholders' equity	295,730	330,038
Total liabilities and stockholders' equity	\$ 418,766	\$ 406,533

The accompanying notes are an integral part of these Consolidated Financial Statements.

QUANTERIX CORPORATION
CONSOLIDATED STATEMENTS OF OPERATIONS
(amounts in thousands, except per share data)

	Year Ended December 31,		
	2025	2024	2023
Revenues:			
Product revenue	\$ 92,941	\$ 79,740	\$ 79,670
Service and other revenue	44,212	51,244	40,089
Collaboration and license revenue	1,501	4,452	1,380
Grant revenue	243	1,985	1,229
Total revenues	138,897	137,421	122,368
Costs of goods sold and services:			
Cost of product revenue	50,981	33,304	29,103
Cost of service and other revenue	22,957	21,013	19,041
Total costs of goods sold and services	73,938	54,317	48,144
Gross profit	64,959	83,104	74,224
Operating expenses:			
Research and development	35,922	31,082	26,064
Selling, general and administrative	138,008	101,618	89,111
Other lease costs	844	3,020	3,712
Impairment and restructuring	15,727	—	1,328
Total operating expenses	190,501	135,720	120,215
Loss from operations	(125,542)	(52,616)	(45,991)
Interest income	8,567	14,655	15,839
Change in fair value of contingent liabilities	4,547	—	—
Other income (expense), net	157	(136)	2,517
Loss before income taxes	(112,271)	(38,097)	(27,635)
Income tax benefit (expense)	5,121	(434)	(719)
Net loss	\$ (107,150)	\$ (38,531)	\$ (28,354)
Net loss per common share, basic and diluted	\$ (2.51)	\$ (1.00)	\$ (0.75)
Weighted-average common shares outstanding, basic and diluted	42,639	38,367	37,594

The accompanying notes are an integral part of these Consolidated Financial Statements.

QUANTERIX CORPORATION
CONSOLIDATED STATEMENTS OF COMPREHENSIVE LOSS
(amounts in thousands)

	Year Ended December 31,		
	2025	2024	2023
Net loss	\$ (107,150)	\$ (38,531)	\$ (28,354)
Other comprehensive income (loss), net of tax:			
Unrealized gains (losses) on marketable securities	(25)	(219)	325
Foreign currency translation	2,382	(1,189)	541
Total other comprehensive income (loss)	2,357	(1,408)	866
Comprehensive loss	\$ (104,793)	\$ (39,939)	\$ (27,488)

The accompanying notes are an integral part of these Consolidated Financial Statements.

QUANTERIX CORPORATION
CONSOLIDATED STATEMENTS OF STOCKHOLDERS' EQUITY
(amounts in thousands)

	Common Stock		Additional paid-in capital	Accumulated other comprehensive income (loss)	Accumulated deficit	Total stockholders' equity
	Shares	Amount				
Balance at December 31, 2022	37,280	\$ 37	\$ 763,629	\$ (2,538)	\$ (403,196)	\$ 357,932
Issuance of common stock under stock plans, including tax effects	734	1	2,690	—	—	2,691
Stock-based compensation expense	—	—	16,823	—	—	16,823
Unrealized gains on marketable securities, net of tax	—	—	—	325	—	325
Foreign currency translation	—	—	—	541	—	541
Net loss	—	—	—	—	(28,354)	(28,354)
Balance at December 31, 2023	38,014	\$ 38	\$ 783,142	\$ (1,672)	\$ (431,550)	\$ 349,958
Issuance of common stock under stock plans, including tax effects	559	1	31	—	—	32
Stock-based compensation expense	—	—	19,987	—	—	19,987
Unrealized losses on marketable securities, net of tax	—	—	—	(219)	—	(219)
Foreign currency translation	—	—	—	(1,189)	—	(1,189)
Net loss	—	—	—	—	(38,531)	(38,531)
Balance at December 31, 2024	38,573	\$ 39	\$ 803,160	\$ (3,080)	\$ (470,081)	\$ 330,038
Issuance of common stock under stock plans, including tax effects	684	1	(134)	—	—	(133)
Common stock issued as consideration in business combinations, net of tax	7,487	7	49,893	—	—	49,900
Stock-based compensation expense	—	—	20,718	—	—	20,718
Unrealized losses on marketable securities, net of tax	—	—	—	(25)	—	(25)
Foreign currency translation	—	—	—	2,382	—	2,382
Net loss	—	—	—	—	(107,150)	(107,150)
Balance at December 31, 2025	46,744	\$ 47	\$ 873,637	\$ (723)	\$ (577,231)	\$ 295,730

The accompanying notes are an integral part of these Consolidated Financial Statements.

QUANTERIX CORPORATION
CONSOLIDATED STATEMENTS OF CASH FLOWS
(amounts in thousands)

	Year Ended December 31,		
	2025	2024	2023
Cash flows from operating activities:			
Net loss	\$ (107,150)	\$ (38,531)	\$ (28,354)
Adjustments to reconcile net loss to net cash used in operating activities:			
Depreciation and amortization expense	15,844	6,463	6,275
Credit losses on accounts receivable	628	588	336
Accretion of marketable securities	(2,078)	(6,833)	(1,964)
Operating lease right-of-use asset amortization	3,081	1,893	2,015
Stock-based compensation expense	20,718	19,987	16,823
Impairment	7,752	—	1,361
Change in fair value of contingent liabilities	(4,547)	—	—
Recognition of off-market liability	(2,726)	—	—
Deferred taxes	(5,867)	(290)	(290)
Other operating activity	(540)	345	140
Changes in assets and liabilities:			
Accounts receivable	10,609	(7,704)	(6,695)
Inventory	4,033	(6,679)	(8,944)
Prepaid expenses and other current assets	3,521	(443)	(2,371)
Other non-current assets	252	(1,215)	(717)
Accounts payable	(528)	723	1,189
Accrued compensation and benefits, accrued expenses, and other current liabilities	(7,914)	1,398	4,410
Net change in other operating assets and liabilities	(12,324)	(4,866)	(2,063)
Net cash used in operating activities	(77,236)	(35,164)	(18,849)
Cash flows from investing activities:			
Purchases of marketable debt securities	(69,757)	(295,606)	(175,613)
Proceeds from sales and maturities of marketable securities	215,829	216,709	31,000
Purchases of property and equipment	(2,612)	(3,368)	(3,841)
Acquisitions, net of cash acquired	(93,229)	—	—
Net cash provided by (used in) investing activities	50,231	(82,265)	(148,454)
Cash flows from financing activities:			
Proceeds from common stock issued under stock plans	738	3,066	2,889
Payments for employee taxes withheld on stock-based compensation awards	(1,446)	(2,610)	(198)
Net cash provided by (used in) financing activities	(708)	456	2,691
Net decrease in cash, cash equivalents, and restricted cash	(27,713)	(116,973)	(164,612)
Effect of exchange rate changes on cash, cash equivalents, and restricted cash	1,574	(734)	301
Cash, cash equivalents, and restricted cash at beginning of period	59,319	177,026	341,337
Cash, cash equivalents, and restricted cash at end of period	\$ 33,180	\$ 59,319	\$ 177,026
Supplemental disclosure of cash flow information:			
Common stock issued in exchange for acquisition of Akoya (Note 3)	\$ 49,900	\$ —	\$ —
Purchases of property and equipment in accounts payable and accruals	\$ 121	\$ 1,230	\$ 419

The accompanying notes are an integral part of these Consolidated Financial Statements.

QUANTERIX CORPORATION
NOTES TO CONSOLIDATED FINANCIAL STATEMENTS

Note 1. Organization and Nature of Business

Quanterix Corporation ("Quanterix" or the "Company") is a life sciences company transforming healthcare innovation by accelerating biomarker breakthroughs from discovery to diagnostics using its ultra-sensitive translational research and spatial biology instruments, consumables, and services. The Company continues to invest in pushing a paradigm shift in healthcare from an emphasis on later-stage treatment to a focus on earlier detection, monitoring, prognosis, and, ultimately, prevention.

Quanterix's proprietary digital "Simoa" detection technology enables customers to reliably detect protein biomarkers at ultra-low concentrations in blood, serum and other fluids that, in many cases, are undetectable using conventional, analog immunoassay technologies. Multi-plexing biomarker analysis in tissue samples with the Company's "Spatial Biology" platforms enables scientists to understand the localized interactions occurring on the cellular level. The Company believes its combination of technologies will enable scientists to help drive diagnostic innovation in the evolving healthcare landscape with data across the tissue to fluid continuum. Currently, the ability of its Simoa platforms to detect proteins in the femtomolar range is enabling the development of novel therapies and diagnostics and has the potential to identify early-stage disease markers before symptoms appear.

The Company sells its proprietary instruments and related consumables worldwide to research laboratories, contract research organizations, academic institutions, and bio-pharmaceutical companies. In addition, the Company provides contract research services and clinical laboratory testing services, including four Laboratory Developed Tests ("LDT"), using its proprietary technology through its Accelerator Laboratory, which is certified under the Clinical Laboratory Improvement Amendments of 1988 ("CLIA") (the "Accelerator Laboratory").

Liquidity

The Company has recognized annual losses from operations since inception and has an accumulated deficit of \$577.2 million as of December 31, 2025. The Company incurred net losses of \$107.2 million, \$38.5 million, and \$28.4 million for the years ended December 31, 2025, 2024, and 2023, respectively. As of December 31, 2025, the Company had cash and cash equivalents of \$29.8 million and marketable securities of \$88.4 million. The Company expects that its current cash, cash equivalents, and marketable securities will be sufficient to fund its operations for a period of at least one year from the date the Consolidated Financial Statements are issued.

Note 2. Significant Accounting Policies

Basis of Presentation

The accompanying Consolidated Financial Statements and Notes to Consolidated Financial Statements have been prepared in accordance with generally accepted accounting principles in the United States of America ("U.S. GAAP") and pursuant to the rules and regulations of the U.S. Securities and Exchange Commission regarding annual financial reporting on Form 10-K.

The Company's fiscal year is the twelve-month period from January 1 through December 31, and all references to "2025," "2024," and "2023," refer to the fiscal year unless otherwise noted. Certain amounts in the prior years' Consolidated Financial Statements have been reclassified to conform to the current year's presentation.

Use of Estimates

The preparation of the Consolidated Financial Statements and Notes to Consolidated Financial Statements in conformity with U.S. GAAP requires management to make estimates and assumptions that affect the reported amounts of assets and liabilities at the dates of the Consolidated Financial Statements, and the reported amounts of revenues and expenses during each fiscal year. Such estimates include, but are not limited to, revenue recognition, valuation of inventory, valuation and impairment of goodwill, intangible and other long-lived assets, valuation of acquired assets and assumed liabilities from acquisitions, valuation of contingent liabilities, recoverability of deferred tax assets, and stock-based compensation expense. The Company bases its estimates on historical experience, known trends, worldwide economic conditions, both general and specific to the life sciences industry, and other relevant factors it believes to be

reasonable under the circumstances. On an ongoing basis, management evaluates its estimates and changes in estimates are recorded in the period in which they become known. Actual results could differ from those estimates.

Principles of Consolidation

The Consolidated Financial Statements include the accounts of Quanterix and its wholly-owned subsidiaries. All intercompany transactions have been eliminated in consolidation.

Business Combinations

The Company accounts for business combinations in accordance with the acquisition method of accounting under ASC 805 - *Business Combinations* ("ASC 805"). The acquisition method of accounting requires the Company to record the acquired assets and liabilities, including identifiable intangibles, at their estimated fair values as of the acquisition date, with any excess of the consideration transferred recorded to goodwill. Goodwill is not amortized; however it is required to be tested for impairment at least annually at the reporting unit level.

If the initial accounting for a business combination is incomplete by the end of a reporting period that falls within a measurement period (not to exceed a year from the date of acquisition), the Company records preliminary amounts. During the measurement period, the Company adjusts the preliminary amounts to reflect new information obtained about facts and circumstances that existed as of the acquisition date that, if known, would have affected the measurement of the amounts recognized as of that date. The Company records these adjustments to the preliminary amounts with a corresponding offset to goodwill. Any adjustments identified after the measurement period are recorded in the Consolidated Statements of Operations.

Acquired intangible assets consist of definite and indefinite-lived intangible assets, including developed technology, customer relationships, know-how, and in process research and development ("IPR&D"). Definite-lived intangible assets are amortized on a straight-line basis over each asset's estimated useful life. The Company accounts for IPR&D as an indefinite-lived intangible asset until the underlying project is completed. Upon completion, the IPR&D asset is accounted for as a finite-lived intangible asset.

For customer contracts acquired in an acquisition, the Company reassesses them under ASC 606 - *Revenue from Contracts with Customers* ("ASC 606") as if it had originated the contract, including assessing differences from the Company's revenue policies and estimates. As part of this reassessment, under ASC 805, the Company elected the practical expedient to reflect the aggregate effect of all modifications that occurred before the acquisition date. Differences from these reassessments, if any, are recorded as contract assets or deferred revenue.

Additionally, the Company assesses acquired customer contracts for off-market terms and records an additional asset or liability for favorable or unfavorable terms. Off-market assets or liabilities are recorded at their estimated fair values as of the acquisition date based on the economic returns that could have been achieved by a market participant in a market transaction.

Refer to Note 3 - *Acquisitions* for discussion of the Company's 2025 acquisitions.

Foreign Currency

The functional currency of the Company's subsidiaries is generally their respective local currencies. These subsidiary financial statements are translated into U.S. dollars using the period-end exchange rates for assets and liabilities, average exchange rates during the corresponding period for revenue and expenses, and historical rates for equity. The effects of foreign currency translation adjustments are recorded in accumulated other comprehensive income (loss), a component of stockholders' equity on the Consolidated Balance Sheets.

Revenue from Contracts with Customers

The Company generates the majority of its revenues from contracts with customers and accounts for them pursuant to ASC 606. Refer to the section titled "Grant Revenue" below for discussion the Company's accounting policy for revenue generated from grant awards.

For contracts with customers, the Company recognizes revenue when a customer obtains control of promised products or services, for an amount that reflects the consideration expected to be received in exchange for those products or services. The Company follows the five step revenue model prescribed by ASC 606 to determine revenue recognition: (i) identify the contract(s) with a customer; (ii) identify the performance obligation(s) in the contract; (iii) determine the transaction price, including variable consideration, if any; (iv) allocate the transaction price to the performance obligations in the contract; and (v) recognize revenue when (or as) the entity satisfies a performance obligation. Revenues are presented net of any sales, value added, or similar taxes collected from customers and remitted to the government.

The Company accounts for a contract when it has approval and commitment from both parties, the fees, payment terms, and rights of the parties regarding the products or services to be transferred are identified, the contract has commercial substance, and it is probable that substantially all of the consideration for the products and services expected to be transferred is collectible. The Company applies judgment in determining the customer's ability and intention to pay for services expected to be transferred, which is based on factors including the customer's payment history, management's ability to mitigate exposure to credit risk (for example, requiring payment in advance of the transfer of products or services, or the ability to stop transferring promised products or services in the event a customer fails to pay consideration when due), and experience selling to similarly situated customers.

The Company's contracts may include either a single promise (referred to as a performance obligation) to transfer a product or service, or a combination of multiple promises to transfer products or services. The Company evaluates the existence of multiple promises within its contracts by using judgment to determine if (1) the customer can benefit from each contractual promise on its own or together with readily available resources and (2) the transfer of each contractual promise is separately identifiable from other promises in a contract. When both criteria are met, each promise is accounted for as a separate performance obligation. Additionally, the Company has elected the practical expedient under ASC 606 to account for shipping and handling as an activity to fulfill a promise to transfer a product, and therefore does not evaluate whether shipping and handling activities are promised services to its customers.

Contracts that include rights to additional products or services that are exercisable at a customer's discretion are generally considered options. The Company assesses if these options provide a material right to the customer and if so, the material right is considered a performance obligation. The identification of material rights requires judgment to determine the value of the option to purchase additional products and services in relation to options that may be provided to, and prices paid by, customers in the normal course of business. Material rights are recognized when exercised by a customer or upon expiration of the right.

The Company determines the transaction price of its contracts based on the amount of consideration it expects to be entitled to, which is generally equal to the contract amount. In some cases, contracts contain variable consideration which primarily relates to (1) sales and usage-based royalties related to the license of intellectual property and (2) contracts with minimum purchase commitments. For sales and usage-based royalties, ASC 606 provides an exception to estimating variable consideration. Under this exception, the Company recognizes revenues from sales or usage-based royalty revenue at the later of when the sales or usage occurs or the satisfaction (or partial satisfaction) of the performance obligation to which the royalty has been allocated. All other variable amounts are constrained to the minimum guaranteed contract amount so that a reversal of cumulative revenue does not occur in future periods. Once there is no longer uncertainty over a variable amount, any incremental fees the Company is entitled to are allocated to the related performance obligation(s).

For contracts that contain multiple performance obligations, the Company allocates the transaction price among the performance obligations on a relative basis according to their standalone selling prices ("SSP"). Determining the SSP for performance obligations requires judgment and is based on factors including prices charged to customers in observable transactions, internal pricing objectives and list prices, pricing of similar products, expected costs to manufacture the Company's products, and estimated margins. The Company has more than one range of standalone selling prices for certain products and services based on the geographic location of customers and sales channel.

Product Revenue

The Company's product revenues are composed of instruments, assay kits, replacement parts, instrument installation fees, and other consumables such as reagents and antibodies. Products are sold directly to customers and are also sold through distributors in EMEA and Asia Pacific regions. Direct instrument sales include installation and an initial year service-type warranty. The Company has determined that the instrument and installation are a combined performance obligation in cases where a customer has a contractual acceptance right that applies through the installation of the instrument. The included service-type warranty is considered a separate performance obligation since a customer could

benefit from it independently with readily available resources and is capable of being sold on its own (with such warranty recorded in services and other revenue on the Consolidated Statements of Operations). Sales of instruments to distributors include a license to import and resell the instruments. The Company has determined these distributor licenses are part of a combined performance obligation with the instrument as the distributor only benefits from the combination of the instrument and ability to resell it.

Instrument sales may also be bundled with assays and other consumables, training, and/or an extended service warranty, each of which is considered a separate performance obligation.

Product revenues for direct instrument sales to customers are recognized when the related performance obligation(s) are satisfied (i.e. when transfer of control of the instrument passes to the customer). For direct instrument sales that require installation prior to contractual acceptance, the combined performance obligation is recognized upon completion of the instrument's installation. For direct instrument sales that do not contain contractual acceptance and instrument sales to distributors, the instrument performance obligation is recognized based on the agreed upon shipping terms (either upon shipment or delivery) as that is when control of the product passes to the customer. When installation is a separate performance obligation, it is recognized upon completion.

Service Revenue

Service revenues are composed of contract research services, initial year of service-type warranties, extended services warranty contracts, repair services, and other services such as training. Contract research services are provided through the Company's Accelerator Laboratory and generally consist of fixed fee contracts.

Revenues from contract research services are generally recognized over time, using an output method based on the number of completed test results, since the work performed does not have an alternative use and the contracts allow for the collection of transaction consideration for costs incurred plus a reasonable margin through the period of performance. For performance obligations that are not recognized over time, they are recognized at the point in time when the Company completes and delivers its research results on each individually completed study. Revenues from warranties and service contracts are recognized ratably over the contract service period since the customer simultaneously receives and benefits from the services.

Collaboration and License Revenue

Collaboration and license revenues are composed of revenue associated with licensing the Company's technology, intellectual property, and know-how associated with the Company's instruments to third parties and for related services. License arrangements consist of sales or usage-based fees and/or future royalties. Revenues are recognized at the point in time the license is delivered as the underlying license is considered functional intellectual property since the customer has the right to use it when it is received. Royalty revenues that are sales or usage-based are recognized at the later of when the sales or usage occurs and the satisfaction (or partial satisfaction) of the performance obligation to which the royalty has been allocated.

Contract Assets and Liabilities

Accounts Receivable and Allowance for Credit Losses

Accounts receivable represent rights to consideration in exchange for products or services that have been transferred by the Company, when payment is unconditional and only the passage of time is required before payment is due. Also included in accounts receivable are unbilled amounts resulting from revenue exceeding the amount billed to the customer, where the right to payment is unconditional. If the right to payment for services performed was conditional on something other than the passage of time, the unbilled amount would be recorded as a separate contract asset.

Generally, the Company's contracts are non-cancellable. For contracts that are cancellable by the customer, the Company records accounts receivable only up to the amount of revenue recognized but not yet collected.

The Company's payment terms vary by customer type and location and the products or services offered. Payment from customers is generally required 30 to 90 days from date of satisfaction of the performance obligation. Accounts receivable do not bear interest and the Company does not provide financing arrangements to its customers.

The Company is exposed to credit losses primarily through accounts receivable from sales of its products and services. The Company's expected loss allowance methodology is developed using historical collection experience, current and future economic and market conditions, a review of the current status of customers' accounts receivable, and management's ability to mitigate exposure to credit risk. Specific allowance amounts are established to record the appropriate provision for customers that have a higher probability of default. The Company monitors changes to the receivables balance on a timely basis, and balances are written off as they are determined to be uncollectible after collection efforts have been exhausted.

Deferred Revenue

The Company refers to contract liabilities as deferred revenue on the Consolidated Balance Sheets. Deferred revenue consists of billings in excess of revenue recognized, primarily related to upfront payments for instruments, consumables, service warranties, and laboratory services.

Costs to Obtain Contracts

The Company capitalizes commissions paid to its sales representatives and related fringe benefit costs that are incremental to obtaining customer contracts. Capitalized commissions are recorded in prepaid expenses and other current assets and other non-current assets on the Consolidated Balance Sheets. These commissions are amortized over the life of the contract and are recorded in cost of goods sold and selling, general and administrative expense on the Consolidated Statements of Operations. The Company has elected the practical expedient allowing commissions with an amortization period of one year or less to be expensed as incurred.

Commissions associated with instrument and consumables sales are earned when the related revenue is recognized. Commissions associated contract research services, warranty, and extended service contracts are earned upon booking.

The Company evaluates potential impairment of these amounts at each balance sheet date, and no related impairments were recorded during the years ended December 31, 2025, 2024, and 2023.

Warranties

For instruments sold directly to customers, the Company provides an initial year of service-type warranty and also sells extended warranty contracts for additional periods beyond the initial year. The service-type warranty guarantees that the Company's instruments are free from material defects in workmanship and materials, excluding normal wear and tear, and includes maintenance services. Under ASC 606, the Company concluded that the initial year service-type warranty is a separate performance obligation since a customer could benefit from it independently with readily available resources and it is capable of being sold on its own. The Company recognizes revenue for these service-type warranties ratably over the service period.

For products sold to distributors, the Company provides an initial year of assurance-type warranty which guarantees that the products conform to the Company's published specifications. Assurance-type warranties do not create a separate performance obligation under ASC 606. In accordance with ASC Topic 460 – *Guarantees*, the Company establishes an accrual for estimated assurance-type warranty expense and records the expense in cost of product revenue on the Consolidated Statements of Operations. Amounts recognized for assurance-type warranties during the years ended December 31, 2025, 2024, and 2023 were not material.

Grant Revenue

Accounting for grants does not fall under ASC 606, as the grantor will not benefit directly from the Company's expansion or product development, and no products or services are transferred to the grantor. As there is no authoritative guidance under U.S. GAAP on accounting for grants to for profit business entities from government entities, the Company accounts for grants by analogy to International Accounting Standards Topic 20, *Accounting for Government Grants and Disclosure of Government Assistance* ("IAS 20") and ASC Topic 958, *Not for-Profit Entities* ("ASC 958"). The decision to account for a grant under IAS 20 or ASC 958 is based on whether the grantor is a government entity.

The Company recognizes grant revenue as the Company performs services under the arrangement when the funding is committed and as each grant's activities are performed. The timing of revenue recognition and receipt of funding

varies by grant and can be independent from performance of the related activities, such as an upfront payment of the award value, or subsequent to the Company's requests for reimbursement for already performed activities (subject to the approval of the granting organization).

Cost of Goods Sold and Services

Cost of Product Revenue

Cost of product revenue consists of manufacturing and assembly costs for instruments, related reagents, other consumables, contract manufacturer costs, personnel costs, allocated overhead, and other direct costs related to product sales. Raw material part costs include inbound shipping and handling costs associated with purchased goods, contract manufacturer costs, personnel costs, overhead and other direct costs related to product sales. Additionally included in cost of product revenue is amortization expense for intangibles that relate to the sale of products and royalty fees due to third parties from revenue generated by collaboration or license deals.

Cost of Service and Other Revenue

Cost of services and other revenue consists of direct costs associated with operating the Company's Accelerator Laboratory on behalf of its customers, including raw materials, personnel costs, allocated overhead costs that include facility and other related costs, and other direct costs. Additional costs include costs related to warranties, and other costs of servicing equipment at customer sites.

Research and Development Expenses

Research and development expense consists primarily of personnel costs, research supplies, third-party development costs for new products, materials for prototypes, quality assurance, amortization expense for intangibles that relate to the development of products, and allocated overhead costs that include facility and other related costs.

Selling, General and Administrative Expenses

Selling, general and administrative expense consists of personnel costs for our sales and marketing, finance, legal, human resources, and general management teams, shipping and handling for product sales, acquisition related costs, other general and administrative costs, as well as professional services costs, such as marketing, advertising, legal and accounting services, and allocated overhead costs that include facility and other related costs. The classification of shipping and handling costs for product sales varies from company to company with some companies recording these as selling, general and administrative expenses and others recording such expenses within costs of goods sold for products. To the extent the classification of the Company's shipping and handling costs differs from the reporting approach used by other companies, the Company's gross margins may not be comparable with those reported by such other companies.

Net Loss Per Share

Basic net loss per common share attributable to common stockholders is calculated by dividing the loss attributable to common stockholders by the weighted-average number of common shares outstanding. Diluted net loss per common share attributable to common stockholders is calculated under the treasury stock method by dividing the loss attributable to common stockholders by the diluted weighted-average number of common shares outstanding. Diluted weighted-average shares outstanding reflect the dilutive effect, if any, of potential common shares issued, such as unvested common stock, unvested restricted stock units ("RSUs"), common stock options, and shares estimated to be purchased under the Company's employee stock purchase plan ("ESPP"). During periods when the Company is in a net loss position, these potential common shares are excluded from the diluted net loss per common share attributable to common stockholders because their effect would be anti-dilutive. Accordingly, basic and diluted net loss per common share attributable to common stockholders were the same for all periods presented.

Cash and Cash Equivalents

Cash and cash equivalents consist of cash deposits and short-term, highly liquid marketable securities that are readily convertible into cash with original maturities of three months or less at the time of purchase. Cash and cash equivalents consist of the following (in thousands):

	As of December 31,	
	2025	2024
Cash	\$ 12,620	\$ 12,283
Money market funds	17,219	44,426
Total cash and cash equivalents	\$ 29,839	\$ 56,709

Restricted Cash

The following table summarizes the period ending cash and cash equivalents as presented on the Consolidated Balance Sheets and the total cash, cash equivalents, and restricted cash as presented on the Consolidated Statements of Cash Flows (in thousands):

	As of December 31,	
	2025	2024
Cash and cash equivalents	\$ 29,839	\$ 56,709
Restricted cash	3,341	2,610
Cash, cash equivalents, and restricted cash	\$ 33,180	\$ 59,319

Restricted cash consists of collateral for letters of credit issued as security for several of the Company's leased facilities and to secure the Company's corporate credit card program. The short-term or long-term classification is determined in accordance with the expiration of the underlying letter of credit and security.

Marketable Securities

The Company's current portfolio of marketable securities is entirely debt securities and may at any time include commercial paper, U.S. Treasuries, corporate notes and bonds, U.S. Government agency bonds, certificates of deposit, and similar types of debt securities. Marketable debt securities with original maturities of three months or less at the time of purchase are recorded in cash and cash equivalents on the Consolidated Balance Sheets as they are considered highly liquid and readily convertible into cash. All other marketable securities, including those with maturities beyond one year, are recorded as current assets on the Consolidated Balance Sheets based on their highly liquid nature and because such securities are available for use in current operations.

The Company classifies its marketable securities as either held to maturity, available-for-sale, or trading at the time of purchase and re-evaluates such classification at each balance sheet date. All of the Company's marketable securities are currently classified as available-for-sale as it may use them in current operations. Available-for-sale securities are recorded at fair value (refer to Note 8 - *Fair Value of Financial Instruments*).

Unrealized gains and losses (other than impairment or credit related losses) are recorded in accumulated other comprehensive income (loss), a component of stockholders' equity on the Consolidated Balance Sheets. Realized gains and losses are determined using the specific identification method and are recorded in other income (expense), net on the Consolidated Statements of Operations.

At least quarterly, the Company monitors its marketable securities for impairment. In the event a security's fair value is less than its amortized cost basis, the Company evaluates whether an impairment exists and if the impairment is a result of credit loss or other factors. For a security in an unrealized loss position, if the Company intends to sell the security in an unrealized loss position, or it is more likely than not that the Company will be required to sell the security before recovery of its amortized cost basis, an impairment loss equal to the difference between the security's fair value and amortized cost basis is recorded in other income (expense), net on the Consolidated Statement of Operations. Additionally, the Company determines if a credit loss exists by considering information about the collectability of the security, current market conditions, and the issuer's financial condition. If a decline in fair value is a result of a credit loss, an allowance for credit losses is recorded in other income (expense), net on the Consolidated Statement of Operations, limited to the portion attributed to the credit loss.

The Company presents accrued interest receivable separately from its marketable securities balance and has elected the practical expedient to exclude such amounts from the assessment and measurement of impairment of its marketable securities. Such accrued interest is recorded in prepaid expenses and other current assets on the Consolidated Balance Sheets.

Inventory

Inventory consists of instruments, assays, and the materials required to manufacture instruments and assays.

Inventory is stated at the lower of cost and net realizable value on a first-in, first-out ("FIFO") basis or average cost basis and includes the cost of materials, labor, and manufacturing overhead. The Company analyzes its inventory levels on each reporting date for slow-moving, excess, and obsolete inventory, and inventory expected to expire prior to being used. These analyses require judgment and are based on factors including, but not limited to, recent historical activity, anticipated or forecasted demand for the Company's products and scientific data supporting the estimated life of materials that expire. If the Company identifies adverse conditions exist, such as unfavorable changes in estimated customer demand, the lives of materials that expire, or actual market conditions that may differ from its projections, the carrying value of the inventory is reduced to its estimated net realizable value.

Property and Equipment

Property and equipment, including leasehold improvements, are stated at cost, net of accumulated depreciation. These assets are depreciated over their estimated useful lives using the straight-line method. Expenditures for maintenance and repairs are charged to expense as incurred, whereas significant expenditures that extend the useful lives of existing assets are capitalized as additions to property and equipment.

Depreciation is calculated based upon the following estimated useful lives:

	Estimated Useful Life
Laboratory and manufacturing equipment	5 years
Demo inventory	3 years
Furniture and fixtures	7 years
Computers and software	3 years
Leasehold improvements	Shorter of asset's life or remaining lease term

The Company develops and modifies internal use software supporting the Company's operations. Certain costs incurred during the application development stage including external direct costs of services used in the development or internal personnel costs for employees directly associated with the development, are capitalized and are recorded in property and equipment on the Consolidated Balance Sheets. The Company begins depreciating these costs over the life of the related asset upon completion of a working model or when it is ready for its intended use. Costs incurred during the preliminary project stage and post-configuration stages are expensed as incurred.

The Company also develops software related to the operation of some of its instruments, which is accounted for under ASC 985 - *Costs of Software to Be Sold, Leased, or Marketed*. The Company capitalizes certain costs incurred after the software reaches technological feasibility and amortizes it once the software is available to customers. The Company did not incur any such costs during the year ended December 31, 2025.

Cloud Computing Costs

For cloud computing arrangements that include a software license or software-as-a-service contracts, certain implementation costs incurred are capitalized and recorded in prepaid expenses and other current assets and other non-current assets on the Consolidated Balance Sheets. Once the software is ready for its intended use, the capitalized costs are amortized over the noncancellable term of the cloud computing service contract, plus any optional renewal periods that are reasonably certain to be exercised.

Leases

The Company primarily enters into operating leases for office, laboratory, manufacturing spaces, and office equipment, and determines whether an arrangement is a lease at inception of the arrangement. The Company accounts for a lease when it has the right to control the leased asset for a period of time, while obtaining substantially all of the assets' economic benefits. Leases are recorded on the Consolidated Balance Sheets as operating lease right-of-use ("ROU") assets and current and non-current operating lease liabilities.

Operating lease ROU assets and lease liabilities are recognized at the lease commencement date based on the estimated present value of the future minimum lease payments over the lease term and any initial direct costs incurred. Initial direct costs are incremental costs of a lease that would not have been incurred had the lease not been executed. The discount rate used to determine the present value of the lease payments is the Company's incremental borrowing rate on a collateralized basis for a similar term and amount, as generally an implicit rate in the lease is not readily determinable. To estimate its incremental borrowing rate, a credit rating applicable to the Company is estimated using a synthetic credit rating analysis since the Company does not currently have an agency-based credit rating.

The Company's lease agreements can contain lease and non-lease components. The Company accounts for the lease and fixed payments for non-lease components as a single lease component under ASC 842 – *Leases*, which increases the amount of the ROU assets and lease liabilities. Most of the Company's lease agreements also contain variable payments, primarily maintenance, utility, and other-related costs, which are not included in the measurement of the ROU assets and lease liabilities and are expensed as incurred.

Some of the Company's leases contain options to extend or terminate the lease. When determining the lease term, these options are included in the measurement and recognition of the Company's ROU assets and lease liabilities when it is reasonably certain that the Company will exercise the option(s). The Company considers various economic factors when making this determination, including, but not limited to, the significance of leasehold improvements incurred in the leased space, the difficulty in replacing the asset, underlying contractual obligations, and specific characteristics unique to a particular lease. Subsequent to entering into a lease arrangement, the Company reassesses the certainty of exercising options to extend or terminate a lease. When it becomes reasonably certain that the Company will exercise an option that was not included in the lease term, the Company accounts for the change in circumstances as a lease modification, which results in the remeasurement of the ROU asset and lease liability as of the modification date.

Leases with a term of 12 months or less upon commencement are not recorded on the Consolidated Balance Sheets and are recorded to expense on a straight-line basis over the lease term.

Impairment of Goodwill and Indefinite-Lived Intangible Assets

The Company assesses goodwill for impairment at the reporting unit level at least annually or whenever events or circumstances change that would more likely than not reduce the fair value of a reporting unit below its carrying amount. Absent an event that indicates a specific impairment may exist, the Company has selected October 1 as the date for performing its annual goodwill impairment test.

A reporting unit is defined as an operating segment or a component of an operating segment to the extent discrete financial information is available that is reviewed by segment management. Prior to the third quarter of 2025, the Company determined it had one reporting unit. After the acquisition of Akoya, the Company determined it had two reporting units, consisting of the legacy Quanterix business and the Akoya business.

An impairment assessment requires evaluating a potential impairment using either a qualitative assessment, to determine if it is more likely than not that the fair value of any reporting unit is less than its carrying amount, or a quantitative analysis, to determine and compare the fair value of each reporting unit to its carrying value, or a combination of both. Judgment is required in determining the use of a qualitative or quantitative assessment, as well as in determining a reporting unit's estimated fair value, as it requires the Company to make estimates of market conditions and operational performance items such as projected financial results, discount rates, control premiums, or valuation multiples for key financial metrics.

During the second quarter of 2025, the Company recorded an impairment charge on its goodwill balance related to the acquisition of Emission. Refer to Note 4 - *Goodwill & Intangible Assets* for further discussion. The Company had no goodwill recorded at December 31, 2024.

At least annually or whenever events or circumstances change, the Company assesses whether IPR&D has been abandoned, in which case it would be written off, or if its estimated fair value is below its carrying value, in which case it could be impaired.

Impairment of Long-Lived Assets

The Company reviews its long-lived assets, including definite lived intangible assets, right of use assets, and property and equipment, for impairment whenever events or circumstances indicate the carrying amount of the asset(s) may not be fully recoverable or that the estimated useful lives may warrant revision. On an ongoing basis, the Company assesses the determination of its asset groups, which primarily focuses on changes in the Company's operating structure, the way in which it expects to deploy its assets, or how the Company intends to recover the cost of its assets.

To assess the recoverability of a long-lived asset or asset group, the Company compares the estimated undiscounted future cash flows for the estimated remaining useful life, or estimated lease term, of the asset (or the primary asset in the asset group) to its carrying value. If the undiscounted cash flows are less than the carrying value, the Company estimates the asset's fair value using the future discounted cash flows associated with the use of the asset. To the extent that the discounted cash flows are less than the carrying value, the asset(s) are impaired and written down to their estimated fair value.

Significant judgment is required to estimate future cash flows, including, but not limited to, estimates about future revenues, expenses, asset disposal value, expected uses of the asset (group), historical customer retention rates, technology roadmaps, customer awareness, trademark and trade name history, contractual provisions that could limit or extend an asset's useful life, market data, discount rates, and potential sublease opportunities including rent and rent escalation rates, time to sublease, and free rent periods.

Refer to Note 15 - *Leases* for a discussion of operating lease impairment charges recorded in 2025.

Fair Value of Financial Instruments

The carrying amount reflected on the Consolidated Balance Sheets for cash, restricted cash, accounts receivable, prepaid expenses and other current assets, accounts payable, accrued compensation and benefits, and accrued expenses and other current liabilities approximate their fair values due to their short-term nature.

Additionally, the Company has certain financial instruments that are required to be measured at fair value on a recurring basis including cash equivalents, marketable securities, and contingent liabilities. The fair values of these financial instruments are classified as Level 1, 2, or 3 within the fair value hierarchy as follows:

- Level 1: Observable inputs based on unadjusted quoted prices in active markets for identical assets.
- Level 2: Inputs, other than Level 1 inputs, that are observable either directly or indirectly, such as quoted prices for similar assets, quoted prices in markets that are not active, other inputs that are observable or can be corroborated by observable market data for substantially the full term of the assets.
- Level 3: Unobservable inputs for which there is little or no market data and such inputs are significant to the fair value of the assets. These inputs reflect the Company's assumptions about the inputs that market participants would use in pricing the asset.

To the extent that the valuation is based on models or inputs that are less observable or unobservable in the market, the determination of fair value requires more judgment.

Income Taxes

The Company recognizes deferred tax assets and liabilities for the expected future tax consequences of events that have been recognized in the Consolidated Financial Statements or tax returns. Deferred tax assets and liabilities are determined based on differences between the carrying amount and the tax basis of the assets and liabilities using the enacted tax rates in effect in the years in which the differences are expected to reverse. A valuation allowance against deferred tax assets is recorded if, based on the weight of the available evidence, it is more likely than not that some or all of

the deferred tax assets will not be realized. On an ongoing basis, the Company reassess the valuation allowance on its deferred tax assets, weighing positive and negative evidence to assess their recoverability.

The Company accounts for uncertain tax positions using a more likely than not threshold for recognizing uncertain tax positions, in accordance with ASC 740 – *Income Taxes*. The evaluation of uncertain tax positions is based on factors that include, but are not limited to, changes in tax law, the measurement of tax positions taken or expected to be taken in tax returns, the effective settlement of matters subject to audit, new audit activity, and changes in facts or circumstances related to a tax position. The Company evaluates uncertain tax positions on an ongoing basis and adjusts any liability recorded to reflect subsequent changes in the relevant facts surrounding the uncertain positions. Amounts recorded for uncertain tax positions, including interest and penalties, are recorded in income tax benefit (expense) on the Consolidated Statement of Operations.

Credit, Product, and Supplier Concentrations, and Off-Balance-Sheet Risk

Financial instruments that potentially expose the Company to concentrations of credit risk primarily consist of cash, cash equivalents, marketable securities, and accounts receivable. The Company limits its risk exposure by having its cash, cash equivalents, and marketable securities held at large commercial banks and investing in highly rated marketable securities. Maximum exposure to losses related to cash, cash equivalents, marketable securities, and accounts receivable are limited to their carrying amounts.

As of both December 31, 2025 and 2024, no customer accounted for greater than 10% of the Company's gross accounts receivable. Customers outside the United States represented 42% and 40% of the Company's gross accounts receivable balance as of December 31, 2025 and 2024, respectively.

For the year ended December 31, 2025 and 2024, no customers accounted for greater than 10% of the Company's total revenue. For the year ended December 31, 2023, one customer of approximately \$14.0 million of revenue accounted for greater than 10% of the Company's total revenue.

The Company is also subject to supply chain risks related to outsourced manufacturing for most of its instruments. Although there are a limited number of manufacturers for its instruments, the Company believes that other suppliers could provide similar manufacturing on comparable terms. A change in suppliers, however, could cause a delay in manufacturing and a possible loss of sales, which would adversely affect operating results. In addition to outsourced manufacturing of some of its instruments, the Company also purchases antibodies through a number of different suppliers. Although a disruption in service from any one of its antibody suppliers is possible, the Company believes that it would be able to find an adequate supply from alternative suppliers.

Stock-Based Compensation

The Company measures and recognizes stock-based compensation expense by calculating the estimated fair value of stock options, RSUs, or purchase rights issued under the Company's ESPP on the grant date. The Company generally issues new common shares upon exercise of options and vesting of RSUs. Awards granted by the Company are routine in nature including new hire, annual, and promotion grants.

The fair value of stock options and purchase rights under the ESPP is estimated using the Black-Scholes option-pricing model. The Black-Scholes model requires the Company to make assumptions about the expected or contractual term of the option or purchase right, the expected volatility, risk-free interest rates, and expected dividend yield. The Company estimates the expected term of options granted to employees utilizing historical exercise data. The expected term is applied to the stock option grant group as a whole, as the Company does not expect substantially different exercise or post-vesting termination behavior among its employee population. The expected volatility is based on the Company's historical volatility. The risk-free interest rate is based on the U.S. Treasury yield curve in effect at the time of grant, commensurate with the expected term. The expected dividend yield is zero as the Company has never paid dividends and has no current plans to pay any dividends on common stock.

The fair value of RSUs is determined using the closing market price of the Company's common stock on the grant date.

The Company recognizes stock-based compensation expense on a straight-line basis over an award's requisite service period, which is the vesting period for stock options and RSUs, and the offering period for purchase rights under the ESPP. The Company recognizes forfeitures as they occur.

Restructuring Costs

The Company records restructuring charges when a restructuring plan is approved and the amounts to be incurred are estimable. Restructuring costs are comprised of employee separation costs, primarily severance and related benefit payments, and any associated costs related to implementing a restructuring plan to reorganize operations.

Refer to Note 19 - *Restructuring* for further discussion on the restructuring plan implemented during the year ended December 31, 2025.

Recent Accounting Pronouncements

Recently Adopted Accounting Standards

In December 2023, the FASB issued ASC Update No. 2023-09, *Income Taxes (Topic 740): Improvements to Income Tax Disclosures* ("ASU 2023-09"). This update enhances income tax disclosure requirements by requiring public entities to provide additional information in its tax rate reconciliation and additional disclosures about income taxes paid. The new standard became effective for the Company's annual financial statements for the period beginning on January 1, 2025. The Company adopted this standard as of January 1, 2025 on a prospective basis and the required disclosures are included in Note 14 - *Income Taxes*.

Recent Accounting Standards To Be Adopted

In December 2025, the FASB issued ASC Update No. 2025-12, *Codification Improvements*. This update provides a variety of language changes and clarity across several topics which are applicable to the Company. The amendments in this update may be applied prospective or retroactively. Additionally, the Company is permitted to elect the transition method for these updates on an issue-by-issue basis. The new standard will be effective for the Company for annual reporting periods beginning after December 15, 2026, and interim reporting periods within those annual reporting periods. The Company is currently evaluating the impact of adoption of the standard on its Consolidated Financial Statements and related disclosures.

In December 2025, the FASB issued ASC Update No. 2025-11, *Interim Reporting (Topic 270): Narrow-Scope Improvements*. This update enhances disclosure of an entity's interim disclosure requirements. The amendments in this update can be applied prospectively or retrospectively. The new standard will be effective for the Company for interim reporting periods within annual reporting periods beginning after December 15, 2027. The Company is currently evaluating the impact of adoption of the standard on its Consolidated Financial Statements and related disclosures.

In November 2025, the FASB issued ASC Update No. 2025-10, *Government Grants (Topic 832): Accounting for Government Grants Received by Business Entities*. This update establishes guidance for recognition, measurement and presentation of government grants received by public business entities. The new standard may be applied using a modified prospective approach, modified retrospective approach, or retrospective approach. The new standard will be effective for the Company for annual reporting periods beginning after December 15, 2028 and interim reporting periods within those annual reporting periods. The Company is currently evaluating the impact of adoption of the standard on its Consolidated Financial Statements and related disclosures.

In September 2025, the FASB issued ASC No. 2025-06, *Goodwill and Other—Internal-Use Software (Subtopic 350-40): Targeted Improvements to the Accounting for Internal-Use Software*. This update enhances disclosure of an entity's internal use software by removing prescriptive and sequential software development stages. The amendments in this update can be applied prospectively or retrospectively. The new standard will be effective for the Company for annual reporting periods beginning after December 15, 2027 and interim reporting periods within those annual reporting periods beginning after December 15, 2027. The Company is currently evaluating the impact of adoption of the standard on its Consolidated Financial Statements and related disclosures.

In July 2025, the FASB issued ASU No. 2025-05, Financial Instruments – Credit Losses (Topic 326): *Measurement of Credit Losses for Accounts Receivable and Contract Assets*. This update provides a practical expedient to assume that current conditions as of the balance sheet date will persist through a reasonable and supportable forecast period for eligible assets when estimating expected credit losses for current accounts receivable and current contract assets arising from transactions accounted for under ASC 606. The update also allows an entity to make an accounting policy election to consider subsequent collections of balances received after the balance sheet date through a date selected by the entity. The amendments in this update are required to be applied prospectively. The new standard will be effective for the Company for annual and interim reporting periods beginning after December 15, 2025. The Company does not expect that the adoption of this standard will have a material impact on its Consolidated Financial Statements and related disclosures.

In November 2024, the FASB issued ASC Update No. 2024-03, *Reporting Comprehensive Income (Topic 220): Expense Disaggregation Disclosures*. This update enhances disclosure of an entity's expenses, primarily through additional disaggregation of income statement expenses. The update also requires entities to disclose qualitative descriptions of the amounts remaining in relevant expense captions that are not separately disaggregated quantitatively. The amendments in this update should be applied prospectively, but entities have the option to apply it retrospectively. The new standard will be effective for the Company's annual reporting periods beginning after December 15, 2026, and interim reporting periods within annual reporting periods beginning after December 15, 2027.. The Company is currently evaluating the impact that the adoption of this standard will have on its Consolidated Financial Statements and related disclosures.

Note 3. Acquisitions

Akoya Biosciences, Inc.

On July 8, 2025 (the "Akoya Closing Date"), the Company completed the transactions contemplated by the Amended and Restated Agreement and Plan of Merger dated as of April 28, 2025, by and among the Company, Wellfleet Merger Sub, Inc., a Delaware corporation and wholly owned subsidiary of the Company, and Akoya Biosciences, Inc., a Delaware corporation. On the Akoya Closing Date, Wellfleet Merger Sub, Inc. merged with and into Akoya (the "Merger"), with Akoya surviving the Merger as a wholly owned subsidiary of the Company.

Akoya is a life sciences technology company based in Marlborough, Massachusetts delivering spatial biology solutions focused on transforming discovery, clinical research and diagnostics. The acquisition is part of the Company's plans to establish an ecosystem to identify and measure biomarkers across tissue and blood, expand its technology offerings into oncology and immunology, and expand its portfolio of laboratory service offerings.

Total Consideration Transferred

The following table presents the fair value of the consideration transferred for the Merger as of the Akoya Closing Date (in thousands, except for exchange ratio and stock price):

Total Akoya common stock and equity instruments outstanding as of July 7, 2025 to be converted		51,136
Exchange ratio		0.147
Total shares of Quanterix common stock to be issued		7,517
Quanterix stock price per share as of the Akoya Closing Date	\$	6.54
Fair value of Akoya common stock and equity instruments converted to Quanterix common stock	\$	49,161
Cash consideration paid (1)	\$	18,942
Cash paid for debt extinguishment (2)	\$	82,131
Fair value of replacement equity awards attributable to pre-acquisition service (3)	\$	739
Total fair value of consideration transferred	\$	150,973

(1) Represents cash consideration paid to Akoya stockholders, including fractional shares, of \$0.37 per share of Akoya common stock.

(2) Represents the repayment of Akoya's long-term debt upon closing of the acquisition, including \$7.0 million of early termination, legal, and prepayment fees.

(3) Represents the fair value of certain equity-based awards held by Akoya employees prior to the acquisition date that have been replaced with Quanterix equity-based awards. The portion of these awards that relates to services performed prior to the acquisition date are included within the purchase price.

Upon completion of the Merger, the Company assumed Akoya's stock incentive plans. All Akoya RSUs that were outstanding immediately prior to the completion of the Merger were automatically adjusted by an exchange ratio and converted into a RSU award covering shares of the Company's common stock on the same terms and conditions, including continuing vesting requirements.

Preliminary Allocation of Purchase Price

The following table summarizes the preliminary allocation of the purchase price to the estimated fair values of the assets acquired and liabilities assumed as of the acquisition date (in thousands):

	As of September 30, 2025	Measurement Period Adjustments (1)	As of December 31, 2025
Assets:			
Cash and cash equivalents	\$ 16,108	\$ —	\$ 16,108
Accounts receivable, net of allowance for expected credit losses	8,616	—	8,616
Inventory	25,800	—	25,800
Prepaid expenses and other assets	5,483	—	5,483
Property and equipment, net	12,087	—	12,087
Intangible assets	121,800	—	121,800
Goodwill (2)	23,460	2,916	26,376
Operating lease right-of-use assets	4,585	—	4,585
Finance lease right-of-use assets	1,041	—	1,041
Total assets acquired	218,980	2,916	221,896
Liabilities:			
Accounts payable	8,266	—	8,266
Accrued expenses and other liabilities	34,206	2,916	37,122
Deferred revenue	18,879	—	18,879
Operating lease liabilities	5,616	—	5,616
Finance lease liabilities	1,040	—	1,040
Total liabilities assumed	68,007	2,916	70,923
Net assets acquired	\$ 150,973	\$ —	\$ 150,973

(1) The Company recorded a measurement period adjustment resulting from new information obtained related to an off-market liability and related tax impacts.

(2) Goodwill represents the estimated fair value of the expected synergies from combining Akoya with Quanterix, as well as the value of the acquired workforce. The goodwill is not deductible for income tax purposes and has been fully assigned to the Akoya reporting unit.

The determination of the fair values the assets acquired and liabilities assumed involves significant judgment in selecting inputs used in the valuation methodologies, including, but not limited to, projected revenues and expenses, future changes in technology, estimated selling prices, replacement costs or margins, customer attrition rates, covenants not to compete, obsolescence of developed technologies, the likelihood and timing of achieving milestones or performance targets, discount rates, and assumptions about the period of time a brand will continue to be used. The use of different estimates could produce different results.

The purchase price allocation set forth above is preliminary as the Company continues to obtain information to complete the purchase price allocation.

Intangible Assets

The preliminary fair value of the intangible assets acquired as of the acquisition date was as follows (in thousands):

	Preliminary Fair Value
Definite-lived intangible assets:	
Developed technology	\$ 99,600
Customer relationships	2,900
Total definite-lived intangible assets	102,500
Indefinite-lived intangible assets:	
In process research and development	19,300
Total intangible assets	\$ 121,800

The total weighted-average useful life of acquired, definite-lived intangible assets was 9.6 years at acquisition.

In determining the fair values, management primarily relied on income based approaches using Level 3 inputs. A multi-period excess earnings valuation methodology was used for the developed technology and IPR&D intangible assets, and a distributor method was used for the customer relationships intangible. These income approaches required the use of estimates including: projected revenues and expenses related to the particular asset, obsolescence rates, customer retention rates, discount rates, and certain published or readily available industry benchmark data. In establishing the estimated useful life of each definite-lived intangible asset, the Company relied primarily on the duration of the cash flows utilized in the valuation model.

Off-Market Customer Contract

The Company assessed the unfavorable terms of an acquired contract between Akoya and a customer and recorded a \$16.7 million off-market liability as of the acquisition date. The Company determined the preliminary fair value of the off-market component based on an income approach using Level 3 inputs. This income approach required the use of estimates including: projected revenue, expected profit margin, and a discount rate. The off-market liability is a non-cash balance that will be recognized into revenue as the Company satisfies the associated performance obligation. During the year ended December 31, 2025, the Company recognized \$2.7 million of non-cash revenue from the amortization of the off-market liability.

Financial Results

The operating results of Akoya have been included in the Company's financial statements since the acquisition date. For the year ended December 31, 2025, the acquisition added \$33.7 million of revenue to total revenues and a loss of \$12.0 million to net loss.

The unaudited pro forma financial information presented below was derived from historical financial records of Quanterix and Akoya and presents the operating results for the periods presented as if the acquisition occurred on January 1, 2024 (in thousands):

	Year Ended December 31,	
	2025	2024
Revenues	\$ 174,269	\$ 216,145
Net loss	\$ (135,454)	\$ (96,919)

The unaudited pro forma financial information has been prepared in accordance with U.S. GAAP and includes adjustments primarily to reflect (1) additional amortization expense for the acquired intangible assets, (2) additional amortization of the fair value increases of acquired inventory and property and equipment, (3) elimination of interest expense from the repayment of Akoya's long-term debt, and (4) recognition of additional non-cash revenue from the amortization of an off-market liability. Accordingly, the unaudited pro forma financial information are presented for

informational purposes only and are not necessarily indicative of the results of operations that would have been achieved if the acquisitions had taken place on January 1, 2024, nor are they indicative of future results.

Emission, Inc.

On January 8, 2025 (the "Emission Closing Date"), the Company acquired all of the issued and outstanding shares of capital stock of Emission, Inc. ("Emission"), a life sciences manufacturing company based in Georgetown, Texas. Emission produces large-scale, highly-uniform dye-encapsulating magnetic beads designed for low and mid-plex assays and a mid-plex platform that reads these proprietary beads. The transaction is part of the Company's plans to secure the use of Emission's highly controlled beads in the Company's next generation platforms and expansion into a new multi-plex market segment targeting third-party original equipment manufacturer customers.

Total Consideration Transferred

The following table summarizes the fair value of the aggregate consideration paid or payable for Emission as of the Emission Closing Date (in thousands):

Cash paid (1)	\$	8,997
Holdback (2)		1,000
Contingent consideration (3)		6,612
Total fair value of consideration transferred	\$	<u>16,609</u>

(1) Cash paid represents the contractual amount paid on the Emission Closing Date. Cash acquired was not material.

(2) The holdback is expected to be paid during the first quarter of 2026 and is subject to applicable adjustments.

(3) The acquisition includes contingent consideration discussed in the section titled "Contingent Payments".

Contingent Payments

The Emission transaction included two arrangements that result in additional cash payments to the seller. An additional \$10.0 million was paid in the fourth quarter of 2025 upon completion of certain technical milestones ("Earnout 1") and up to \$50.0 million could be payable based on the amount and timing of certain performance targets over a five year period ending December 31, 2029 ("Earnout 2").

Under ASC 805, the Company determined Earnout 1 was compensation expense and was recognized separately from the business combination. In accordance with ASC 710 - *Compensation*, Earnout 1 was recognized over the period certain technical requirements were transferred and certain milestones were achieved. During the year ended December 31, 2025, the Company recorded \$5.1 million in research and development and \$4.9 million in selling, general and administrative on the Consolidated Statements of Operations related to Earnout 1.

The fair value of Earnout 2 on the acquisition date was \$6.6 million, which represents purchase price and is included in the accounting for the business combination. Monte-Carlo simulations were used to determine the fair value, including the following significant unobservable inputs: projected revenue, a risk adjusted discount rate, and revenue volatility. Refer to Note 8 - *Fair Value of Financial Instruments* for further discussion.

Allocation of Purchase Price

The following table summarizes the allocation of the purchase price to the estimated fair values of the assets acquired and liabilities assumed as of the acquisition date (in thousands):

Assets:	
Cash and cash equivalents	\$ 43
Accounts receivable, net of allowance for expected credit losses	49
Inventory	307
Intangible asset (1)	12,900
Goodwill (2)	6,374
Total assets acquired	19,673
Liabilities:	
Accounts payable	57
Deferred tax liability (3)	3,007
Total liabilities assumed	3,064
Net assets acquired	<u>\$ 16,609</u>

- (1) The acquired intangible asset is finite-lived, represents developed technology, and has an estimated useful life of 14 years. The determination of the fair value of the definite-lived intangible asset required management judgment and the consideration of a number of factors. In determining the fair value, management primarily relied on a multi-period excess earnings valuation methodology. This methodology required the use of estimates including: projected revenues related to the particular asset, its obsolescence rate, royalty, margin, and discount rates, and certain published or readily available industry benchmark data. In establishing the estimated useful life of the acquired intangible asset, the Company relied primarily on the duration of the cash flows utilized in the valuation model.
- (2) Goodwill represents the estimated fair value of the expected synergies from combining Emission with Quanterix, as well as the value of the acquired workforce. The goodwill is not deductible for income tax purposes. During the second quarter of 2025 the Company recorded an impairment charge related to this goodwill (refer to Note 4 - *Goodwill & Intangible Assets*).
- (3) Recorded in other non-current liabilities on the Consolidated Balance Sheets.

The operating results of Emission have been included in the Company's financial statements since the acquisition date and are not material to the Company's consolidated financial results. No measurement period adjustments have been recorded since the Emission Closing Date.

Call Option Agreement

In connection with the closing of the acquisition of Emission, the Company entered into a call option agreement (the "Option Agreement"), in which the Emission selling shareholders have the right to repurchase all of the outstanding capital stock of Emission for \$10.0 million after five years if Emission's revenues do not exceed \$5.0 million in any one year during such five-year period. If the Emission selling shareholders exercise the right to repurchase Emission, the Company will retain a perpetual, fully-paid, irrevocable license to all Emission intellectual property required to continue to manufacture and commercialize the Company's products. The Company determined that the call option is embedded in the purchased shares of Emission and does not require separate accounting unless exercised.

Acquisition Costs

Acquisition costs related to the Company's business combinations were \$12.4 million and \$1.1 million for the year ended December 31, 2025 and 2024, respectively, and are recorded in selling, general and administrative in the Consolidated Statements of Operations.

Note 4. Goodwill & Intangible Assets***Goodwill and Impairment***

During the second quarter of 2025, the Company assessed several events and circumstances, including a larger than expected decline in the Company's revenue and bookings primarily due to the rapidly changing macro-economic conditions resulting from reductions in U.S. federal research funding, reductions in research and development spending by larger pharmaceutical customers, and new import tariffs. As a result, the Company concluded it was more likely than not that the fair value of its then single reporting unit was less than its carrying value and performed a quantitative impairment test for its goodwill. The Company estimated the implied fair value of its reporting unit using a market valuation approach, which included inputs such as the Company's quoted stock price. As a result of the quantitative test, the Company concluded that the goodwill arising from the acquisition of Emission was fully impaired and recorded an impairment charge of \$6.4 million.

At September 30, 2025 and December 31, 2025, the Company performed additional quantitative interim impairment tests as result of continued events and circumstances that indicated its remaining goodwill could be impaired. The Company estimated its reporting units' implied fair values based on an income valuation approach using Level 3 inputs, which used estimates including: projected revenues and expenses, obsolescence rates, customer retention rates, a discount rate, and certain published or readily available industry benchmark data. As a result of these quantitative tests, the Company determined that its goodwill was not impaired as of either date.

The Company performed its annual impairment test as of October 1, 2025 and concluded its goodwill was not impaired. The annual impairment test utilized a qualitative analysis since an interim impairment test had been performed immediately prior to the annual test and no additional events or changes in circumstances occurred between the interim test and annual test.

Changes in the carrying amount of goodwill are as follows (in thousands):

	Total Goodwill
Balance as of December 31, 2024	\$ —
Acquisition of Emission (1)	\$ 6,374
Goodwill impairment	\$ (6,374)
Acquisition of Akoya, including measurement period adjustments (1)	\$ 26,376
Balance as of December 31, 2025	<u>\$ 26,376</u>

(1) Refer to Note 3 - *Acquisitions*.

Intangible Assets, Long-Lived Assets, and Impairment

As a result of the acquisitions of Akoya and Emission, the Company reassessed its asset groups. Primarily due to management's updated expectations of how it plans to deploy and recover the costs of its assets, the Company determined the assets acquired in the Akoya and Emission acquisitions would each become their own asset group, along with vacant leased facilities which are each evaluated as separate asset groups for the purposes of assessing recoverability.

Prior to performing each of its goodwill impairment tests, the Company tested the recoverability of its long-lived assets. The Company utilized an undiscounted cash flow analysis to determine if the cash flows expected to be generated by each of its asset groups over the remaining estimated useful lives of each group's primary asset were sufficient to recover the carrying value of each asset group. Significant assumptions that form the basis of the forecasted results utilized to calculate undiscounted cash flows include: projected revenues and expenses and market conditions related to these assets. The Company also estimated the disposal value of certain asset groups by relying on a market based valuation approach using Level 2 inputs. Collectively, these assumptions and estimates are based on a complex series of judgments about future events and rely heavily on estimates and assumptions that have been deemed reasonable by the Company. Changes in the estimates or assumptions used could materially affect the determination of recoverability. Potential events and circumstances that could have an adverse impact on the Company's estimates and assumptions include, but are not limited to, lower than expected bookings growth, increases in costs, and other macroeconomic factors.

As of December 31, 2025, the Company concluded that none of its intangible or other long-lived assets were impaired, other than a portion of its operating lease right of use assets as described in Note 15 - *Leases*.

Should economic conditions deteriorate further or remain depressed for a prolonged period of time, estimates of future cash flows for each of the Company's asset groups may be insufficient to support their carrying value, requiring an impairment. Impairment charges, if any, may be material to the results of operations and financial position.

Acquired intangible assets consisted of the following (in thousands, except useful life and weighted average life amounts):

	Estimated Useful Life (in years)	As of December 31, 2025				
		Gross Carrying Value	Accumulated Amortization	Cumulative Translation Adjustment	Net Carrying Value	Weighted Average Life Remaining (in years)
Definite-lived intangible assets:						
Developed technology	7.0 - 14.0	\$ 114,150	\$ (7,596)	\$ —	\$ 106,554	9.6
Know-how	8.5	13,000	(8,445)	(1,470)	3,085	2.0
Customer relationships	8.5 - 10.0	4,260	(1,408)	(4)	2,848	8.5
Non-compete agreements	5.5	340	(340)	—	—	—
Trade names	3.0	50	(50)	—	—	—
Total definite-lived intangible assets		131,800	(17,839)	(1,474)	112,487	
Indefinite-lived intangible assets:						
In-process research and development		19,300	—	—	19,300	
Total intangible assets		\$ 151,100	\$ (17,839)	\$ (1,474)	\$ 131,787	

	Estimated Useful Life (in years)	As of December 31, 2024				
		Gross Carrying Value	Accumulated Amortization	Cumulative Translation Adjustment	Net Carrying Value	Weighted Average Life Remaining (in years)
Definite-lived intangible assets:						
Know-how	8.5	\$ 13,000	\$ (7,057)	\$ (2,093)	\$ 3,850	3.0
Developed technology	7.0	1,650	(1,645)	—	5	0.1
Customer relationships	8.5 - 10.0	1,360	(1,166)	(18)	176	3.1
Non-compete agreements	5.5	340	(285)	(55)	—	—
Trade names	3.0	50	(50)	—	—	—
Total intangible assets		\$ 16,400	\$ (10,203)	\$ (2,166)	\$ 4,031	

The Company recorded amortization expense of \$7.6 million, \$1.6 million, and \$1.6 million for the years ended December 31, 2025, 2024, and 2023, respectively. Amortization of know-how and developed technology is recorded in cost of goods sold, and amortization of customer relationships, non-compete agreements, and trade names is recorded in selling, general and administrative on the Consolidated Statements of Operations.

Future estimated amortization expense is as follows (amounts in thousands):

	As of December 31, 2025	
2026	\$	13,258
2027		13,236
2028		11,686
2029		11,656
2030		11,656
Thereafter		50,995
Total amortization expense	\$	<u>112,487</u>

Note 5. Revenue and Related Matters

Revenue from Contracts with Customers

The Company's customers primarily consist of entities engaged in the life sciences research that pursue the discovery and development of new drugs for a variety of neurologic, oncologic, cardiovascular, infectious disease, and other protein biomarkers associated with diseases. The Company's customer base includes pharmaceutical, biotechnology, contract research organizations, academic, and government institutions.

Disaggregated Revenue

When disaggregating revenue, the Company considers all of the economic factors that may affect its revenues. The following tables disaggregate the Company's revenue by geography, based on the location products and services are consumed, and revenue type (in thousands):

	Year Ended December 31, 2025			
	North America	EMEA	Asia Pacific	Total
Product revenue:				
Instruments	\$ 7,039	\$ 5,007	\$ 5,814	\$ 17,860
Consumable and other products	42,049	23,661	9,371	75,081
Total product revenue	<u>\$ 49,088</u>	<u>\$ 28,668</u>	<u>\$ 15,185</u>	<u>\$ 92,941</u>
Service and other revenue:				
Research services	\$ 23,677	\$ 1,799	\$ 505	\$ 25,981
Service-type warranties	6,270	3,844	1,028	11,142
Other services	4,508	2,405	176	7,089
Total service and other revenue	<u>\$ 34,455</u>	<u>\$ 8,048</u>	<u>\$ 1,709</u>	<u>\$ 44,212</u>
Collaboration and license revenue:				
Collaboration and license revenue	\$ 1,571	\$ —	\$ (70)	\$ 1,501
Total collaboration and license revenue	<u>\$ 1,571</u>	<u>\$ —</u>	<u>\$ (70)</u>	<u>\$ 1,501</u>

	Year Ended December 31, 2024			
	North America	EMEA	Asia Pacific	Total
Product revenue:				
Instruments	\$ 3,924	\$ 3,857	\$ 2,699	\$ 10,480
Consumable and other products	39,794	20,760	8,706	69,260
Total product revenue	\$ 43,718	\$ 24,617	\$ 11,405	\$ 79,740
Service and other revenues:				
Research services	\$ 31,252	\$ 5,833	\$ 913	\$ 37,998
Service-type warranties	6,319	3,490	798	10,607
Other services	1,618	976	45	2,639
Total service and other revenue	\$ 39,189	\$ 10,299	\$ 1,756	\$ 51,244
Collaboration and license revenue:				
Total collaboration and license revenue	\$ 3,681	\$ —	\$ 771	\$ 4,452

	Year Ended December 31, 2023			
	North America	EMEA	Asia Pacific	Total
Product revenue:				
Instruments	\$ 6,374	\$ 4,384	\$ 4,947	\$ 15,705
Consumable and other products	35,122	21,216	7,627	63,965
Total product revenue	\$ 41,496	\$ 25,600	\$ 12,574	\$ 79,670
Service and other revenue:				
Research services	\$ 24,706	\$ 2,000	\$ 1,122	\$ 27,828
Service-type warranties	6,265	3,001	613	9,879
Other services	1,436	951	(5)	2,382
Total service and other revenue	\$ 32,407	\$ 5,952	\$ 1,730	\$ 40,089
Collaboration and license revenue:				
Total collaboration and license revenue	\$ 1,380	\$ —	\$ —	\$ 1,380

The following table disaggregates the Company's revenue by technology type (in thousands):

	Year Ended December 31,		
	2025	2024	2023
Product revenue:			
Simoa	\$ 71,034	\$ 79,740	\$ 79,670
Spatial Biology	21,907	—	—
Total product revenue	\$ 92,941	\$ 79,740	\$ 79,670
Service and other revenue:			
Simoa	\$ 32,388	\$ 51,244	\$ 40,089
Spatial Biology	11,824	—	—
Total service and other revenue	\$ 44,212	\$ 51,244	\$ 40,089

All of the Company's collaboration and license revenue was generated by Simoa technology.

Contract Assets

There were no contract assets as of December 31, 2025 or 2024.

Deferred Revenue

During the years ended December 31, 2025 and 2024, the Company recognized \$6.5 million and \$7.9 million of revenue, respectively, related to its deferred revenue balance at January 1 of each such period.

Remaining Performance Obligations

As of December 31, 2025, the aggregate amount of transaction prices allocated to performance obligations that were not yet satisfied, or were partially satisfied, was \$26.6 million. Of this amount \$20.7 million is expected to be recognized as revenue in the next 12 months, with the remaining \$5.8 million expected to be recognized thereafter. The \$5.8 million primarily consists of amounts billed for undelivered services related to extended service-type warranties and research services.

Note 6. Allowance for Credit Losses

The change in the allowance for credit losses on accounts receivable is summarized as follows (in thousands):

	2025	2024	2023
Balance at December 31 of prior year	\$ 1,042	\$ 454	\$ 118
Provision acquired through acquisition	1,015	—	—
Provision for expected credit losses	1,465	712	729
Write-offs and recoveries collected	(1,154)	(124)	(393)
Balance at December 31	<u>\$ 2,368</u>	<u>\$ 1,042</u>	<u>\$ 454</u>

Note 7. Marketable Securities

The amortized cost, gross unrealized gains, gross unrealized losses, and fair value of the Company's marketable securities by major security type were as follows (in thousands):

	As of December 31, 2025			
	Amortized cost	Unrealized Gains	Unrealized Losses	Fair Value
Commercial paper	\$ 12,875	\$ 2	\$ —	\$ 12,877
U.S. Treasuries	29,572	78	—	29,650
U.S. Government agency bonds	13,588	7	(1)	13,594
Corporate bonds	32,279	13	(20)	32,272
Total marketable securities	<u>\$ 88,314</u>	<u>\$ 100</u>	<u>\$ (21)</u>	<u>\$ 88,393</u>

	As of December 31, 2024			
	Amortized Cost	Unrealized Gains	Unrealized Losses	Fair Value
Commercial paper	\$ 1,494	\$ —	\$ —	\$ 1,494
U.S. Treasuries	61,891	19	(53)	61,857
U.S. Government agency bonds	93,987	89	(98)	93,978
Corporate bonds	74,937	148	(1)	75,084
Total marketable securities	<u>\$ 232,309</u>	<u>\$ 256</u>	<u>\$ (152)</u>	<u>\$ 232,413</u>

The following table shows the fair value and gross unrealized losses of the Company's marketable securities aggregated by major security type and length of time that the individual securities have been in a continuous unrealized loss position (in thousands):

As of December 31, 2025	Less Than 12 Months		Greater Than 12 Months	
	Fair Value	Unrealized Losses	Fair Value	Unrealized Losses
U.S. Government agency bonds	\$ 450	\$ —	\$ 3,589	\$ (1)
Corporate bonds	19,759	(20)	—	—
Total	\$ 20,209	\$ (20)	\$ 3,589	\$ (1)

As of December 31, 2024	Less Than 12 Months	
	Fair Value	Unrealized Losses
U.S. Treasuries	\$ 35,085	\$ (53)
U.S. Government agency bonds	32,148	(98)
Corporate bonds	7,415	(1)
Total	\$ 74,648	\$ (152)

As of December 31, 2024, the Company did not have any individual securities in a continuous loss position for greater than 12 months, and there were no individual securities that were in a significant unrealized loss position. For marketable securities in an unrealized loss position, the Company does not intend to sell them, it is not more likely than not that the Company will be required to sell them before recovery of their amortized cost bases, and the unrealized losses are not credit related.

During the year ended December 31, 2025, the Company sold \$35.9 million of marketable securities. Realized gains related to the sale were not material. The Company did not sell any marketable securities or record any realized gains or losses for the year ended December 31, 2024 and 2023. Additionally, the Company has not recorded any impairment losses or a credit loss allowance on its marketable securities.

At December 31, 2025 and 2024, the Company had \$0.7 million and \$1.6 million of accrued interest receivable on its marketable securities, respectively.

The following table summarizes the contractual maturities of the Company's marketable securities (in thousands):

	As of December 31, 2025		As of December 31, 2024	
	Amortized cost	Fair Value	Amortized Cost	Fair Value
Due within one year	\$ 71,054	\$ 71,141	\$ 197,141	\$ 197,306
Due in one to two years	17,260	17,252	35,168	35,107
Total marketable securities	\$ 88,314	\$ 88,393	\$ 232,309	\$ 232,413

Note 8. Fair Value of Financial Instruments

Recurring Fair Value Measurements

The following tables present the Company's fair value hierarchy for its financial instruments that are measured at fair value on a recurring basis (in thousands):

As of December 31, 2025	Total	Quoted prices in active markets (Level 1)	Significant other observable inputs (Level 2)	Significant unobservable inputs (Level 3)
Financial assets:				
Cash equivalents: (1)				
Money market funds	\$ 17,219	\$ 17,219	\$ —	\$ —
Total cash equivalents	17,219	17,219	—	—
Marketable securities:				
Commercial paper	12,877	—	12,877	—
U.S. Treasuries	29,650	—	29,650	—
U.S. Government agency bonds	13,594	—	13,594	—
Corporate bonds	32,272	—	32,272	—
Total marketable securities	88,393	—	88,393	—
Total financial assets	\$ 105,612	\$ 17,219	\$ 88,393	\$ —
Financial liabilities:				
Contingent liabilities (2)	\$ 5,684	\$ —	\$ —	\$ 5,684
Total financial liabilities	\$ 5,684	\$ —	\$ —	\$ 5,684

As of December 31, 2024	Total	Quoted prices in active markets (Level 1)	Significant other observable inputs (Level 2)	Significant unobservable inputs (Level 3)
Financial assets:				
Cash equivalents: (1)				
Money market funds	\$ 44,426	\$ 44,426	\$ —	\$ —
Total cash equivalents	44,426	44,426	—	—
Marketable securities:				
Commercial paper	1,494	—	1,494	—
U.S. Treasuries	61,857	—	61,857	—
U.S. Government agency bonds	93,978	—	93,978	—
Corporate bonds	75,084	—	75,084	—
Total marketable securities	232,413	—	232,413	—
Total financial assets	\$ 276,839	\$ 44,426	\$ 232,413	\$ —

(1) Included in cash and cash equivalents on the Consolidated Balance Sheets.

(2) The Company's recurring fair value measurements using Level 3 inputs related to the Company's contingent consideration liability from the acquisition of Emission and the contingent liability assumed in the acquisition of Akoya.

Cash equivalents and marketable securities classified as Level 2 financial assets are initially valued at their purchase price and subsequently valued at the end of each reporting period utilizing third party pricing services or other observable data. The pricing services utilize industry standard valuation methods, including both income and market-based approaches, and observable market inputs to determine the fair value. These observable market inputs include reportable trades, benchmark yields, credit spreads, broker/dealer quotes, bids, offers, current spot rates, and other industry and economic events.

Level 3 Financial Instruments

The following table presents the changes in the Company's Level 3 financial instruments measured at fair value on a recurring basis, which consist of contingent liabilities:

	Level 3 Liabilities		
	Emission (1)	PKI License (2)	Total
Balance as of December 31, 2024	\$ —	\$ —	\$ —
Acquisition	6,612	3,619	10,231
Change in fair value of contingent liabilities	(4,624)	77	(4,547)
Balance as of December 31, 2025	\$ 1,988	\$ 3,696	\$ 5,684

(1) Earnout 2 requires additional consideration to be paid to the selling shareholders based on the amount and timing of certain performance targets.

Earnout 2 is measured and paid over a five year period ending December 2029.

(2) As part of Akoya's 2018 acquisition of the Quantitative Pathology Solutions division of Perkin Elmer, Inc., subsequently known as Revvity, Inc. ("PKI"), Akoya entered into a license agreement with PKI (the "PKI License") (refer to Note 16 - *Commitments and Contingencies*). The Company recognizes the assumed contingent liability at fair value in accordance with ASC 805. The PKI License is measured and paid over the remaining eight year period ending March 2033.

Monte-Carlo simulations and discounted cash flow analyses were used to determine the fair values of the Level 3 liabilities, including the following significant unobservable inputs: projected revenue, a risk adjusted discount rate, and revenue volatility. Changes in fair value subsequent to the acquisition date were due to updated valuation inputs and the passage of time. Increases or decreases in the inputs would have resulted in higher or lower fair value measurements.

The range of outcomes payable for Earnout 2 is zero to \$50.0 million. It is not possible to estimate a range of outcomes payable for the PKI License as there is no cap on the amount that could be earned.

The fair value of the contingent liabilities are recorded in accrued expenses and other current liabilities and non-current portion of contingent liabilities on the Consolidated Balance Sheets. Changes in fair value are recorded in change in fair value of contingent liabilities on the Consolidated Statements of Operations.

Other Fair Value Disclosures

During the years ended December 31, 2025 and 2024, the Company did not transfer financial assets between levels of the fair value hierarchy. Additionally, there have been no changes to the valuation techniques for Level 2 or Level 3 financial assets.

Note 9. Inventory

Inventory, net of inventory reserves, consisted of the following (in thousands):

	As of December 31,	
	2025	2024
Raw materials	\$ 13,727	\$ 7,215
Work in process	11,030	7,980
Finished goods	30,006	17,580
Total inventory	\$ 54,763	\$ 32,775

Note 10. Property and Equipment

Property and equipment consisted of the following (in thousands):

	As of December 31,	
	2025	2024
Laboratory and manufacturing equipment	\$ 20,211	\$ 14,648
Demo inventory	2,650	—
Furniture and fixtures	7,332	1,999
Computers and software	3,174	6,036
Leasehold improvements	16,132	13,291
Total cost	49,499	35,974
Less: accumulated depreciation	(25,827)	(18,824)
Property and equipment, net	\$ 23,672	\$ 17,150

The Company incurred depreciation expense of \$7.0 million, \$4.8 million, and \$4.7 million for the years ended December 31, 2025, 2024, and 2023, respectively. Substantially all of the Company's property and equipment is located in North America.

There were no material disposals during the years ended December 31, 2025, 2024, and 2023. For the years ended December 31, 2025, and 2024 the Company did not record any impairments related to property and equipment. Refer to Note 15 - *Leases* for discussion related to the impairment of leasehold improvements in 2023.

As of December 31, 2025 and 2024, total gross capitalized software development costs were \$4.4 million and \$4.3 million, respectively. Depreciation of capitalized software costs was \$1.2 million for the year ended December 31, 2025, and not material for the years ended December 31, 2024, and 2023.

Note 11. Accrued Expenses and Other Current Liabilities

Accrued expenses and other current liabilities consisted of the following (in thousands):

	As of December 31,	
	2025	2024
Accrued professional services	\$ 2,766	\$ 4,897
Accrued royalties	1,784	1,361
Accrued tax liabilities	1,125	1,018
Acquisition holdback (1)	1,000	—
Off-market liability (2)	6,869	—
Other accrued expenses	4,027	1,575
Total accrued expenses and other current liabilities	\$ 17,571	\$ 8,851

(1) Represents the holdback associated with the Emission acquisition. Refer to Note 3 - *Acquisitions*.

(2) Represents the current portion of an off-market component of an acquired customer contract in the acquisition of Akoya. Refer to Note 3 - *Acquisitions*.

Note 12. Stock-Based Compensation

Stock-Based Compensation Plans

In December 2017, the Company adopted the 2017 Employee, Director and Consultant Equity Incentive Plan (the "2017 Plan"), under which it may grant incentive stock options, non-qualified stock options, RSUs, and other stock-based awards. As of December 31, 2017, the 2017 Plan allowed for the issuance of (1) up to 1.0 million shares of common stock and (2) up to 2.5 million shares of common stock represented by awards granted under the 2007 Stock Option and Grant Plan (which was terminated upon completion of the Company's initial public offering) that were forfeited, expired, or cancelled without delivery of shares or which result in the forfeiture of shares of common stock back to the Company on or

after the date the 2017 Plan became effective. The 2017 Plan contains an “evergreen” provision, which allows for an annual increase in the number of shares of common stock available for issuance under the 2017 Plan on the first day of each fiscal year during the period beginning in 2019 and ending in 2027. The annual increase is equal to the lowest of (1) 4% of the number of shares of common stock outstanding as of such date and (2) an amount determined by the Company’s board of directors or compensation committee. As of December 31, 2025, 6.2 million shares were outstanding and there were 4.9 million shares available for grant under the 2017 Plan.

In December 2017, the Company adopted the 2017 Employee Stock Purchase Plan (the “2017 ESPP”). As of December 31, 2019, the 2017 ESPP allowed for the issuance of up to 0.6 million shares of common stock. The 2017 ESPP contains an “evergreen” provision, which allows for an increase in the number of shares under the plan on the first day of each fiscal year beginning with 2018 and ending in 2027. The increase is equal to the lowest of: (1) 1% of the number of shares of common stock outstanding on the last day of the immediately preceding fiscal year and (2) an amount determined by the Company’s board of directors or compensation committee. As of December 31, 2025, 1.9 million shares were available for grant under the 2017 ESPP. The 2017 ESPP provides for six-month offering periods commencing and ending on March 1 through August 31, and September 1 through February 28. During the years ended December 31, 2025, 2024, and 2023, employees purchased 0.2 million, 0.1 million, and 0.1 million shares, respectively, of the Company’s common stock pursuant to the 2017 ESPP.

In July 2025, the Company adopted the Quanterix Corporation Restricted Stock Unit Inducement Awards Plan (the "Inducement Plan"). The Inducement Plan allows for issuance of up to 0.5 million shares of Quanterix common stock, consisting of: (i) up to 0.2 million shares of Quanterix common stock that may be become available for issuance again under the 2021 Akoya Equity Incentive Plan pursuant to the terms of such plan and (ii) 0.3 million shares of Quanterix common stock issuable upon vesting of restricted stock units granted to Akoya employees as an inducement to employment with Quanterix following the Merger.

In November 2025, the Company adopted the 2025 Inducement Plan which allows for issuance of up to 0.9 million shares of Quanterix Common Stock, par value \$0.001 per share.

In accordance with Nasdaq listing rules, equity awards issued under the Inducement Plan and the 2025 Inducement Plan are restricted to individuals who are not already employees or directors of the Company.

Stock Options

Under the 2017 Plan, stock options may not be granted with exercise prices of less than fair market value on the date of the grant. Options generally vest ratably over a four-year period with 25% vesting on the first anniversary and the remaining 75% vesting ratably on a monthly basis over the remaining three years. These options expire ten years after the grant date.

Stock option activity for the year ended December 31, 2025 is presented below (in thousands, except per share and contractual life amounts):

	Number of options	Weighted-average exercise price per share	Weighted-average remaining contractual life (in years)	Aggregate intrinsic value
Outstanding at December 31, 2024	3,563	\$ 19.94	7.7	\$ 678
Granted	2,676	8.21		
Exercised	(4)	3.12		
Forfeited/expired	(1,355)	15.04		
Outstanding at December 31, 2025	4,880	\$ 14.89	7.9	\$ 170
Exercisable at December 31, 2025	2,042	\$ 20.06	6.5	\$ 12
Vested and expected to vest at December 31, 2025	4,880	\$ 14.89	7.9	\$ 170

The weighted average grant-date fair value per share of awards granted was \$5.85 in 2025, \$14.43 in 2024, and \$10.63 in 2023. The total intrinsic value of stock options exercised was not material in 2025, \$1.9 million in 2024, and \$1.9 million in 2023.

Restricted Stock Units

RSUs represent the right to receive shares of common stock upon meeting specified vesting requirements. Shares are delivered to the grantee upon vesting, less shares for the payment of withholding taxes. RSU activity for the year ended December 31, 2025 is presented below (in thousands, except per share amounts):

	Number of shares	Weighted-average grant date fair value per share
Unvested as of December 31, 2024	1,115	\$ 18.55
Assumed through acquisition of Akoya (1)	253	6.54
Granted	1,619	7.76
Vested	(677)	10.44
Forfeited	(735)	13.58
Unvested as of December 31, 2025	1,575	\$ 11.33
Expected to convert at December 31, 2025	1,575	\$ 11.33

(1) Refer to Note 3 - *Acquisitions*.

The weighted average grant-date fair value per share of awards granted was \$7.76 in 2025, \$21.51 in 2024, and \$15.90 in 2023. The total fair value of shares that vested was \$13.0 million in 2025, \$10.3 million in 2024, and \$10.4 million in 2023.

Stock-Based Compensation Expense

Stock-based compensation expense was recorded in the following categories on the Consolidated Statements of Operations (in thousands):

	Year Ended December 31,		
	2025	2024	2023
Cost of product revenue	\$ 901	\$ 1,132	\$ 839
Cost of service and other revenue	970	1,142	1,124
Research and development	2,178	2,117	1,713
Selling, general and administrative	16,669	15,596	13,147
Total stock-based compensation expense	\$ 20,718	\$ 19,987	\$ 16,823

As of December 31, 2025, there was \$30.7 million of total unrecognized stock-based compensation expense related to unvested RSUs and stock options, which is expected to be recognized over the remaining weighted-average vesting period of 2.5 years.

The fair value of the Company's stock options granted and purchase rights to the ESPP were estimated using the Black-Scholes valuation model with the following assumptions:

	Year Ended December 31,		
	2025	2024	2023
Stock Options:			
Risk-free interest rate	3.7% - 4.4%	3.4% - 4.6%	3.5% - 4.7%
Expected dividend yield	None	None	None
Expected term (in years)	5.4 - 5.7	5.3 - 5.4	5.0 - 5.2
Expected volatility	82.1% - 84.5%	81.4% - 83.1%	71.1% - 83.1%
Weighted-average grant date fair value per share	\$ 5.85	\$ 14.43	\$ 10.63
Employee Stock Purchase Plan:			
Risk-free interest rate	3.8% - 4.0%	4.3% - 4.9%	5.2% - 5.5%
Expected dividend yield	None	None	None
Expected term (in years)	0.5	0.5	0.5
Expected volatility	72.3% - 91.1%	54.9% - 66.0%	72.8% - 82.5%
Weighted-average grant date fair value per share	\$ 1.39	\$ 3.84	\$ 3.19

Note 13. Net Loss Per Share

The following table presents the computation of basic and diluted net loss per share (in thousands, except per share data):

	Year Ended December 31,		
	2025	2024	2023
Numerator:			
Net loss	\$ (107,150)	\$ (38,531)	\$ (28,354)
Denominator:			
Weighted average common shares outstanding, basic and diluted	42,639	38,367	37,594
Net loss per share, basic and diluted	\$ (2.51)	\$ (1.00)	\$ (0.75)

As the Company was in a net loss position for all periods, the following common share equivalents were excluded from the calculation of diluted net loss per share because their effect would be anti-dilutive (in thousands):

	Year Ended December 31,		
	2025	2024	2023
Stock options	5,407	3,561	2,783
Common stock and RSUs	1,954	1,341	1,513
Estimated ESPP purchases	36	8	23
Total dilutive shares	7,397	4,910	4,319

Note 14. Income Taxes

The following table presents the components of loss before income taxes (in thousands):

	Year Ended December 31,		
	2025	2024	2023
United States	\$ (112,811)	\$ (39,968)	\$ (30,355)
Foreign	540	1,871	2,720
Loss before income taxes	\$ (112,271)	\$ (38,097)	\$ (27,635)

The following table presents the components of the Company's income tax benefit (expense) (in thousands):

	Year Ended December 31,		
	2025	2024	2023
Current taxes:			
United States - State	\$ (93)	\$ (37)	\$ (161)
Foreign	(416)	(683)	(850)
Total current taxes	(509)	(720)	(1,011)
Deferred taxes:			
United States			
Federal	4,447	—	—
State	802	—	—
Foreign	381	286	292
Total deferred taxes	5,630	286	292
Income tax benefit (expense)	\$ 5,121	\$ (434)	\$ (719)

Effective Tax Rate Reconciliation

In accordance with the Company's prospective adoption of ASU 2023-09 effective for its annual financial statements beginning January 1, 2025, the following table is a reconciliation of the Company's income taxes and its effective tax rate as compared to the United States ("U.S.") federal statutory income tax rate of 21.0% for the year ended December 31, 2025 (in thousands, except percentages):

	Year Ended December 31, 2025	
	Amount	Rate
U.S. federal statutory income tax	\$ 23,577	21.0 %
U.S. state and local income tax, net of federal income tax effect (1)	555	0.5 %
Foreign tax effects	(55)	— %
Effect of cross border transactions	(72)	(0.1)%
Nontaxable or nondeductible items		
Transaction costs	(2,187)	(1.9)%
Goodwill impairment	(1,339)	(1.2)%
Stock compensation	(2,281)	(2.0)%
Limitation on executive compensation	(710)	(0.6)%
Other nondeductible items	(1,235)	(1.1)%
Tax credits	293	0.3 %
Change in valuation allowance	(11,411)	(10.3)%
Other items	(14)	— %
Income tax benefit and effective tax rate	\$ 5,121	4.6 %

(1) The state that contributes to the majority of the state and local tax effect is Massachusetts.

The Company's effective income tax rate for 2025 differs from the U.S. federal statutory rate primarily as a result of the valuation allowance maintained against the Company's net deferred tax assets as well as non-deductible items including stock-based compensation, transactions costs, and goodwill impairment.

A reconciliation of the Company's effective tax rate as compared to the U.S. federal statutory income tax rate of 21.0% for the years ended December 31, 2024 and 2023 in accordance with the guidance prior to the adoption of ASU 2023-09, was as follows:

	Year Ended December 31,	
	2024	2023
U.S. federal statutory tax rate	21.0 %	21.0 %
U.S. state income taxes, net of federal benefit	2.7 %	(0.3)%
Change in valuation allowance	(22.5)%	(24.3)%
Non-deductible stock compensation	(2.4)%	(1.6)%
Permanent items	(0.3)%	(0.6)%
Non-deductible executive compensation	(3.2)%	(2.6)%
Tax credits	4.0 %	4.9 %
Other	(0.4)%	0.9 %
Effective tax rate	(1.1)%	(2.6)%

Deferred Tax Assets and Liabilities

Deferred tax assets and liabilities reflect the net tax effects of temporary differences between the carrying amounts of assets and liabilities for financial reporting and the amounts used for income tax purposes. The Company records deferred tax assets in other non-current assets and deferred tax liabilities in other non-current liabilities on the Consolidated Balance Sheets. Significant components of the Company's deferred tax assets and liabilities were as follows (in thousands):

	As of December 31,	
	2025	2024
Deferred tax assets:		
Net operating loss carryforwards	\$ 147,973	\$ 80,541
Tax credits	10,929	10,178
Deferred revenue	6,322	2,412
Amortization	—	645
Stock-based compensation	4,539	3,808
Inventory	4,092	1,069
Capitalized R&D costs	16,784	13,753
Lease liabilities	8,802	9,093
Other deferred tax assets	653	497
Non deductible interest - 163(j)	3,839	—
Accrual and reserves	5,107	1,876
Gross deferred tax assets	209,040	123,872
Less: valuation allowances	(177,108)	(118,663)
Total deferred tax assets	31,932	5,209
Deferred tax liabilities:		
Right-of-use assets	(3,927)	(3,970)
Depreciation	(1,474)	(1,137)
Amortization acquired intangibles	(26,702)	(831)
Other deferred tax liabilities	(273)	(71)
Total deferred tax liabilities	(32,376)	(6,009)
Net deferred tax liabilities	\$ (444)	\$ (800)

The change in the Company's valuation allowance was as follows (in thousands):

	2025	2024
Balance at December 31 of prior year	\$ 118,663	\$ 110,082
Change in valuation allowance	58,445	8,581
Balance at December 31	\$ 177,108	\$ 118,663

The increase in the valuation allowance during the year ended December 31, 2025 was primarily due to U.S. operating losses incurred. The increase was partially offset by the release of a portion of the valuation allocation related to taxable temporary differences recorded as part of the Emission and Akoya acquisitions. These temporary differences are treated as a source of income for pre-existing U.S. federal and state deferred tax assets.

In determining the need for a valuation allowance, the Company considers the cumulative book income and loss positions of each of its entities, as well as its worldwide cumulative loss position. The Company has assessed, on a jurisdictional basis, the available means of recovering deferred tax assets, including the ability to carry back net operating losses ("NOLs"), the existence of reversing taxable temporary differences, the availability of tax planning strategies, and forecasted future taxable income. At December 31, 2025, the Company maintained a full valuation allowance against its worldwide net deferred tax assets, as it concluded that it was more likely than not that the deferred assets will not be utilized.

Tax Attributes

As of December 31, 2025, the Company had U.S. federal NOLs of approximately \$597.7 million. U.S. federal NOLs generated through December 31, 2017 of approximately \$111.1 million expire at various dates through 2037, and U.S. federal NOLs generated after December 31, 2017 of approximately \$486.6 million do not expire. As of December 31, 2025, the Company had U.S. federal tax credit carryforwards of approximately \$13.2 million that expire at various dates through 2045.

As of December 31, 2025, the Company had \$374.4 million of state NOLs, approximately \$348.4 million of which expire at various dates through 2045, and approximately \$26.0 million of which do not expire. As of December 31, 2025, the Company had U.S. state tax credit carryforwards of approximately \$6.6 million that expire at various dates through 2040.

Under Sections 382 and 383 of the U.S. Internal Revenue Code, if a corporation undergoes an ownership change, the corporation's ability to use its pre-ownership change NOLs and other pre-ownership change tax attributes, such as research tax credits, to offset its post-change income and taxes may be limited. Generally, an ownership change occurs if there is a cumulative change in an entity's ownership by 5% stockholders that exceeds 50 percentage points over a rolling three-year period. Similar rules may apply under U.S. state tax laws. Under the Tax Cuts and Jobs Act of 2017 ("TCJA"), the use of federal NOLs arising in taxable years beginning after December 31, 2017 is limited to 80% of current year taxable income and NOLs arising in taxable years ending after December 31, 2017 may not be carried back (though any such NOLs may be carried forward indefinitely).

The Company may have experienced an ownership change in the past and may experience ownership changes in the future as a result of future transactions in its share capital, some of which may be outside of the control of the Company. As a result, if the Company earns net taxable income, its ability to use its pre-ownership change NOLs, or other pre-ownership change tax attributes, to offset U.S. federal and state taxable income and taxes may be subject to significant limitations. As of December 31, 2025, the Company has not completed a Section 382 analysis.

Unrecognized Tax Benefits

Liabilities for unrecognized tax benefits related to uncertain tax positions are classified as other unrecognized tax benefits. These unrecognized benefits are reviewed on an ongoing basis and are adjusted for changing facts and circumstances, including management's judgment in the interpretation of applicable tax law, regulation of tax rulings, the progress of tax audits, and the closing of statute of limitations.

For the year ended December 31, 2025, the Company increased its gross unrecognized tax benefits \$9.8 million, primarily due to tax benefits relating to federal and state research and development tax credits. These unrecognized tax benefits were related to tax positions taken by Akoya in prior years. The unrecognized tax benefits were recorded as a

reduction to the related deferred tax asset. If recognized in the future, the entire \$9.8 million would not impact the Company's effective tax rate due to the current valuation allowance recorded. As a result, it is estimated that any future release would increase the valuation allowance and not generate an income rate benefit. Prior to 2025, the Company did not have any unrecognized tax benefits.

At December 31, 2025, the Company had no accrued interest or penalties related to uncertain tax positions. Interest and penalty charges, if any, would be classified as income tax benefit (expense) in the Consolidated Statements of Operations.

Other Tax Disclosures

The Company is subject to taxation in the United States as well as the Netherlands, Sweden, China, and the United Kingdom. At December 31, 2025, the Company is generally no longer subject to examination by taxing authorities in the United States for years prior to 2022. However, NOLs and tax credits in the United States may be subject to adjustments by taxing authorities in future years in which they are utilized. The Company's foreign subsidiaries remain open to examination by taxing authorities from 2018 onward.

On July 4, 2025, the One Big Beautiful Bill Act ("OBBBA") was enacted in the United States. The OBBBA includes significant provisions, such as the permanent extension of certain expiring provisions of the Tax Cuts and Jobs Act, modifications to the international tax framework, and the restoration of favorable tax treatment for certain business provisions. The legislation has multiple effective dates, with certain provisions effective in 2025 and others implemented through 2027. These changes are reflected in our results for the year ended December 31, 2025 and did not have a material impact.

Cash Paid for Taxes

The following table summarizes income taxes paid in 2025 (in thousands):

	Year Ended December 31, 2025
U.S. state	
Massachusetts	\$ —
Other states	44
Foreign	
Sweden	1,220
Other foreign jurisdictions	140
Total income taxes paid	\$ 1,404

Cash paid for taxes for the years ended December 31, 2024 and 2023 was \$0.9 million and \$0.8 million, respectively.

Note 15. Leases

During the year ended December 31, 2025, the Company acquired certain leases through the acquisition of Akoya and did not enter into any material leases. At December 31, 2025, the Company had no additional leases which had not yet commenced. Substantially all of the Company's leases are located in North America.

The Company has leased facilities, including leases acquired in acquisitions, that are vacant and are not used in its operations. At the point a decision is made to stop use of a leased facility, the Company moves the related ROU asset and leasehold improvements into their own asset group and performs quarterly impairment assessments. The impairment analyses evaluate the present value of net cash flows under the original leases and the estimated cash flows under an estimated sublease and considered industry and economic factors such as rental rates, interest rates, and recent real estate activities to estimate the net cash flows analysis and impairment amount. These assessments indicated the ROU asset for certain facilities were fully impaired, which resulted in the Company recording an impairment charge of \$1.4 million during the year ended December 31, 2025. During the year ended December 31, 2024, the Company did not record any

lease related impairments. During the year ended December 31, 2023, the Company recorded a \$1.3 million impairment charge related to unused ROU assets and leasehold improvements.

The components of the lease costs and supplemental cash flow information relating to the Company's leases were as follows (in thousands):

	Year Ended December 31,		
	2025	2024	2023
Operating lease cost	\$ 5,970	\$ 4,949	\$ 5,209
Short-term and variable lease cost	5,692	4,359	3,996
Total operating lease cost	\$ 11,662	\$ 9,308	\$ 9,205

	Year Ended December 31,	
	2025	2024
Cash paid for amounts included in the measurement of operating lease liabilities	\$ 9,887	\$ 8,554
Weighted average remaining lease term - operating leases (in years)	4.4	5.8
Weighted average discount rate - operating leases	7.71%	7.92%

The undiscounted future lease payments for non-cancelable operating leases as of December 31, 2025 were as follows (in thousands):

Maturity of lease liabilities as of December 31, 2025	Operating Leases
2026	\$ 10,205
2027	8,952
2028	8,395
2029	8,570
2030	7,028
Thereafter	695
Total lease payments	43,845
Less imputed interest	(6,606)
Total operating lease liabilities	\$ 37,239

Note 16. Commitments and Contingencies

Purchase Commitments

The Company's non-cancellable purchase commitments primarily consist of purchases of raw materials for manufacturing operations under annual and multi-year agreements, some of which have minimum quantity requirements. The Company's total purchase commitments under these agreements as of December 31, 2025 was \$1.1 million, most of which the Company expects to incur in the year ending December 31, 2026.

License Agreements

Eli Lilly and Company

In February 2022, the Company and Eli Lilly and Company ("Lilly") entered into a Technology License Agreement (the "Lilly License") under which Lilly granted a non-exclusive license to Lilly's proprietary p-Tau 217 antibody technology for use by the Company in RUO products, services, and future in vitro diagnostics ("IVD") applications within the field of Alzheimer's disease. Pursuant to the Lilly License, the Company paid an upfront fee, is required to make milestone payments based on the achievement of predetermined regulatory and commercial events, and will pay royalties on net sales of licensed products.

Harvard University

In August 2022, the Company and Harvard University ("Harvard") entered into an exclusive license agreement (the "Harvard License Agreement") for certain intellectual property owned by Harvard. Under the Harvard License Agreement, the Company paid an upfront fee, and is required to pay Harvard low single-digit royalties on net sales of products and services using the licensed technology as well as a portion of its applicable sublicense revenues. The Company incurred no royalty expense under the Harvard License Agreement for the years ended December 31, 2025, 2024 and 2023, respectively. Refer to Note 17 - *Related Party Transactions* for a discussion of a related party relationship with Harvard.

Tufts University

In June 2007, the Company and Tufts University ("Tufts") entered into a license agreement (the "Tufts License Agreement") for certain intellectual property owned by Tufts. The Tufts License Agreement, which was subsequently amended, is exclusive and sub-licensable, and will continue in effect on a country by country basis as long as there is a valid claim of a licensed patent in a country. The Company is contractually obligated to pay license and maintenance fees that are creditable against royalties, in addition to low single-digit royalties on direct sales and services, and a royalty on sublicense income. The Company incurred royalty expenses related to the Tufts License Agreement of \$1.5 million, \$2.1 million, and \$1.7 million, during the years ended December 31, 2025, 2024, and 2023, respectively, which was recorded in cost of product revenue on the Consolidated Statements of Operations. Refer to Note 17 - *Related Party Transactions* for a discussion of a related party relationship with Tufts.

Stanford University

In November 2015, Akoya and Stanford University ("Stanford") entered into a license agreement (the "Stanford License Agreement") to use certain patent rights owned by Stanford to make and sell products and services (the "Stanford Licensed Products"). The Stanford License Agreement, which was subsequently amended, is exclusive and sublicensable, and the Company is obligated to use commercially reasonable efforts to develop, manufacture, sell, and develop markets for Stanford Licensed Products. The Stanford License Agreement will continue until the expiration, revocation, invalidation or abandonment of the last licensed patent or patent application, the last of which is set to expire in 2036, unless terminated earlier in accordance with its terms. Subject to Stanford's approval, the Company controls the prosecution and maintenance of the licensed patents.

Under the Stanford License Agreement, the Company is required to pay Stanford annual license maintenance fees, make certain milestone payments based on specified patent issuance and sales milestone events, and pay Stanford low single-digit royalty on net sales of Stanford Licensed Products and a portion of any sublicensing income. Royalty expenses related to the Stanford License Agreement incurred during the year ended December 31, 2025 were not material.

University of Washington

In June 2018, Akoya and the University of Washington (the "University") entered into an license agreement (the "University License Agreement") to use certain patent rights owned by the University to make and sell products and services (the "University Licensed Products"). The University License Agreement is exclusive and sublicensable, and the Company is obligated to use commercially reasonable efforts to develop, manufacture, sell, and develop markets for University Licensed Products. The University License Agreement will continue until the expiration, revocation, invalidation or abandonment of the last licensed patent or patent application, the last of which is set to expire in 2032, unless terminated earlier in accordance with its terms. Subject to the University's approval, the Company controls the prosecution and maintenance of the licensed patents.

Under the University License Agreement, the Company is required to make certain milestone payments based on the achievement of certain commercial milestones and pay low single-digit royalty on net sales of University Licensed Products, subject to certain minimum annual royalty payments and potential reductions based on a royalty-stacking allowance for certain third-party rights, as well as sharing a portion of any non-royalty sublicensing income. Royalty expenses related to the University License Agreement incurred during the year ended December 31, 2025 were not material.

Perkin Elmer

Under the PKI License acquired as part of the acquisition of Akoya, PKI granted Akoya an exclusive, nontransferable, sublicensable license under certain patent rights to make, use, import, and commercialize certain products and services. The Company is required to pay single digit royalties through March 2033 on net sales of products and services that are covered by patent rights under the PKI License. Royalty expense under the PKI License for the year ended December 31, 2025 was not material. Refer to Note 8 - *Fair Value of Financial Instruments* for further discussion.

Legal Contingencies

The Company is subject to claims in the ordinary course of business; however, the Company is not currently a party to any pending or threatened litigation, the outcome of which would be expected to have a material adverse effect on its financial condition or results of operations. The Company accrues for contingent liabilities when losses are probable and estimable. If an estimate of a probable loss is a range and no amount within the range is more likely than any other amount in the range, the Company accrues the minimum amount of the range.

Note 17. Related Party Transactions

Akoya, which the Company acquired in July 2025, is party to a diagnostic development agreement with a biopharmaceutical customer of Akoya's. A member of the Company's board of directors also serves on the board of directors of the biopharmaceutical customer. Revenue recorded from sales of products and services to this customer was \$4.8 million for the year ended December 31, 2025. Cost of product revenue for goods and services delivered to this customer was \$4.6 million for the year ended December 31, 2025. At December 31, 2025, the Company had \$8.2 million of deferred revenue related to this customer and did not have any open payables to or open receivables from the customer.

In the first quarter of 2025, the Company entered into agreements with two entities owned by selling shareholders of Emission (refer to Note 3 - *Acquisitions*) to continue development work on certain future products for Quanterix. At December 31, 2025, the Company did not have any open payable balances with these entities. During 2025, the Company paid Earnout 1 (refer to Note 3 - *Acquisitions*), a portion of which was paid to these entities.

In August 2022, the Company entered into the Harvard License Agreement for certain intellectual property owned by Harvard (refer to Note 16 - *Commitments and Contingencies*). Harvard is obligated to pay a portion of the payments received from the Company under the Harvard License Agreement to a former member of the Company's board of directors. This individual is also affiliated with Harvard and Mass General Brigham. Revenue recorded from sales of products and services to Harvard and its affiliates and to Mass General Brigham and its affiliates totaled \$1.5 million, \$2.2 million, and \$1.3 million for the years ended December 31, 2025, 2024, and 2023, respectively. Cost of product revenue and operating expenses with Harvard and its affiliates and Mass General Brigham and its affiliates were not material for the years ended December 31, 2025, 2024, and 2023, respectively. At December 31, 2025 and 2024, open payables to and receivable balances from Harvard and Mass General Brigham were not material.

In June 2007, the Company entered into the Tufts License Agreement for certain intellectual property owned by Tufts (refer to Note 16 - *Commitments and Contingencies*). A former member of the Company's board of directors who served on the board of directors during the year ended December 31, 2025, was previously affiliated with Tufts. This Board member continued to receive compensation from Tufts on a formulaic basis based on royalties and license payments the Company makes to Tufts while on the Board of Directors of the Company. At December 31, 2025 and 2024, open payable balances to Tufts were not material.

Note 18. Segment Reporting

Operating segments are defined as components of an enterprise about which separate discrete information is available for evaluation by the chief operating decision-maker ("CODM"), in deciding how to allocate resources and assess performance. The Company's CODM is the chief executive officer.

As a result of the acquisition of Akoya, the Company reassessed its operating segments and concluded that the Company continues to operate as one reportable segment as of December 31, 2025. This operating segment is focused on development and commercialization of comprehensive protein biomarker solutions that identify signatures in blood and tissue to provide insights to providers, patients, and research organizations.

The Company utilizes consolidated net loss as the measure of segment profitability (loss) as required by ASU 2023-07. The CODM uses this measure, along with the significant revenue and expense lines included in the table below, when analyzing the Company's operations and performance and determining how to allocate resources. These measures are consistently used by the CODM in comparing budgeted results versus actuals, in determining when or where to invest resources into the business, and for decisions on strategic initiatives, all of which is assessed at the consolidated level.

The following table presents the reconciliation of significant segment information reviewed by the CODM to consolidated net loss:

	Year Ended December 31,		
	2025	2024	2023
Revenues:			
Revenue from contracts with customers (Note 5)	\$ 138,654	\$ 135,436	\$ 121,139
Grant revenue	243	1,985	1,229
Total revenues	138,897	137,421	122,368
Less:			
Costs of goods sold and services, including shipping and handling costs	79,519	62,430	56,290
Certain operating expenses, excluding shipping and handling costs (1)	168,349	124,587	107,029
Other segment items (2)	(1,821)	(11,065)	(12,597)
Consolidated net loss	\$ 107,150	\$ 38,531	\$ 28,354

(1) Expenses consist of research and development and selling, general and administrative from the Consolidated Statements of Operations and exclude shipping and handling costs.

(2) Other segment items represent discrete events, non-recurring transactions, or insignificant items that are not used by the CODM to evaluate the Company's performance or allocate resources, and include:

- a. Impairment and restructuring costs – impairment charges for goodwill and other long-lived assets, and costs associated with approved restructuring plans, including costs to implement restructurings and employee separation costs;
- b. Change in fair value of contingent liabilities – changes in the fair value of contingent payments as a result of updated valuation inputs;
- c. Other lease costs – amortization of operating lease right-of-use assets and other facility operating expenses from leased facilities not in use;
- d. Interest income – interest earned on cash, cash equivalents, and marketable securities, and the accretion of discounts on marketable securities;
- e. Other income (expense), net – gains and losses on foreign currency, and other non-recurring items that are not a part of the Company's core business operations; and
- f. Income tax benefit (expense) – income taxes related to federal, state, and foreign jurisdictions in which the Company conducts business.

The CODM also reviews consolidated balance sheet accounts and activity including cash usage and other working capital changes using the balances as reported on the Consolidated Balance Sheets

There have been no changes to the methods used to determine segment profit or loss, or the significant segment captions, across any of the periods presented.

Note 19. Restructuring

In May and July 2025, the Company announced actions to reduce operating costs, preserve cash, and, specific to the July action, realize anticipated synergies and other benefits of the Akoya acquisition. These actions included reductions in force and elimination of duplicate corporate positions and were completed as of December 31, 2025.

During the year ended December 31, 2025, the Company incurred approximately \$8.0 million of expenses related to the restructuring, which also represents the total amount expected to be incurred. These costs were recorded in impairment and restructuring on the Consolidated Statements of Operations and were substantially for cash payments of severance and employee benefits, \$7.7 million of which was paid by December 31, 2025.

The Company did not have any additional restructuring activities during the years ended December 31, 2024 and 2023.

Note 20. Employee Benefit Plans

The Company sponsors a 401(k) savings plan for employees and may make discretionary contributions. During the years ended December 31, 2025, 2024, and 2023, the Company made contributions of \$2.2 million, \$2.4 million, and \$0.8 million, respectively.

Note 21. Subsequent Events

Termination of Diagnostic Development Agreement

As part of the acquisition of Akoya, the Company assumed a diagnostics development agreement (the "Development Agreement") with a biopharmaceutical customer (the "Biopharma Customer"). On February 25, 2026, the Development Agreement was terminated by mutual agreement of the parties and, in connection with such termination, Quanterix will transfer certain know-how to the Biopharma Customer and grant a non-exclusive, sub-licensable, fully paid license of the related intellectual property. No further consideration is due to either party for the know-how transfer or license. The Company is still assessing the impact to its Consolidated Financial Statements related to the off-market liability (refer to Note 3 - *Acquisitions*) of \$14.0 million, deferred revenue of \$8.2 million, and IPR&D intangible asset of \$19.3 million associated with the Development Agreement as of December 31, 2025. The contract termination could result in the potential for recognition of one-time revenue and/or an impairment charge.

Quanterix
900 Middlesex Turnpike
Billerica, MA 01821

tel: 617.301.9400
fax: 617.301.9401

www.quanterix.com

November 20, 2025

Re: Employment Agreement

Dear Bill:

This Employment Agreement (the "Agreement") is entered into between you, William P. Donnelly, and Quanterix Corporation (the "Company"), and is effective as of November 20, 2025 (the "Effective Date").

1. **Role.** As of the Effective Date, you will be employed by the Company as its Executive Chair (which position shall include, but not be limited to, the role and responsibilities of the Chair of the Board of Directors). As the Company's Executive Chair, you will report directly to the lead independent director of the Board of Directors of the Company (the "Board"), and you will be responsible for providing the Company with strategic and other advice and services upon reasonable request by the Board, including, but not limited to, advice and services related to investor relations, employee relations and execution on operational and strategic initiatives. You and the Board agree that this role will not be full-time. The Board acknowledges that your principal location for such work will remain at your home office. During and after the Term (as defined below), your continuing Board membership will remain subject to the rules and requirements of the Company's charter and by-laws, applicable law (including any rules or regulations of any exchange on which the common stock of the Company is listed), and any other requirements of the Board and/or the Nominating and Corporate Governance Committee of the Board.
2. **Term.** This Agreement will begin on the Effective Date and continue until the earlier of (i) December 31, 2026, and (ii) the date that it is terminated in accordance with the provisions of Section 5 (the "Term").
3. **Salary.** Commencing on the Effective Date, the Company will pay you a salary at the rate of \$250,000 per year ("Annual Salary"), payable in accordance with the Company's standard payroll schedule. This salary will not be subject to increase during the Term. You will not be entitled to any compensation under the Company's Non-Employee Director Compensation Policy while employed by the Company, and you will not be entitled to any cash bonus compensation as an employee. You agree that you will not be eligible to participate in any of the Company employee benefit plans, programs, or arrangements during the Term, and you waive your right to participate in any such plans, programs, or arrangements.
4. **Equity Incentive Compensation.** Subject to approval by the Compensation Committee of the Board, you will be entitled to an annual grant of non-qualified stock options equal to 0.15% of the shares of common stock of the Company outstanding on the date of grant, which shall vest ratably on a monthly basis for 24 months following grant. The first grant will occur on or promptly following the Effective Date and, if the Term has not earlier terminated, the second grant will occur on the first anniversary of the Effective Date. Each of your equity grants outstanding as of the Effective Date will continue to vest during the Term, for so long as you continue to provide services to the Company (in your capacity as Executive Chair or as a Board director of the Company), subject to the terms of the applicable equity award agreement and equity plan. You will not be entitled to future equity grants under the Company's Non-Employee Director Compensation Policy while employed by the Company as Executive Chair or in any other role as an employee.
5. **Termination.** Your employment is "at will," meaning you or the Company may terminate it at any time for any or no reason, provided that your employment hereunder automatically will conclude on December 31, 2026, unless it is sooner terminated pursuant to this Section 5. You and the Company agree that (a) you may resign for any reason at any time and (b) the Company may terminate your employment for any reason at any time. Additionally, if you shall fail to be

reelected as a director of the Company at the Company's 2026 Annual Meeting of Stockholders, your employment hereunder shall automatically terminate as of the date of such meeting (the "Meeting Date"). In such case, 75% of any unvested options held by you as of the Meeting Date shall vest automatically on the Meeting Date and any remaining unvested options shall expire.

6. Withholding. All forms of compensation referred to in this Agreement are subject to applicable withholding and payroll taxes and other deductions required by law. You hereby acknowledge that the Company does not have a duty to design its compensation policies in a manner that minimizes your tax liabilities, and you will not make any claim against the Company or its Board related to tax liabilities arising from your compensation.
7. Expenses. You shall be entitled to reimbursement of reasonable expenses incurred by you in connection with your employment in accordance with the Company's Travel and Entertainment Policy.
8. Confidential Information. In connection with your entry into this Agreement, you will enter into agreement(s) with the Company regarding obligations of non-competition, non-solicitation, confidentiality and assignment of intellectual property (the "Confidentiality Agreement").
9. Interpretation, Amendment, Waiver and Enforcement. This Agreement, the Confidentiality Agreement, any equity award agreements entered into between you and the Company, and any agreements concerning your rights to defense and indemnification constitute the complete agreement between you and the Company, contain all of the terms of your employment with the Company and supersede any prior agreements, representations or understandings (whether written, oral or implied) between you and the Company. This Agreement may not be modified or amended, and no breach will be deemed to be waived, unless agreed to in writing by you and a duly authorized officer or the lead independent director of the Board of the Company. No delay or omission by the Company in exercising any right under this Agreement shall operate as a waiver of that or any other right. A waiver or consent given by the Company on one occasion shall be effective only in that instance and shall not be construed as a bar to or waiver of any right on any other occasion. The terms of this Agreement and the resolution of any disputes as to the meaning, effect, performance or validity of this Agreement or arising out of, related to, or in any way connected with, this Agreement, your employment with the Company or any other relationship between you and the Company (the "Disputes") will be governed by Massachusetts law, excluding laws relating to conflicts or choice of law. You and the Company submit to the exclusive personal jurisdiction of the federal and state courts located in the Commonwealth of Massachusetts in connection with any Dispute or any claim related to any Dispute, and the prevailing party shall be awarded its attorneys' fees and costs.
10. Return of Company Property. Upon the end of the Term, you shall promptly return to the Company any keys, passes, computer equipment, or other property belonging to the Company, and, unless your service as a director of the Company shall continue following the Term, you shall also return all writings, files, records, correspondence, notebooks, notes and other documents and items (including any copies thereof) containing confidential information or relating to the business or proposed business of the Company or its Affiliates or containing any trade secrets relating to the Company or its Affiliates. For purposes of the preceding sentence, the term "trade secrets" shall have the meaning ascribed to it under the Uniform Trade Secrets Act. If requested, you agree to represent in writing to the Company that you have complied with the foregoing provisions of this Section.
11. Assignment. Neither you nor the Company may make any assignment of this Agreement or any interest in it, by operation of law or otherwise, without the prior written consent of the other; provided, however, that the Company may assign its rights and obligations under this Agreement without your consent to one of its Affiliates or to any person with whom the Company will hereafter effect a reorganization, consolidate with, or merge into or to whom it transfers all or substantially all of its properties or assets. This Agreement will inure to the benefit of and be binding upon you and the Company, and each of our respective successors, executors, administrators, heirs and permitted assigns.
12. Miscellaneous. The headings and captions in this Agreement are for convenience only and in no way define or describe the scope or content of any provision of this Agreement. The words "include" "includes" and "including" when used herein will be deemed in each case to be followed by the words "without limitation." The term "Affiliates" means all persons and entities directly or indirectly controlling, controlled by or under common control with the Company, where control may be by management authority, equity interest or otherwise. This Agreement may be executed in two or more

counterparts, each of which will be an original and all of which together will constitute one and the same instrument. As with all employees, our offer to you is contingent on your submission of satisfactory proof of your identity and your legal authorization to work in the United States.

We appreciate your service to the Company.

Sincerely,

QUANTERIX CORPORATION

Jeffrey T. Elliot
Lead Independent Director

AGREED TO AND ACCEPTED

/s/ Bill Donnelly

William P.
Donnelly

Date:
11/22/2025

Amended and Restated Quanterix Corporation Non-Employee Director Compensation Policy

Effective as of January 1, 2025

I. Overview

The Board of Directors (the “Board”) of Quanterix Corporation (the “Company”) has approved this Amended and Restated Non-Employee Director Compensation Policy (the “Policy”) to provide an inducement to attract and retain the services of qualified persons to serve as directors.

II. Eligibility

This Policy shall apply to each director of the Board who is not an employee of, or compensated consultant to, the Company or any of its Affiliates (as defined in the 2017 Employee, Director and Consultant Equity Incentive Plan, or any successor plan (“the Plan”)) (a “Non-Employee Director”). Employees of the Company and their affiliates are not eligible to receive compensation under this Policy.

III. Director Compensation

The following is a description of the compensation arrangements under which our Non-Employee Directors are compensated for their service as directors, including as members of the various committees of our Board, consisting of the cash retainers described in Section III.A and the equity awards described in Section III.B.

A. Cash Compensation

1. *Terms for Cash Payment*

Subject to Section III.A.2, each Non-Employee Director shall receive the following annual cash compensation for his or her service on the Board and committees of the Board:

Base Board Retainer	\$50,000
Additional Lead Director/Non-Employee Board Chairman Retainer	\$45,000
Additional Audit Committee Chairman Retainer	\$20,000
Additional Compensation Committee Chairman Retainer	\$15,000
Additional Nominating and Governance Committee Chairman Retainer	\$10,000
Additional Audit Committee Member Retainer	\$10,000
Additional Compensation Committee Member Retainer	\$7,500
Additional Nominating and Governance Committee Member Retainer	\$5,000

Cash payments to Non-Employee Directors shall be paid quarterly in arrears on the first Company payroll date, or as soon as administratively practicable, following the end of the fiscal quarter to which service relates (each, a “Payment Date”).

Each Non-Employee Director: (i) who is elected or appointed to the Board after the date hereof or (ii) ceases to be a Non-Employee Director during a fiscal quarter, shall receive a prorated cash retainer for the portion of such partial fiscal quarter during which he or she served on the Board or a committee of the Board (the "Prorated Retainer"). The Prorated Retainer shall be an amount equal to the product of (A) the aggregate amount payable in respect of such Non-Employee Director's service for a full fiscal quarter multiplied by (B) a fraction, the numerator of which is (x) the number of days during such fiscal quarter which the Non-Employee Director served on the Board or committees, and the denominator of which is (y) the total number of days during such fiscal quarter. The Prorated Retainer shall be paid on first Payment Date following such fiscal quarter.

2. Election for Equity in Lieu of Cash Retainers

Prior to the end of each calendar year, each Non-Employee Director shall make an annual election by delivery to the Company of an election form, substantially in the form attached hereto as Exhibit A (the "Election Form"), with respect to cash retainers for the following calendar year, indicating whether he or she elects to receive the retainers in cash, as described in Section III.A.1, or in the Company's common stock, \$0.001 par value per share ("Common Stock"), in lieu of the cash retainers. If no election has been made as of the first day of the year, the Non-Employee Director shall receive all retainers in cash as set forth in Section III.A.1 or, if a previous election has been made to receive Common Stock in lieu of the cash retainers, such election shall remain in effect for subsequent calendar years until such election is changed by the completion, signature and delivery to the Company of a new Election Form, in accordance with the terms of this Policy. Each newly elected or appointed Non-Employee Director shall make an election prior to, or within 30 days of, his or her initial appointment or election to the Board, for the remainder of the year of such appointment or election, whether to receive the retainers in cash or in Common Stock.

In the event an election is made to receive Common Stock in lieu of cash retainers, such director shall automatically be granted, without any further action by the Board, on the first trading day following each fiscal quarter a number of shares of Common Stock having an aggregate fair market value equal to the aggregate amount of such Non-Employee Director's cash retainer for such fiscal quarter, determined by dividing (A) the aggregate amount of the retainers by (B) the Fair Market Value (as defined in the Plan) of the Common Stock on such trading day.

All Common Stock granted to Non-Employee Directors under this Policy shall be granted under the Plan and will be subject to the terms and conditions set forth in the Plan.

B. Equity Compensation

1. Annual Equity Awards

Each Non-Employee Director will automatically be granted, without any further action by the Board, on the first trading day of each fiscal year, an annual equity award valued at \$200,000, rounded to the nearest whole share (the "Annual Award"). The Annual Award is comprised of (A) 60% non-qualified stock options to purchase Common Stock at an exercise price equal to

the Fair Market Value as of such grant date (the “Options”) and (B) 40% restricted stock units, (each RSU relating to one (1) share of Common Stock (the “RSUs”). The number of Options shall be determined on a value basis using the Company’s standard Black-Scholes valuation methodology. The number of RSUs shall be determined by dividing total value of the RSUs by the Fair Market Value (as defined in the Plan) of the Common Stock on such grant date. The Annual Awards shall become vested in full on December 31st of the year in which such awards were granted, provided that the Non-Employee Director is a director of the Company on the applicable vesting date.

2. Initial Equity Awards for Newly Elected Directors

Upon initial election or appointment of a Non-Employee Director to the Board, such Non-Employee Director will automatically be granted, on his or her election or appointment date, without any further action by the Board, an award valued at \$400,000, rounded to the nearest whole share (the “Initial Award”). The Initial Award is comprised of (A) 60% Options at an exercise price equal to the Fair Market Value as of such grant date and (B) 40% RSUs. The number of Options shall be determined on a value basis using the Company’s standard Black-Scholes valuation methodology. The number of RSUs shall be determined by dividing total value of the RSUs by the Fair Market Value (as defined in the Plan) of the Common Stock on such grant date. The Options granted pursuant to Initial Awards shall vest over a three-year period, with one-third vesting on the first anniversary of the applicable grant date, and the remainder vesting over the following two years in 24 successive equal monthly installments at the end of each month until the third anniversary of such grant date, provided that the Non-Employee Director is a director of the Company on the applicable vesting date. The RSUs granted pursuant to Initial Awards shall vest over a three-year period, with one-third vesting on each of the first, second, and third anniversaries of the applicable grant date, provided that the Non-Employee Director is a director of the Company on the applicable vesting date.

All Annual Awards and Initial Awards granted to Non-Employee Directors under this Policy shall be granted under the Plan, and will be subject to the terms and conditions set forth in the Plan, and the form of Stock Option Agreement and form of Restricted Stock Unit Agreement, each as approved by the Board.

C. Expense Reimbursement

Upon presentation of documentation of such expenses reasonably satisfactory to the Company, each Non-Employee Director shall be reimbursed for his or her reasonable out-of-pocket business expenses incurred in connection with attending meetings of the Board and its committees or in connection with other business related to the Board. Each Non-Employee Director shall also be reimbursed for his or her reasonable out-of-pocket business expenses authorized by the Board or one of its committees that are incurred in connection with attendance at meetings with the Company’s management. Each Non-Employee Director shall abide by the Company’s travel and other policies applicable to Company personnel.

IV. Policy Review / Amendments

The Compensation Committee or the Board shall review this Policy from time to time to assess whether any amendments in the type and amount of compensation provided herein should be adjusted in order to fulfill the objectives of this Policy. This Policy may only be amended by the Board.

EXHIBIT A

QUANTERIX CORPORATION

NON-EMPLOYEE DIRECTOR COMPENSATION ELECTION FORM

In accordance with the Amended and Restated Director Compensation Policy (the "Policy"), of Quanterix Corporation (the "Company"), the undersigned hereby makes the following election for the period from January 1, ____ through December 31, ____ (the "Period") with respect to the non-employee director board and committee cash retainers (the "Retainers") to be earned by the undersigned during the Period:

I elect to receive my Retainers during the Period (please check one of the following):

_____ in cash

_____ in the Company's common stock, \$0.001 par value per share ("Common Stock")

In accordance with the Policy, Common Stock shall be granted quarterly in arrears on the first trading day following the end of the fiscal quarter to which service relates and the undersigned shall be granted automatically and without any further action required by the Board of Directors ("Board") under the Company's 2017 Employee, Director and Consultant Equity Incentive Plan or any successor plan (the "Plan") a number of shares of Common Stock having an aggregate fair market value equal to the aggregate amount of the Retainers for such fiscal quarter, determined by dividing (A) the aggregate amount of the Retainers by (B) the Fair Market Value (as defined in the Plan) of the Common Stock on such trading day (rounded down to the nearest whole share), in lieu of the aggregate amount of the Retainers that would otherwise be paid in cash in respect of such fiscal quarter.

Signature

Print Name

Date

Manufacturing Agreement

Whereas, Columbia Electrical Contractors, Inc. d/b/a Columbia Tech, a Massachusetts corporation with a principal place of business of 27 Otis Street, Westborough, MA 01581 (“Columbia Tech”), provides contract manufacturing services;

Whereas, Akoya Biosciences, Inc., a Delaware corporation with a place of business at 68 Elm Street, Hopkinton, MA 01748 (“Customer”), desires to have products they designed manufactured by another party; and

Whereas, Columbia Tech and Customer (collectively “Parties” or individually “Party”) wish to enter an agreement whereby Columbia Tech provides contract manufacturing services to Customer.

Now, therefore, Columbia Tech and Customer agree as follows, effective as of July 17, 2019 (“Effective Date”):

1. ORDERS:

- a. Quotations. Customer may from time to time request that Columbia Tech provide pricing for certain goods (the “Products”) based on minimum order sizes in amounts designated by Customer (“Order Sizes”). In any such request, Customer shall also designate the applicable quantity, if any, of each material that it intends to provide Columbia Tech pursuant to Addendum 1 for use in manufacturing such Products (the “Consigned Materials”). Columbia Tech shall thereafter provide a quotation for the manufacture of Products at each Order Size when using the Consigned Materials (the “Quotation”). Customer may request changes to the Quotation based on different Order Sizes, different amounts of Consigned Materials or other changes, and Columbia Tech will use reasonable efforts to provide new Quotations. All Quotations shall (i) list any requirements regarding Consigned Materials, Order Size, lead time for initial delivery and any other similar requirements to achieve the price stated therein, and (ii) show any costed bill of materials reflecting Markup Percentages, as may be updated pursuant to this paragraph, that Columbia Tech uses to arrive at the Price. The Quotation shall clearly identify the document by a unique number (the “Quotation Number”) and list the expiration date for the Quotation. The initial “Markup Percentages” shall be included in the Initial Quotation. Thereafter, Columbia Tech may change the Markup Percentages once every three months by providing written notice to Customer. All such changes shall apply to new quotations and not previously accepted orders.
- b. Orders. For any Quotation that has not expired, Customer may from time to time submit one or more purchase orders to Columbia Tech, each of which includes the Quotation Number, the total number of Products to be ordered, and a schedule of when to deliver Products pursuant to the order, provided that no such schedule

shall exceed any lead time needed to acquire any necessary materials and build the Product, as such Lead time may change from time to time. Any other terms and conditions appearing on the purchase order shall be null and void, placed for convenience only, and the terms and conditions of this Agreement and the terms shown on the Quotation shall control (collectively the “Requested Order”).

- c. Provided that: (i) the Requested Order conforms to the Quotation and the requirements of Section 1.b above and does not otherwise create confusion on the obligation of the parties; and (ii) the schedule of deliveries (each applicable date of delivery, a “Delivery Date”) does not have a Delivery Date more than one (1) year after the date of the Requested Order (collectively, the “Order Requirements”), then Columbia Tech shall accept the Requested Order. Should any Requested Order not conform to the Order Requirements, Columbia Tech may reject the Requested Order by providing notice describing the non- conformity in reasonable detail, provided that Columbia Tech shall thereafter work diligently with Customer to agree as soon as practicable on a modified Requested Order that conforms to the Order Requirements. Any Requested Order not rejected within ten (10) days shall be deemed to be accepted. Any accepted Requested Order shall hereinafter be referred to as an “Order”.

2. SPECIFICATIONS: Prior to Columbia Tech’s manufacture of any Product, Customer shall provide Columbia Tech with all material instructions, drawings, specifications, tests, procedures, processes, and other information, including packaging requirements and storage instructions, necessary to manufacture the Products pursuant to any Order (collectively the “Specifications”). All Specifications shall be clear and allow the manufacture of the Products without reference to any outside resources, other than manufacturing processes or other know-how generally known by contract manufacturers in the industry or specifically known by Columbia Tech. Columbia Tech will not be responsible for performing any action or have any duty to Customer except: (a) what expressly appears in the Specifications, or (b) appears in this Agreement; and then only such actions that are capable of reasonably being performed by a similarly situated contract manufacturer and that are not contrary to any law, rule or regulation.

3. DELIVERY:

- a. The Delivery Dates on the Order may not be changed except as described in this Section 3.a. Customer shall have the right, upon notice to Columbia Tech, to push out the Delivery Date for any unit of Product, one time, to a later date that is not more than a maximum of thirty (30) days after the original Delivery Date for such unit as set forth in the Order, and thereupon such later date shall be deemed the Delivery Date for such unit for all purposes hereunder. In addition, Customer shall have the right, upon notice to Columbia Tech, to cause Columbia Tech to store any Product (“Stored Product”) at Columbia Tech’s facility after the applicable Delivery Date until such time as Customer orders actual delivery of such Stored Product (the date for which Customer orders actual delivery of

such Stored Product, the “Stored Product Delivery Date”), with the applicable charges for such storage set forth in Attachment C. In such event, Columbia Tech shall store the Stored Product in accordance with any applicable Specifications and shall keep Customer’s title to the Stored Product free and clear of all liens and encumbrances.

- b. Columbia Tech may not push out any Delivery Date or Stored Product Delivery Date, however, Columbia Tech shall be granted a grace period of fifteen (15) days after any Delivery Date (or, for delivery of Stored Product, 5 days after any Stored Product Delivery Date) to deliver the applicable Product(s) without penalty or breach of this Agreement (the “Grace Period”). If Columbia Tech fails to deliver any Product or Stored Product within the applicable Grace Period, then, unless such failure is due to late delivery to Columbia Tech, for reasons that are beyond Columbia Tech’s control, of materials that (i) are on allocation by the applicable supplier, (ii) are, to Columbia Tech’s knowledge after reasonable inquiry, only available from one source, or (iii) required by Customer to be supplied from a supplier designated by Customer (any such failure, an “Excused Late Delivery”), Columbia Tech shall provide an invoice credit of one percent (1%) for each week that the delivery is late, up to a maximum of five percent (5%) credit (the “Late Delivery Credit”). Subject to Customer’s additional termination rights and obligations in Section 19, the Late Delivery Credit shall be Customer’s sole remedy for any late delivery by Columbia Tech.
- c. All delivery of Products by Columbia Tech shall be made EX Works (EXW Incoterms 2010) Columbia Tech’s dock in Worcester County, MA. For Products other than Stored Products, title and risk of loss shall pass to Customer upon such delivery. For Stored Products, title shall pass to Customer upon the earlier of such delivery (based on the Stored Product Delivery Date) or Customer’s payment therefor, and risk of loss shall pass to Customer upon such delivery (based on the Stored Product Delivery Date). In addition and without limiting the previous statement, after Columbia Tech has placed the Products with a freight carrier, Columbia Tech not carry any insurance on the Products for delivery or storage. Customer shall maintain all such insurance. Should any change occur in Incoterms, then the 2010 version shall still control.
- d. Notwithstanding the previous section, if Customer provides a freight carrier account number that is billed to the Customer and the choice of freight carrier on the Order (“Delivery Number”), then Columbia Tech will arrange for shipment using Customer’s Delivery Number.
- e. Customer or Customer’s designee shall have five (5) days after any shipment of Product has been delivered by the freight carrier to its listed destination to accept such Product. After such five (5) day period, the Product will be deemed to have been accepted, and such acceptance may not be revoked. The five (5) day period for acceptance shall not increase the warranty term, which shall begin to run on

actual delivery pursuant to Section 3.c based on the Delivery Date or Stored Product Delivery Date, as applicable. "Accepted" shall have the same meaning of acceptance pursuant to the Uniform Commercial Code.

- f. For any commercial invoice prepared by Customer, which involves any shipment that originates from Columbia Tech, Customer shall include the following statement on the commercial invoice:

These items are controlled by the U.S. Government and authorized for export only to the country of ultimate destination for use by the ultimate consignee or end-user(s) herein identified. They may not be resold, transferred, or otherwise disposed of, to any other country or to any person other than the authorized ultimate consignee or end-user(s), either in their original form or after being incorporated into other items, without first obtaining approval from the U.S. government or as otherwise authorized by U.S. law and regulation.

- g. Each Party agrees to abide by any and all applicable laws or regulations relating to the export of products, including, but not limited to, the Export Arms Regulations and International Traffic in Arms Regulations. Customer shall promptly provide Columbia Tech with any applicable Export Control Classification Number upon Columbia Tech's request.

4. PRICE:

- a. The purchase price to be paid for the Products shall be determined using the applicable Quotation referenced on the Order ("Invoice Price"). The Credit Amount for any Consigned Materials provided by Customer shall be provided on a monthly basis based on materials actually used and at the price shown on an SMI Order (as such term is defined in this Agreement).
- b. Columbia Tech may arrive at prices based on reasonable criteria of its choosing; provided, however, that in the event that cost savings are identified based on lower costs from suppliers, Columbia Tech shall apply fifty percent (50%) of any such cost savings to reducing the applicable components of its pricing formula, with any such price reduction applying to any invoice for any Product that uses or incorporates the applicable lower-cost materials; and any change in tariffs, duties or other material costs may be immediately passed on to the Customer.

5. PAYMENTS/SECURITY INTEREST:

- a. Customer shall pay to Columbia Tech a deposit not greater than Columbia Tech's estimated invoice costs for any materials for such Order; provided, however, that the Credit Amount for any Consigned Materials provided by

Customer to Columbia Tech in accordance with a delivery schedule on the SMI Order based on the dates Columbia Tech would reasonably need such Consigned Materials for use in manufacturing Products under such Order (up to the full amount of such deposit requirement) shall be credited against such deposit requirement (and Customer shall not be required to pay any part of the deposit in cash so long as (1) such delivery schedule contemplates eventual delivery of Consigned Materials having an aggregate Credit Amount at least equal to the amount of the deposit requirement and (2) Customer delivers Consigned Materials in accordance with such delivery schedule). Columbia Tech shall invoice Customer for the remaining balance of the Invoice Price, upon (i) for Products other than Stored Products, the later of (x) the Delivery Date for such quantity or (y) actual delivery of such quantity pursuant to Section 3.c, and (ii) for Stored Products, the Delivery Date for such quantity. Customer shall pay any undisputed amount set forth in any such invoice within forty-five (45) days after the date of such invoice. Notwithstanding the above, invoices shall only contain credits for cash deposit amounts, if any, and Customer may use any credit on its account based on any consumed Consigned Materials by providing written notice to Columbia Tech at the time of payment.

- b. Columbia Tech shall invoice Customer for any storage fees with respect to Stored Product monthly in arrears. Customer shall pay any undisputed amount set forth in any such invoice within forty-five (45) days after the date of such invoice.
- c. Any undisputed amount which is not paid when due shall bear interest at the rate of one and one half percent (1.5%) per month (or if that amount is prohibited by law, then at the maximum interest rate allowable by law) or portion thereof from the date such amount became due through the date on which payment is received by Columbia Tech. All payments shall be made in US Dollars to Columbia Tech at 27 Otis Street, Westborough, MA 01581.

6. [Reserved]

7. WARRANTIES:

- a. Workmanship. Subject to the limitations included herein and within the Warranty Period, Columbia Tech warrants to the Customer that (i) all Products shall conform to the Specifications and free from defects in workmanship and (ii) Columbia Tech's manufacture of the Products shall conform to the Specifications.
- b. Materials. Subject to the limitations included herein and within the Warranty Period, Columbia Tech shall on behalf of the Customer use all reasonable efforts to enforce warranties Columbia Tech obtains from its vendors and suppliers on the materials in the Products ("Vendor Warranties"). Columbia Tech shall use reasonable efforts to notify Customer in writing prior to ordering any material for use in manufacturing Products if such material is not covered by a warranty

from the applicable vendor or supplier against defects in materials or workmanship for a period of at least 13 months after date on which delivery of the applicable Product to Customer is expected to be made; provided that any breach of this notification shall not subject the Products to any additional warranty remedy. Columbia Tech shall have no liability on the materials except as set forth in this Section 7.b.

- c. Title. Columbia Tech warrants that, upon delivery, title to the Products shall pass to the Customer free and clear of any liens or encumbrances. Nothing in this Section 7.c shall affect Customer's title to the Consigned Materials as provided in Addendum 1.
- d. Intellectual Property. Columbia Tech warrants that its manufacturing processes will not infringe any third party's intellectual property rights. Columbia Tech makes no warranty on any intellectual property right of the customer in any assembly, process, design, product, expression or article of manufacture supplied by the Customer, contained in the Specifications, or arising specifically in result of compliance with the Specifications.
- e. Preconditions to Warranty Repair. Customer shall be responsible for identifying, and communicating to Columbia Tech the cause of failures prior to Columbia Tech performing any warranty work. If the defect is reported during the Warranty Period, Customer shall immediately thereafter request return, repair, replace, refund or credit instructions in accordance with Section 7.h (the "RMA instructions").
- f. [Reserved]
- g. Warranty Period. Subject to the limitations included herein, Columbia Tech's warranties pursuant to Section 7.a shall survive as to any Product for a period of thirteen (13) months after delivery of such Product (for clarity, in case of Stored Product, for a period of thirteen (13) months after actual delivery based on the Stored Product Delivery Date). Columbia Tech will assist with enforcing Vendor Warranties pursuant to Section 7.b for the remaining duration of any Vendor Warranties (collectively this section defines the "Warranty Period").
- h. Remedies for Warranty. For any Product not conforming to the warranty set forth in Section 7.a, promptly when notified of the nonconformity by the Customer within the Warranty Period, Columbia Tech shall issue a return material authorization (an "RMA") for the repair or replacement of such non-conforming Product by Columbia Tech. Upon receiving such non-conforming Product pursuant to the RMA, Columbia Tech shall repair or replace such non-conforming Product, and re-deliver such repaired or replacement Product to its applicable destination, as soon as possible, but subject to material availability. Customer shall bear all costs of shipping and returning any Products to Columbia Tech. If Customer follows all RMA instructions for warranty repair, the warranty

repair was due to Columbia Tech's breach of warranty during the warranty period (and not on Columbia Tech providing assistance on Vendor Warranties), and the shipping is from a location in the contiguous forty-eight states, Columbia Tech will refund Customer's cost of shipping the product, and pay for return shipping to a location in the contiguous forty-eight states.

- i. Exclusions and Limitations. The warranties in Section 7.a do not cover defects caused by materials, services, designs, specifications, instructions, tests or information supplied by the Customer or failures which result from accident, misuse, abuse, neglect, mishandling, alteration or modification by anyone other than Columbia Tech or its employees, agents or subcontractors. , CUSTOMER'S SOLE REMEDY FOR ANY WARRANTY CLAIM UNDER SECTION 7.a SHALL BE THE REMEDIES SET FORTH IN SECTION 7.h. THE WARRANTIES SET FORTH HEREIN ARE EXCLUSIVE AND IN LIEU OF ALL OTHER WARRANTIES (WHETHER EXPRESS OR IMPLIED) INCLUDING, WITHOUT LIMITATION THE WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE CUSTOMER MAY NOT ASSIGN THIS WARRANTY TO ANY OTHER PARTY WITHOUT THE EXPRESS WRITTEN PERMISSION OF COLUMBIA TECH. COLUMBIA TECH OFFERS NO WARRANTY WHATSOEVER ON OTHER SERVICES, INCLUDING, BUT NOT LIMITED TO SERVICES FOR NON-RECURRING ENGINEERING, RECURRING COSTS OR COBUILDS. **UNDER NO CIRCUMSTANCES SHALL COLUMBIA TECH BE RESPONSIBLE FOR THE RECALL OF ANY PRODUCT.**
- j. Each Party represents and warrants: (i) that such Party has the requisite power and authority to enter into and perform this Agreement and all necessary rights to grant any rights or licenses purported to be granted by it herein; (ii) that this Agreement has been duly authorized by all necessary action on the part of such Party; (iii) that the execution, delivery and performance of this Agreement by such Party shall not conflict with, breach or result in a violation of such Party's organizational documents or any judgment, order or decree of any court or arbiter or contract or agreement to which such Party is bound; (iv) that this Agreement constitutes the valid and binding obligation of such Party and is enforceable against such Party in accordance with its terms, subject to applicable bankruptcy, insolvency, reorganization, creditors' rights and other similar laws; and (v) it has all governmental approvals required to enter into this Agreement and to perform its obligations hereunder.

8. INVENTIONS/PATENT RIGHTS: Unless modified by a separate addendum, as between the parties, (a) Customer shall own any enhancements to Customer's intellectual property created by Columbia Tech or its subcontractors performing assembly-related activities hereunder (but not other third parties) pursuant to this Agreement ("Inventions"), and (b) Columbia Tech shall own any enhancement to Columbia Tech's manufacturing processes, other than any such enhancement that qualifies as an Invention. Nothing herein shall create any obligation for Columbia Tech to obtain title on behalf of Customer to any Inventions created by third parties related to the manufacture of Customer's products; provided, however, that Columbia Tech shall ensure that any subcontractor performing assembly-related activities hereunder assigns to Akoya (or to Columbia Tech for subsequent assignment by Columbia Tech to Akoya) any rights that such subcontractor may have in or to any Inventions or Customer Records created by or on behalf of such subcontractor in connection with such activities. At Customer's request, sole cost and expense, Columbia Tech will perform all lawful and necessary acts, sign all patent, trademark and copyright applications, oaths, assignments and other papers necessary to apply for, obtain and assign to Customer and enforce the patents, trademarks and copyrights for any and all such Inventions. Columbia Tech's duty to enforce any intellectual property right shall only include responding to reasonable requests from Customer based upon Customer's actions in policing Customer's own intellectual property. Columbia Tech shall not have any independent responsibility or duty to police any patent or initiate any legal action against infringers of any intellectual property. Customer shall conduct and pay for all costs, expenses and other aspects of any patent, trademark or copyright application, prosecution, renewal, or policing process, including but not limited to, the assignment of intellectual property to Customer, all searches of intellectual property, attorney's fees and all other expenses incident thereto. Columbia Tech shall reasonably assist Customer in all such activities provided that, in addition to the any other fees or amounts owed pursuant to this Agreement, Customer shall reimburse Columbia Tech for any out-of-pocket costs incurred by Columbia Tech in connection with such assistance.

9. RELEASE/INDEMNIFICATION:

- a. Customer shall indemnify and defend Columbia Tech and its shareholders, directors, officers, employees, Affiliates and agents from and against any and all claims, actions, suits, proceedings, obligations, liabilities, orders, damages, government inquiries and investigations of any nature (including, but not limited to subpoenas, expressions of interest, audits and all other phases of inquiries or investigations), and costs and expenses (including attorneys' and paralegals' fees and expenses) (collectively "Losses"), in connection with any third party claim or government inquiry or investigation ("Third Party Claim"), regardless of whether such Loss(es) are based on contract (including indemnification regardless of whether such claim is due to Columbia Tech providing express or implied indemnification to any third party including third party testing or certification companies), tort (including negligence, regardless of whether the negligence is by Columbia Tech or its Affiliates acting alone, Columbia Tech or its Affiliates

acting with others, the negligence of Customer, or the negligence of third parties), strict liability, or any other theory or form of action, even if such Party has been advised of the possibility thereof, arising out of, in connection with, or incidental to: (a) Customer's breach of this Agreement, (b) Customer's negligence, willful misconduct, gross negligence or unlawful acts whether through action or inaction or (c) any service, manufacture, sale, or use of any Product by or on behalf of Customer or its customers, or (d) the Consigned Materials, all of which includes, but is not limited to: (i) any violation or infringement upon the rights of any individual or entity, whether based on patents, trademarks, trade designs, copyrights, trade secrets or otherwise; and (ii) any bodily injury, property damage, loss of use of a product, incidental damages and consequential damages; provided, however, that Customer shall have no such obligation with respect to any Losses to the extent arising out of, in connection with, or incidental to any gross negligence or willful misconduct of Columbia Tech or its employees or agents or any infringement to the extent based on Columbia Tech's manufacturing process (excluding any manufacturing process contained in the Specifications or arising specifically as a result of compliance with the Specifications). Additionally, if the indemnified event is based on Columbia Tech's breach of this Agreement or the negligence (other than gross negligence) of Columbia Tech, Customer shall be responsible for such indemnification only for the amount of such Losses, if any, that exceeds the Cap.

- b. Columbia Tech shall indemnify and defend Customer and its shareholders, directors, officers, employees, Affiliates and agents from and against any and all Losses in connection with Third Party Claims arising out of, in connection with, or incidental to (a) Columbia Tech's breach of this Agreement, (b) Columbia Tech's negligence, willful misconduct, gross negligence or unlawful acts whether through action or inaction or (c) infringement of a third party's intellectual property right based on Columbia Tech's manufacturing process (excluding any manufacturing process contained in the Specifications or arising specifically as a result of compliance with the Specifications), except that, for any Losses arising from Columbia Tech's negligence or any breach of the Agreement by Columbia Tech, Columbia Tech shall not be responsible for such indemnification to the extent such Losses exceed the Cap; provided, however, that Columbia Tech shall have no such obligation with respect to any Losses to the extent arising out of, in connection with, or incidental to any breach, negligence, gross negligence or willful misconduct of Customer or its employees or agents.
- c. Any indemnification claim under Section 9.a or 9.b in respect of a Party or its affiliated persons shall be asserted solely by such Party ("Indemnitee") in accordance with this paragraph. Indemnitee shall provide the other Party ("Indemnitor") with prompt written notice of the applicable Third Party Claim; provided, however, that failure to provide such notice shall not relieve Indemnitor of its obligations hereunder except to the extent that Indemnitor can demonstrate

prejudice attributable to such failure. If Indemnitor acknowledges its indemnification obligation in writing, then defense of the applicable Third Party Claim shall be provided by legal counsel chosen by the Indemnitor and reasonably acceptable to Indemnitee; provided, however, that the Indemnitor shall not have the right to assume or continue control of the defense of such Third Party Claim, if (i) such Third Party Claim seeks non-monetary relief (in whole or in part) or relates to or arises in connection with any criminal proceeding, (ii) the named parties in any such action (including any impleaded parties) include both the Indemnitee and the Indemnitor (or their respective Affiliates) and the representation of both parties by the same counsel would be inappropriate due to actual or potential differing or conflicts of interests between them, or (iii) the Indemnitor fails to actively and diligently conduct the defense of such Third Party Claim; provided, further, that, if Indemnitor assumes the defense of such Third Party Claim, Indemnitee shall have the right to retain separate counsel at its own expense to participate in such defense. Without the prior written consent of Indemnitee, Indemnitor may not settle any such third party claim unless the sole relief is money damages that are paid in full by Indemnitor and such settlement includes a release of Indemnitee (or its applicable affiliated person) from all liability for such claim or a dismissal of litigation with prejudice.

- d. "Affiliates" shall mean any corporation, limited liability company, trust, partnership, or any other form of business that is under the same common control of a Party, including but not limited to any company having the same shareholder(s) as such Party.

10. SAFETY/WARNINGS /COMPLIANCE WITH LAWS/QUALITY:

- a. Subject to Columbia Tech's warranties and obligations hereunder, Customer accepts and assumes all liability and responsibility for any and all safety testing, labeling and warnings in connection with any of the Products manufactured by Columbia Tech pursuant to any Order. Subject to Columbia Tech's warranties and obligations hereunder, Customer agrees to take any and all actions necessary or desirable actions to ensure that the Products manufactured by Columbia Tech pursuant to any Order and the use thereof comply with all federal, state or local laws, including any regulatory requirements. Columbia Tech's performance of any regulatory duty or function shall not waive or change the duties and obligations of the Parties pursuant to this section.
- b. Columbia Tech shall perform its activities hereunder in compliance with (i) all applicable laws, rules or regulations of any governmental authority having authority over Columbia Tech or its activities under this Agreement (for clarity, excluding any laws, rules or regulations applicable to the marketing, sale or distribution of the Products to end users), and (ii) if the parties enter into a written quality agreement relating to their activities under this Agreement ("Quality Agreement"), any applicable requirements set forth in the Quality Agreement.

- c. For the term of this Agreement and for 5 years thereafter, Columbia Tech shall keep, complete and accurate records with respect to the manufacture of Products and related activities in sufficient detail to demonstrate compliance with their obligations under this Agreement. At reasonable times and upon reasonable notice, Columbia Tech shall make such records available for inspection, examination and copying by Customer or its representatives. Columbia Tech agree that all records specifically related to activities for Customer in connection herewith (“Customer Records”) shall be owned by Customer and delivered to Customer upon its request.
 - d. At reasonable times and upon reasonable notice, Columbia Tech shall permit Customer or its representative to observe and inspect Columbia Tech’s applicable facilities and processes and the performance of its activities hereunder, and ensure that its suppliers and subcontractors permit Customer or its representatives to observe and inspect their applicable facilities and process and the performance of their activities in connection herewith, to verify compliance with the terms of this Agreement. Columbia Tech shall reasonably cooperate, with any such inspection, including by making personnel reasonably available to provide any information or assistance reasonably requested by Customer or its representative in connection therewith.
 - e. The Parties shall enter into the Customer Consigned Materials Agreement attached hereto as Addendum 1 (“CCMA”). Without limiting any other provisions of this Agreement, each Party shall timely perform all of its obligations under the CCMA and the Quality Agreement (if any).
11. INSURANCE: Customer shall maintain in full force and effect necessary insurance to operate the Customer. Columbia Tech will maintain general liability insurance in the amount of the Cap. Without limiting the foregoing, Customer shall maintain insurance policies set forth below, at its sole cost and expense, written by companies with an A.M. Best’s Insurance Guide rating of A or better, (i) Product Liability Insurance in the amount of two million dollars (\$2,000,000) (ii) General Liability in the amount of Two million dollars (\$2,000,000). Customer’s Commercial General Liability and Product Liability Insurance policies shall include Columbia Tech, its directors, officers, and employees as Additional Insureds thereunder. Each such policy shall waive or otherwise prohibit insurer subrogation against Columbia Tech and all such other Additional Insureds. Each such policy shall also include a severability of interests (or “separation of insureds”) provision. The Additional Insured status under Customer's Commercial General Liability Policy shall be provided pursuant to an Insurance Services Office (ISO) form CG 20 10 11 85 Additional Insured Endorsement, by a combination of ISO forms CG 20 10 and CG 20 37, or by other equivalent Additional Insured endorsement form(s) that provide both Premises and Operations and Completed Operations liability coverage that is at least as broad as that afforded by the above-referenced forms. With respect to liability of Customer under this Agreement, all insurance required of Customer hereunder shall

respond on a primary (not excess or contributory) basis with respect to any similar insurance maintained by Columbia Tech and/or any other party required to be included as an Additional Insured hereunder, notwithstanding any policy language to the contrary.

Prior to the commencement of any performance under this Agreement, Customer will provide Columbia Tech with evidence that the insurance coverage required of Customer hereunder is in full force and effect. In the event any such insurance renews or is terminated during the course of Customer's performance, Customer will promptly provide Columbia Tech with evidence that such coverage will be renewed or replaced upon termination with insurance that complies with these provisions. Such evidence of insurance will be in the form of a standard Certificate of Insurance or other form of evidence of insurance acceptable to Columbia Tech, and shall contain sufficient information to allow Columbia Tech to determine whether there is compliance with these provisions. Such evidence of insurance shall be accompanied by copies of any Additional Insured endorsements or automatic Additional Insured policy provisions necessary to achieve compliance with the Additional Insured requirements of this Agreement. All such evidence of insurance shall require that the insurer provide at least thirty (30) day written notice to Columbia Tech prior to the effective date of policy cancellation (ten (10) day notice in the event of cancellation due to nonpayment of premium). All policies shall remain in effect for at least as long as the expected useful life of the Product plus an additional two (2) years. Nothing in this section shall serve as a limit to Customer's liability, and the failure of Customer to obtain insurance or of Customer to enforce the obligations of this section shall also not limit Customer's liability pursuant to any section of this Agreement.

12. WAIVER: Any waiver by any Party of its rights under this Agreement shall be in writing and signed by the Party waiving such right. The failure of either Party to enforce any of the provisions of this Agreement or any rights in respect thereto, or to exercise any election herein provided, shall not waive such provisions, rights or elections or subsequent breaches thereof.
13. RELATIONSHIP OF PARTIES: The relationship of the Parties shall be that of independent contractors and not as partners or joint ventures. Each Party is, and is intended to be, engaged in its own and entirely separate business.
14. ASSIGNMENT/BENEFIT: This Agreement and the rights and obligations in connection herewith may not be assigned by either Party without the prior written consent of the other Party, except that Columbia Tech may subcontract pursuant to the terms of this Agreement, and either Party may assign this Agreement and all of its rights and obligations hereunder pursuant to a merger, sale of all or substantially all of its assets, reorganization or other similar business combination transaction. Without limiting any applicable provisions of the Quality Agreement (if any), upon any subcontract by Columbia Tech of any assembly-related activities, including any assembly that must be produced in a clean room, Columbia Tech will notify Customer within 90 days after the date of the subcontract. Any assignment in contravention hereof shall be null and void.

This Agreement shall be binding upon and inure to the benefit of the Parties and their respective representatives, successors and assigns.

15. NOTICES: All notices and other communications given hereunder shall be in writing and deemed to have been given when (i) personally delivered, (ii) one (1) business day after delivery to a nationally recognized overnight courier service, (iii) upon the written confirmation of receipt following the transmission of a telecopy or email or (iv) three (3) days after being mailed by certified mail, postage prepaid, to the addresses of Columbia Tech or Customer as set forth in the preamble of this agreement or to such other addresses as either Party hereto may provide for such purpose.
16. GOVERNING LAW: This Agreement shall be governed by and construed in accordance with the laws of the Commonwealth of Massachusetts.
17. JURISDICTION AND VENUE: COLUMBIA TECH AND CUSTOMER IRREVOCABLY AGREE THAT ALL ACTIONS OR PROCEEDINGS IN ANY WAY, MANNER OR RESPECT, ARISING OUT OF OR FROM OR RELATED TO THIS AGREEMENT, THE SERVICES, THE PRODUCTS OR ANY DOCUMENT, INSTRUMENT OR TRANSACTION IN CONNECTION HERewith OR THEREWITH SHALL BE HEARD OR LITIGATED EXCLUSIVELY IN COURTS HAVING SITUS WITHIN THE CITY OF BOSTON, COMMONWEALTH OF MASSACHUSETTS. COLUMBIA TECH AND CUSTOMER CONSENT AND SUBMIT TO THE JURISDICTION OF ANY LOCAL, STATE OR FEDERAL COURT LOCATED WITHIN SAID CITY, COUNTY AND STATE AND IRREVOCABLY WAIVE ANY RIGHT TO TRANSFER OR CHANGE VENUE OF ANY SUCH ACTION OR PROCEEDING OR OBJECT TO THE JURISDICTION OF ANY SUCH COURT OVER THE PARTY HERETO.
18. PERSONNEL: Customer shall not solicit or employ the employees of Columbia Tech, for a period of twelve (12) months following the termination of this Agreement; provided, however, that nothing in this Section 18 shall prohibit or restrict Customer from (a) soliciting or employing any person who has ceased to be an employee of Columbia Tech for any reason (other than as a result of solicitation activities of Customer in violation of this Section 18), or (b) responds to any public or general solicitation or otherwise voluntarily contacts Customer regarding employment (other than as a result of solicitation activities of Customer in violation of this Section), or (c) conducting any public or general solicitation activities (including through a bona fide search firm), so long as such activities are not specifically targeted at Columbia Tech employees. Additionally, after termination, if Columbia Tech has hired any person that was employed by Customer or PerkinElmer Health Sciences, Inc. (or any affiliate thereof) on or within 3 months prior to the Effective Date, then Customer may hire also rehire such persons without violating this Section 18.
19. TERM:

- a. This Agreement shall remain in effect until terminated. (i) Either Party may terminate this Agreement for the other Party's material breach by providing the other Party thirty (30) days written notice specifying such material breach in reasonable detail, unless the other Party cures such breach within such thirty (30) day period, and (ii) Customer may terminate this Agreement immediately upon written notice to Columbia Tech, if Columbia Tech fails to deliver any Product under any Order within 30 days after the end of the applicable Grace Period, or fails on 3 or more occasions to deliver any Product under any Order within 15 days after the end of the applicable Grace Period, in either case excluding any Excused Late Delivery ((i) and (ii), collectively, "Termination for Breach"). Additionally, either Party may terminate this Agreement for convenience by providing the other Party ninety (90) days written notice ("Termination for Convenience").
- b. Upon any Termination for Convenience by Columbia Tech, if requested by Customer, Columbia Tech shall, prior to the effective date of such Termination for Convenience, prepare one or more Quotations as requested by Customer and provide Customer with a reasonable opportunity to effect one or more final Orders, all in accordance with all applicable provisions of this Agreement.
- c. Upon any Termination for Convenience, or any Termination for Columbia Tech's Breach by Customer, Columbia Tech shall, to the extent requested by Customer, continue to process and fulfill any Orders accepted and any Requested Orders delivered (and, if applicable, any Orders accepted based on Requested Orders delivered) prior to the effective date of such termination, including any Orders effected pursuant to Section 19.b, all in accordance with all applicable provisions of this Agreement, which shall survive to the extent necessary therefor. Upon a Termination for Customer's Breach by Columbia Tech, Columbia Tech shall have the right to stop processing all orders.
- d. Customer may not alter an order except pursuant to the express terms of this Agreement. In the event of cancellation or termination of any Order except if such Termination of the Order is made accompanying the termination of the Agreement by Customer for Columbia Tech's breach, subject to any obligation to continue to process open orders as may apply pursuant to Section 19.c, Customer shall pay for all finished products, work in process, materials on hand and all costs of cancelling commitments reasonably made by Columbia Tech that are specific to Customer ("Customer-Specific Commitments"), as follows. Customer shall pay finished products at the Invoice Price. Customer shall pay for work in process at Columbia Tech's cost of materials plus thirteen and one half percent (13.5%), and the cost of labor at Columbia Tech's then applicable labor rate plus a markup of three percent on the sum of the labor, material, and material markup ("Interim Amount"), plus an additional markup of seven percent on the Interim Amount, but not more than the Invoice Price. Customer shall pay for all materials on hand and the cost of cancelling any Customer-Specific Commitment at

Columbia Tech's actual cost plus thirteen and one half percent (13.5%), plus a markup of three percent ("Second Interim Amount"), plus an additional markup of seven percent on the Second Interim Amount, but not more than the Invoice Price. In the event of cancellation or termination of any Order that is made in connection with the Termination of this Agreement for breach, subject to any obligation to continue to process open orders as may apply pursuant to Section 19.c, Customer shall pay for all Customer-Specific Commitments as follows. Customer shall pay finished products at the Invoice Price. Customer shall pay for all materials on hand that cannot be returned and all materials that are on order that cannot be cancelled at Columbia Tech's cost. (collectively all the above, the "Termination Payments"). For avoidance of doubt, but without limiting any other rights or remedies that may be available to Customer, the Termination Payments shown above are independent covenants and not conditioned on the occurrence of any event or discharged through any failure of Columbia Tech's performance.

- e. Upon any termination of this Agreement, Columbia Tech shall (i) reasonably cooperate and assist in Customer's transition to another contract manufacturer, which shall only consist of, upon request, informing all suppliers and subcontractors used by Columbia Tech in connection with manufacturing Products of such transition and using good faith efforts to cause each such supplier or subcontractor to support Customer's new contract manufacturer, and by cooperating with Customer to effect the transfer to Customer's new contract manufacturer of any Consigned Materials in the possession or under the control of Columbia Tech, (ii) except as required to comply with any surviving obligations, return and deliver to Customer or, at Customer's option, destroy all tangible embodiments of the Customer's Confidential Information or other intellectual property in the possession or under the control of Columbia Tech, and (iii) except as required to comply with its surviving obligations, discontinue in all respects its use of Customer's intellectual property.
- f. Without limiting Section 19.c, Sections 5 (as to amounts accrued as of termination or arising pursuant to any post-termination activities, 6, 7, 8, 9, 10, 11, 12, 13, 14, 15, 16, 17, 18, 19 (as applies to post-termination activities), 20, 22, 24, 27, 29, 31, 32 and 33 shall survive any termination of this Agreement for the period stated therein or, if no period is stated therein, indefinitely. Termination of this Agreement for any reason shall not affect any rights or obligations that may have accrued prior to such termination or limit any rights or remedies that may otherwise be available to a Party at law or in equity.
- g. Upon Termination, Columbia Tech may apply any remaining Deposit to any amount owed by Customer pursuant to this section or due to any other agreement.

- h. Nothing herein shall restrict Columbia Tech's rights to terminate pursuant to Section 14 or place any additional obligation on Columbia Tech in connection with such termination. Customer shall pay all Termination Payments upon any such exercise of the option by Columbia Tech.
20. JIGS, FIXTURES AND TOOLING: Any jigs, fixtures or tooling purchased by Customer shall remain the property of Customer after the termination of this Agreement and shall be delivered by Columbia Tech to Customer or its designee in accordance with Customer's instructions. All such jigs, fixtures and tooling supplied to Columbia Tech shall be properly labeled by Customer.
21. SUPPLIERS AND SUBCONTRACTORS.
- a. Columbia Tech shall not incorporate any materials that have been identified on the Specification or other approved vendor list as originating from a specific manufacturer in connection with the manufacture of Products unless materials have been approved in writing by Customer; provided, however, that: (a) any supplier or subcontractor specifically identified in the Specifications or an AVL shall be deemed approved by Customer as to the applicable activities; (b) nothing shall limit Columbia Tech's ability to obtain materials from distributors, secondary sources, or other suppliers provided that the requirements from item (a) above are met; and (c) if the specification or bill of materials list a generic item, then Columbia Tech may obtain such generic item from any supplier.
 - b. When ordering materials or using subcontractors, Columbia Tech will send its standard purchase order terms or operation under agreements it has established with those suppliers. Except as provided in Section 8, nothing herein shall require Columbia Tech to obtain any terms with any supplier or contractors, provided, however, that no use of subcontractors shall relieve Columbia Tech of any of its obligations under this Agreement.
22. HEADINGS: Section headings are not to be considered a part of this Agreement and are not intended to be a full and accurate description of the contents hereof.
23. CHANGE ORDERS: Any changes to any Order or Specifications, in whole or in part, must be approved in writing by both Parties prior to implementation. Subject to Section 4.b, Columbia Tech shall have the right to change prices in conjunction with any Change Order, and may also change any production schedule as reasonably necessary to implement the change order. Columbia Tech shall continue to work on all active orders until a change order is signed by both Parties.
24. LIMITATION OF LIABILITY: In addition to any other exclusion that may appear in this Agreement, under no circumstances shall either Party be responsible for any incidental damages, consequential damages, indirect damages, lost profits of any nature, loss of reputation, or loss of orders, even if due to breach of contract or breach of warranty; provided, however, that such limitation shall not apply to Customer's indemnification

obligations hereunder, a Party's breach of confidentiality obligations or use of the other Party's intellectual property other than as expressly authorized hereby or any Late Delivery Credit. The maximum remedy that Customer may have, and the maximum liability of Columbia Tech pursuant to this Agreement, whether due to contract, tort or any other theory, shall be an aggregate amount of one million dollars (\$1,000,000.00) (the "Cap"); provided, however, that such limitation shall not apply to Columbia Tech's breach of confidentiality obligations or use of Customer's intellectual property other than as expressly authorized hereby, any Late Delivery Credit or any loss, damage or destruction of Consigned Materials for which Columbia Tech is responsible under Addendum 1. The Cap shall not increase any other liability that is otherwise limited or liquidated in this Agreement and all such liabilities (except as excluded pursuant to the preceding sentence) shall contribute to the aggregate amount of the Cap.

25. PREVIOUS ORDERS. Both Parties agree that the previous Orders shown at Attachment A shall be subject to the terms and conditions recited in this Agreement.
26. EMPLOYEES AND ACCESS TO BUSINESS PREMISES: Columbia Tech and Customer both acknowledge that the other may from time to time have access to the other Company's business premises. Columbia Tech and Customer covenant and warrant that they have adequate general liability insurance and worker's compensation insurance to cover any employee, agent, or representative that is given access to the business premises of the other. Columbia Tech and Customer also covenant they will follow any rules established by the other for access to the business premises. Columbia Tech and Customer also expressly warrant they will pay all employment taxes, benefits taxes and salary of their respective employees, and nothing in this agreement shall confer employment status on any person. Upon request, either Party shall deliver to the other certificates of insurance evidencing their insurance coverage. Except as expressly provided herein, neither Columbia Tech nor Customer is given any license to enter the other's business premises by virtue of this Agreement.
27. CONFIDENTIALITY.
- a. "Confidential Information" means any technical, business or other information that is disclosed by or on behalf of a Party (the "Disclosing Party") to the other Party (the "Receiving Party") under or in connection with this Agreement; provided, however, that (i) the terms of this Agreement shall be deemed the Confidential Information of each Party and (ii) the Specifications, Inventions and Customer Records shall be deemed the Confidential Information of Customer only.
 - b. During the term of this Agreement and for a period of seven (7) years thereafter, the Receiving Party shall not (i) use Confidential Information of the Disclosing Party for any purpose other than exercising the Receiving Party's rights or performing the Receiving Party's obligations under this Agreement or (ii) disclose Confidential Information of the Disclosing Party to any person or entity other than the Receiving Party's employees, agents, suppliers and contractors on a

need-to-know basis and under commercially reasonable obligations of confidentiality; provided, however, that the Receiving Party shall be responsible to the Disclosing Party for any unauthorized use or disclosure of the Disclosing Party's Confidential Information by any such person. The Receiving Party shall use the same degree of care to prevent unauthorized use or disclosure of the Disclosing Party's Confidential Information as it uses to prevent unauthorized use or disclosure of its own confidential information, but no less than a reasonable degree of care. The Receiving Party shall notify the Disclosing Party in writing promptly after becoming aware of any unauthorized use or disclosure of the Disclosing Party's Confidential Information and cooperate with the Disclosing Party as the Disclosing Party may reasonably request in responding to such unauthorized use or disclosure.

- c. Notwithstanding Section 27.b, the Receiving Party's obligations hereunder shall not apply to any Confidential Information of the Disclosing Party that (i) is generally known to the public through no wrongful act or omission of the Receiving Party, (ii) is received by the Receiving Party, without confidentiality restrictions, from a third party without breach of any confidentiality obligations owed by such third party to the Disclosing Party, as evidenced by the Receiving Party's written records, (iii) was in the Receiving Party's possession, without confidentiality restrictions, before it was disclosed to the Receiving Party by the Disclosing Party, as evidenced by the Receiving Party's written records, or (iv) is independently developed by the Receiving Party, without use of any Confidential Information of the Disclosing Party, as evidenced by the Receiving Party's written records; provided, however, that clauses (iii) and (iv) above shall not limit Columbia Tech's obligations with respect to the Specifications, Inventions and Customer Records.
- d. Notwithstanding Section 27.b, the Receiving Party may disclose Confidential Information of the Disclosing Party (i) as required by applicable law, regulation or court order, so long as the Receiving Party provides the Disclosing Party with reasonable advance notice (to the extent possible and permitted by law) and reasonably cooperates with any efforts of the Disclosing Party to challenge, limit or seek a protective order for any such disclosure; (ii) to the extent reasonably necessary in filing or prosecuting patent applications or preparing or submitting regulatory filings; or (iii) in communications with existing or bona fide prospective acquirers, merger partners, lenders, investors, licensees, sublicensees or collaborators, and consultants and advisors of Receiving Party in connection with transactions or bona fide prospective transactions with any of the foregoing, in each case on a need to know basis and under commercially reasonable obligations of confidentiality; provided, however, that the Receiving Party shall be responsible to the Disclosing Party for any unauthorized use or disclosure of the Disclosing Party's Confidential Information by any such person under this clause (iii).

- e. The Parties acknowledge and agree that any breach or threatened breach of the Parties' obligations under Section 27 may result in irreparable injury to the non-breaching Party. Accordingly, in the event of any such breach or threatened breach, in addition to any other rights or remedies that may be available to the non-breaching Party, the non-breaching Party shall be entitled to seek preliminary and permanent injunctive relief, without the necessity of proving irreparable injury or actual damages and without the necessity of posting a bond.

28. **SUSPENSION:** Columbia Tech may suspend performance under this Agreement, if at any time:

- a. Columbia Tech reasonably believes that the introduction of the Product into commerce shall cause imminent danger; or
- b. Columbia Tech reasonably believes that the Products in their current form are misbranded, adulterated, or otherwise in violation of any law, rule or regulation.

Immediately after becoming aware of any such issue, and prior to initiating any such suspension, Columbia Tech shall provide written notice to Customer describing the issue in appropriate detail. Thereafter, Columbia Tech shall use its best efforts to agree with Customer on any required changes to address the issue and resume performance as soon as possible.

- 29. **SEVERABILITY:** If any provision (or sub-provision) of this Agreement is declared invalid by a court of competent jurisdiction, such provision (or sub-provision) shall be ineffective only to the extent of such invalidity, so that the remainder of the provision and all remaining provisions of this Agreement will continue in full force and effect.
- 30. **FORCE MAJEURE:** Columbia Tech shall not be liable for delays in or non-performance of its obligations under this Agreement as a result of strikes, lockouts, fires, war conditions, accidents, foreign or domestic governmental controls or other actions, embargoes or other causes beyond Columbia Tech's control, including but not limited to, any delay from receiving material from one of Columbia Tech's suppliers beyond Columbia Tech's control.
- 31. **INAPPROPRIATE AND DANGEROUS USES:** Customer hereby represents and warrants that no Products will be used for an inherently dangerous purpose, including, but not limited to, the operation of rail systems, or other motorized devices; demolition or activities involving explosive discharges; or control of artillery, armaments or other instruments of warfare. Customer further represents and warrants that no Products, goods, services or any delivery made from Columbia Tech to Customer shall at any time be used in, installed on or in, become spare parts for, be relied upon for, be used in the construction of, or be used in any way whatsoever in conjunction with aircraft, missiles, or spacecraft (collectively "Aircraft"), including any ground handling tools or equipment,

training aids, test equipment, instructions, manuals, blueprints, engineering tools, or other services and labor relating in any way to Aircraft or the aerospace industry.

32. ENTIRE AGREEMENT: This Agreement and its Attachments and Addendums and the Quality Agreement (if any) constitute the entire agreement and understanding between the Parties with respect to the subject matter hereof and supersedes all previous negotiations, agreements and representations between the Parties, written or oral, all of which shall be deemed to be merged into this Agreement.
33. DISCLAIMERS: Columbia Tech makes no representations or warranties (i) regarding the intellectual property rights of Customer in any invention, discovery, design or Products, (ii) regarding any actual or potential infringement of the Products on any intellectual property or other rights of any person or entity, and (iii) regarding the prior development or current existence of any invention, discovery, design or product similar to the Products. Columbia Tech expressly disclaims all liability and responsibility regarding (A) safety testing or warnings necessary or desirable in connection with any of the Products and (B) compliance with any federal, state or local laws with respect to the Products. Columbia Tech shall have no liability or responsibility to conduct any investigation or inquiry with respect to the foregoing; provided, however, Columbia Tech shall disclose to Customer all infringements upon the rights or patents of others and all violations of federal, state or local laws for which Columbia Tech has actual knowledge. Notwithstanding the foregoing, nothing in this Section 33 shall be deemed to limit any warranties or obligations of Columbia Tech that are expressly set forth in any other provision of this Agreement.
34. TRAINING: Upon Columbia Tech's request, Customer shall provide training regarding any process, know how, skill, or other technique known or possessed by Customer and reasonably needed for Product manufacture to any persons designated by Columbia Tech.
35. NON-RECURRING ENGINEERING; RECURRING COSTS, CO-BUILD: any quotation, time and materials costs for non-recurring engineering, the participation in co-builds or recurring charges related to the manufacture of Products; provided, however, that Customer shall not be responsible for any such amounts unless specifically included in the applicable Quotation.

THE REMAINDER OF THIS PAGE IS INTENTIONALLY BLANK.

Signed Under Seal, effective as of the Effective Date.

Columbia Electrical Contractors, Inc.

/s/ Jason Bourque

Jason Bourque
Sr. Vice President of Finance and
Administration

Akoya Biosciences, Inc.

/s/ Joseph Driscoll

Name: Joseph Driscoll

Title: Chief Financial Officer

ATTACHMENT A- ADDITIONAL ORDERS SUBJECT TO THIS AGREEMENT NONE

ATTACHMENT B – This Attachment is intentionally blank

ATTACHMENT C – STORAGE FEES FOR STORED PRODUCT DISCUSSION ON THIS
ATTACHMENT IS STILL OPEN.

ADDENDUM 1: CUSTOMER CONSIGNED MATERIALS AGREEMENT

This Customer Consigned Materials Agreement, effective as of the date last signed below, is by and between Columbia Electrical Contractors d/b/a **Columbia Tech** (“Columbia Tech”) with a place of business at 27 Otis Street, Westborough, MA 01581, and Akoya Biosciences, Inc., a Delaware Corporation with a place of business at 68 Elm Street, Hopkinton, MA 01748, in connection with that certain Manufacturing Agreement between the Parties (the “Manufacturing Agreement” or “MA”). Capitalized terms used but not defined herein shall have the meanings given to them in the Manufacturing Agreement.

The parties agree:

1. **Consigned Materials.** To the extent contemplated by any Quotation, or otherwise with the consent of Columbia Tech, Customer may from time to time provide Columbia Tech certain materials that Columbia Tech will then use when manufacturing goods for Customer (the “Consigned Materials”). Columbia Tech, at its election, may: issue a purchase order indicating such order is an SMI Order to receive the Consigned Materials (an “SMI Order”). Should Columbia Tech agree to receive such Consigned Materials, and Customer does not provide the Materials on the date stated by Columbia Tech on any such document, Columbia Tech may charge Customer for all such costs associated with the delay in receiving the Consigned Materials, including but not limited to any amount due to Customer for Late Delivery or costs associated with obtaining similar materials.

2. **Storage Area.** Columbia Tech will store Consigned Materials at a facility owned or leased by Columbia Tech (a “Facility”). Columbia Tech will make available a physically segregated area within a Facility (the “Storage Area”). Columbia Tech shall store, handle and transport the Consigned Materials in accordance with of the same care used for its own materials. Columbia Tech shall not use the Consigned Materials with customers other than Customer.

3. **Title to Consigned Materials.** Title to Consigned Materials will remain with Customer at all times. Columbia Tech shall keep Customer’s title to the Consigned Materials free and clear of all liens and encumbrances. If any Consigned Materials are lost, damaged or destroyed due to any breach, negligence or other factors under the control of Columbia Tech, then Columbia Tech shall be responsible for the replacement cost of the affected Consigned Materials. Columbia Tech shall be responsible for insuring the Consigned Materials against all risk and all perils while they are in the possession or under the control of Columbia Tech. Columbia Tech will provide Customer with certificate(s) of insurance from insurers reasonably satisfactory to Customer. Such certificate(s) will be endorsed to contain a provision requiring the insurers to provide Customer with thirty (30) days written notice of any cancellation or adverse material change in such insurance. Within thirty (30) days from the execution of this document, and throughout the term hereof at appropriate intervals, Columbia Tech will deliver satisfactory certificate(s) evidencing such insurance coverage to Customer.

Notwithstanding anything to the contrary, the insurance provisions set forth in this Section 3 will survive any termination of this Addendum.

4. **Inspection of Storage.** Upon reasonable notice, Customer's personnel or agent will have the right to inspect the inventory of Consigned Materials located in the Storage Area and perform a full or partial physical count of all Consigned Materials. The inspection will be scheduled at times mutually agreed upon by the parties so as to minimize any impact to Columbia Tech's operations. Columbia Tech will provide an authorized representative to review such physical inventory or cycle count and initial the resulting physical inventory or cycle count report.

5. **Credit for Consigned Materials.** Columbia Tech will continue to invoice Customer for any Products manufactured by Columbia Tech at the price set forth on the applicable purchase order for the Products related to the invoice. Each month, Columbia Tech shall supply a report of any Consigned Materials used in the Build of the Products. Columbia Tech will place on the account of customer a credit for the Consigned Materials equal to the unit value of the materials shown on the SMI Order; ("Credit Value"). The Credit Value shall not include any material handling fees, material markup, profit, SG&A or other charges Columbia Tech may impose for the build of Products or the services shown hereunder.

6. **No Purchase Obligation.** Columbia Tech shall not have any purchase obligation for the Consigned materials under any circumstances. Any listing of credits, prices, or other units of value on quotations, invoices, purchase orders, SMI orders or any similar documents is for recordkeeping purposes only. Under no circumstances shall Columbia Tech be responsible for payment to Customer for any Consigned Materials (excluding credit or replacement for loss, damage or destruction for which Columbia Tech is responsible hereunder, subject to paragraph 16 below) or have title and the risk of loss to the Consigned Materials.

7. **Storage and Management Fees.** Columbia Tech may charge and invoice Customer fees for the management and storage of Consigned Materials in amounts shown in Columbia Tech quotations or SMI Orders.

8. **Term And Termination** This Customer Consigned Materials Agreement shall become effective once signed by both parties and shall remain effective until terminated as provided for herein. This Customer Consigned Materials Agreement shall terminate upon the termination of the MA (as defined below). Should Customer have any credit remaining in its account based on Section 5, and such credit cannot be applied to any future invoice (whether pursuant to this Agreement, the MA or otherwise), Columbia Tech shall refund such credit amount.

9. **No Warranties.** Columbia Tech offers no warranties whatsoever on the Consigned Materials. In the event Columbia Tech has made any warranty on materials to Customer for the products containing Consigned Materials, then such warranty shall not apply to the Consigned Materials. Customer agrees to receive all Consigned Materials "AS IS" and "WHERE IS".

10. Order of Precedence. This Agreement deals with the procedures for Consigned Materials. It shall not replace any agreement regarding the purchase of products except as it relates to the subject matter herein. For all such matters, this Agreement is completed integrated, and except for such documents contemplated and expressly shown to be issued by the parties, contains the full and complete agreement of the parties. In the event of any conflict between the terms of this document and any other agreement, then this Agreement shall control. The parties do acknowledge the Manufacturing Agreement.

11. **[Reserved]**

12. Return of Consigned Materials. Once materials have been consigned, Columbia Tech may keep such materials until the earlier of: (a) Columbia Tech uses the materials on orders from Customer; (b) Columbia Tech sends Customer 90 days' prior written notice that it is returning Consigned Materials, provided that Columbia Tech must hold any such Consigned Materials for at least ninety (90) days; (c) Customer sends Columbia Tech a notice to return Consigned Materials; or (d) termination of this Agreement (a "Return Event"). Customer agrees to receive all Consigned Materials identified in a Return Event. In the event Customer does not accept any such return delivery or expressly states in writing that it will refuse to accept delivery of such return, Columbia Tech may at its election, at the end of a 90-day period following written notice to Customer, sell or dispose of the applicable Consigned Materials and charge the Customer for all such reasonable costs and expenses related to such sale or disposition net of sales proceeds. All deliveries of Consigned Materials to Columbia Tech shall be DDP. Columbia Tech's premises in Westborough, Massachusetts or its suburban area. All deliveries from Columbia Tech to Customer shall be EXW. Columbia Tech's premises. For avoidance of doubt, Customer shall bear all costs of delivery, all costs for clearing an item for export and all risk of loss to items in transit.

13. Payment/Security Interest/Enforcement Expenses. Columbia Tech may invoice Customer for any amount due under this agreement at any time after such charge is authorized. Customer agrees to pay all such charges as shown in the MA and subject to any security interests, collection expenses or other items shown in the MA.

14. Delivery. All deliveries of Consigned Materials to Columbia Tech shall be DDP Columbia Tech's dock (with Customer acting as the Supplier). All deliveries from Columbia Tech to Customer shall be EXW Columbia Tech's dock (with Columbia Tech acting as the Supplier). For avoidance of doubt, Customer shall bear all costs of delivery, all costs for clearing an item for export and all risk of loss to items in transit.

15. Other Terms. All terms related to Waiver, Relationship of the Parties, Assignment, Notices, Governing Law, Jurisdiction, Severability and Force Majeure shall be as listed in the MA and such terms shall be considered integrated herein.

QUANTERIX CORPORATION

INSIDER TRADING POLICY

TABLE OF CONTENTS

	<u>Page</u>
I. The Need for an Insider Trading Policy	3
II. What is Material Non-Public Information?	3
III. The Consequences of Insider Trading	4
IV. Our Policy	5
General Prohibition on Trading.	5
Transactions by Family Members, Others in Your Household and Entities You Control.	5
Other Companies' Non-public Information.	1
Personal or Independent Reasons Are Not Exceptions.	1
Policy Administrator	1
When Information Becomes Public	1
Prohibited Trading Periods	1
Exceptions for Certain Transactions	9
Pre-Clearance of All Acquisitions, Sales and Other Transfers by Certain Company Personnel	10
V. Individual Responsibility	10
VI. Additional Prohibited Transactions	11
VII. Post-Termination Transactions	11
VIII. Company Assistance	11
IX. Certifications	11

Quanterix Corporation (the “Company”) has adopted the following policy regarding trading by Company personnel in the Company’s securities (the “Insider Trading Policy,” or this “Policy”). This Policy applies to *all* Company personnel, including directors, officers, employees and consultants of the Company and its subsidiaries. This Policy also applies to certain family members, other members of a person’s household and entities controlled by Company personnel, as described in Section IV below.

I. The Need for an Insider Trading Policy

This Policy has been developed:

- to educate all Company personnel as to the federal securities laws and the rules of the Securities and Exchange Commission (the “SEC”) on insider trading in public company securities;
- to set forth requirements that apply to Company personnel and other persons covered by this Policy who seek to trade in the Company’s securities;
- to protect the Company and its personnel from legal liability; and
- to preserve the reputation of the Company and its personnel for integrity and ethical conduct.

Because the Company is a public company, transactions in the Company’s securities are subject to the federal securities laws and regulations adopted by the SEC. These laws and regulations make it illegal for an individual to buy or sell securities of the Company while aware of *material non-public information*. The SEC takes insider trading very seriously and devotes significant resources to uncovering the activity and to prosecuting offenders. Liability may extend not only to the individuals who trade while in possession of material non-public information but also to their “tippers,” people who leak material non-public information to individuals who then trade based on that information. The Company and “controlling persons” of the Company may also be liable for violations by Company employees.

II. What is Material Non-Public Information?

Definition.

Material non-public information is any information (positive or negative) that:

- is not generally known to the public, and
- which, if publicly known, would likely affect either the market price of the Company’s securities or a person’s decision to buy, sell or hold the Company’s securities.

Examples. Common examples of information that will frequently be regarded as material include, but are not limited to:

- quarterly or annual earnings results;
- projections of future financial results;
- the gain or loss of a significant customer or supplier;
- news regarding timing of development or launch of new products;
- news of regulatory approvals or failure to ascertain required regulatory approvals;
- news of a pending or proposed merger, acquisition or tender offer;
- news of a pending or proposed acquisition or disposition of a significant asset;
- news of a pending or proposed joint venture;
- a company restructuring;
- significant transactions with officers, directors or greater than 5% shareholders;
- financing transactions;
- the declaration of a stock split or the offering of additional securities;
- establishment of a stock repurchase program;
- changes in management;
- changes in auditors, or a notification that the auditor's reports may no longer be relied upon;
- pending or threatened significant litigation, or the resolution of such litigation;
- impending bankruptcy or financial liquidity problems;
- the gain or loss of a significant contract, license, registration or collaboration; or
- the entry into, amendment or termination of a material contract.

Twenty-Twenty Hindsight. The foregoing list is illustrative and is not exhaustive of the type of information that may be deemed to be material. In determining whether information is material, the SEC and other regulators will view the information after-the-fact with the benefit of hindsight. As a result, in determining whether any information is material, we will and you should carefully consider whether regulators and others might view the information as being material in hindsight, with the benefit of all relevant information that later becomes available. For example, if there is a significant change in the Company's stock price following release of certain information, that information will likely be determined to have been material when viewed with the benefit of hindsight.

In addition to addressing the relevant statutes and regulations in this area, we are adopting this Policy to avoid even the appearance of improper conduct on the part of anyone employed by or associated with the Company and certain related persons, not just members of senior management.

III. The Consequences of Insider Trading

The consequences of insider trading violations can be severe:

For individuals who trade while in possession of material non-public information (or tip information to others):

- a civil penalty of up to three times the profit gained or loss avoided;

- a criminal fine (no matter how small the profit) of up to \$5 million; and
- a jail term of up to 20 years.

These penalties can apply even if the individual is not a member of the Board of Directors or an officer of the Company. **Moreover, if an employee violates this Policy, he or she may also be subject to Company-imposed sanctions, including termination for cause.**

For a Company (as well as possibly any supervisory person) that fails to take appropriate steps to prevent illegal trading:

- a civil penalty of the greater of \$1 million or three times the profit gained or loss avoided as a result of the employee's violation; and
- a criminal penalty of up to \$25 million.

Any of the above consequences, including an SEC investigation that does not result in prosecution, can tarnish the Company's or an individual's reputation and irreparably damage a career.

IV. Our Policy

General Prohibition on Trading. Company personnel and Related Persons (as defined below in this Section IV) may not buy or sell securities of the Company while in possession of material non-public information or engage in any other action to take advantage of, or pass on to others, that information, subject to the specific exceptions noted below in this Section IV under the caption "Exceptions for Certain Transactions."

Transactions by Family Members, Others in Your Household and Entities You Control. The restrictions in this Policy also apply to (1) immediate family members who reside with you, (2) others living in your household (whether or not related to you), (3) family members who do not live in your household but whose transactions in the Company's securities are directed by you or are subject to your influence or control (e.g., parents or children who consult with you before they trade in the Company's securities) and (4) any entities that you influence or control, including any corporations, limited liability companies, partnerships or trusts (each person or entity identified in clauses (1) – (4), a "Related Person"). SEC regulations specifically provide that any material non-public information about the Company communicated to any spouse, parent, child or sibling is considered to have been communicated under a duty of trust or confidence; and that any trading in the Company's securities by such family members while they are aware of such information may, therefore, violate insider trading laws and regulations. Company personnel are expected to be responsible for the compliance of all Related Persons with this Policy. This means that, to the extent such Related Persons of Company personnel intend to trade in the Company's securities, the Related Persons need to comply with the black-out periods and all other restrictions in this Policy. Furthermore, you should not participate in any investment club (i.e., groups of people who pool their money to make investments) that may invest in the Company's securities.

Other Companies' Non-public Information. This Policy also applies with equal force to information relating to any other company, including our collaborators, vendors and customers, obtained by Company personnel during the course of their service to or employment by the Company. Specifically, no Company personnel who, in the course of work on behalf of the Company, learns of material non-public information about a company with which the Company does business may trade in the other company's securities until the information becomes public or is no longer material.

Personal or Independent Reasons Are Not Exceptions. Transactions in the Company's securities that may be necessary or justifiable for independent reasons (such as the need to raise money for an emergency expenditure) are no exception. Even the appearance of an improper transaction must be avoided to preserve our reputation for adhering to the highest standards of conduct.

Policy Administrator. This Policy shall be administered by the "Policy Administrator," who shall be the Company's General Counsel, and if such person is not available, then the Company's Vice President, Senior Legal Counsel or Chief Financial Officer, shall serve as the alternate Policy Administrator. The Policy Administrator may, however, change from time to time; to confirm the name of the then-current Policy Administrator, please check the Company's internal HR website.

When Information Becomes Public. This Policy applies to material *non-public* information about the Company, which means that trading is permitted once the information becomes known to the public (unless some other Company policy or legal obligation restricts trading at that time). Because the Company's shareholders and the investing public should be afforded time to receive and absorb information, as a general rule you should not engage in any transactions until the beginning of the second business day after material information has been released. Thus, if an announcement is made on a Monday, Wednesday generally would be the first day on which you may trade. If an announcement is made a Friday, Tuesday generally would be the first day on which you may trade. However, if the information released is complex, such as a major financing or other significant transaction, it may be necessary to allow additional time for the information to be absorbed by the investing public. In such circumstances, you will be notified by the Policy Administrator regarding a suitable waiting period before trading. In addition, we have established specified black-out periods, as described below.

Prohibited Trading Periods. While it is never permissible to trade based on material non-public information, we are implementing the following procedures to help prevent inadvertent violations of this Policy and avoid even the appearance of an improper transaction (which could result, for example, where Company personnel engage in a trade while unaware of a pending major development).

(1) Company Wide Black-Out Periods Applicable to All Company Personnel. All Company personnel and Related Persons are prohibited from trading in any of the Company's securities during the following periods:

- from the time each such individual becomes aware of the material information (the black-out start times often vary), until the beginning of the second business day after the day the Company has made a public announcement of material information, including earnings releases, unless the information released is complex, in which case it may be necessary to extend this period and the Policy Administrator will notify you of any such extension of the black-out period; and
- during other specified periods when significant developments or announcements are anticipated, as notified by the Policy Administrator.

You will be notified by e-mail when you may not trade in the Company's securities during periods when significant developments or announcements are anticipated, in which event you will also be notified when trading restrictions are lifted. ***Of course, even during periods when trading is permitted, no one, including persons or entities who do not fall within the definition of Related Persons, should trade in the Company's securities if he or she possesses material non-public information.***

(2) Additional Black-Out Periods Applicable to the Board of Directors, Senior Management, Financial Team Members and Designated Employees. In addition to being subject to the trading procedures applicable to all Company personnel (above), members of the Company's Board of Directors, Senior Management, Financial Team Members, Designated Employees (each as defined below) and Related Persons of such individuals are also subject to additional trading procedures and restrictions during the following periods:

- the periods from 15 days prior to the close of each fiscal quarter until the beginning of the second business day after the release of the Company's financial results for each quarter and, in the case of the fourth quarter, financial results for the year end; and
- any other periods as determined by the Company.

The following members of management constitute the "Senior Management" of the Company: all Executive (Section 16) Officers and Executive Team members.

The following individuals constitute the "Financial Team Members" of the Company: all members of the Company's financial team.

The following individuals constitute other "Designated Employees" of the Company: all members of the Company's sales and marketing team, and certain additional members of Company personnel, as listed on Exhibit A hereto, which list shall be amended from time to time to reflect the then-current group of such individuals.

The Policy Administrator may, from time to time, amend the list of and/or designate other employees as Senior Management, Financial Team Members or Designated Employees, in which case the Policy Administrator shall notify the affected individuals.

Exceptions for Certain Transactions.

(1) **Gifts.** *Bona fide* gifts are not transactions that are subject to this Policy, unless the person making the gift (the donor) has reason to believe that the recipient of the gift intends to sell the Company's securities while the donor is in possession of material non-public information.

(2) **Mutual Funds.** Transactions in mutual funds that are invested in the Company's securities are not transactions subject to this Policy.

(3) **Transactions Involving Company Equity Plans.** Except as otherwise noted below, this Policy does not apply to the following transactions:

- ***Stock Option Exercises.*** This Policy does not apply to the exercise of an employee stock option acquired pursuant to the Company's equity plans, or to the exercise of a tax withholding right pursuant to which a person has elected to have the Company withhold shares subject to an option to satisfy tax withholding requirements. This Policy does apply, however, to any sale of stock as part of a broker-assisted cashless exercise of an option, or any other market sale of stock for the purpose of generating the cash needed to pay the exercise price and or taxes upon the exercise of an option.
- ***Restricted Stock Awards and Restricted Stock Unit Awards.*** This Policy does not apply to the vesting of restricted stock or restricted stock units, or the exercise of a tax withholding right pursuant to which a person elects to have the Company withhold shares of stock to satisfy tax withholding requirements upon the vesting of any restricted stock or restricted stock unit. This Policy does apply, however, to any market sale of restricted stock or shares received upon vesting of restricted stock units.
- ***Employee Stock Purchase Plan.*** This Policy will not apply to purchases of the Company's securities under an employee stock purchase plan. This Policy does apply, however, to subsequent sales or other transfers of such securities.
- ***Other Transactions with the Company.*** Any other purchase of the Company's securities from the Company or sales of the Company's securities to the Company are not subject to this Policy.

(4) **Rule 10b5-1 Trading Plans.** Notwithstanding the restrictions and prohibitions on trading in the Company's securities set forth in this Policy, persons subject to this Policy are permitted to effect transactions in the Company's securities pursuant to approved trading plans established under Rule 10b5-1 of the Securities Exchange Act of 1934, as amended ("**Trading Plans**"), which may include transactions during the prohibited periods discussed above. Rule 10b5-1 requires that these transactions be made pursuant to a plan that was established while the person was not in possession of material non-public information, and the SEC requires that these plans not be entered into during any applicable Company-imposed black-out period. In order to comply with this Policy, the Company must pre-approve any such Trading Plan prior to its effectiveness. After a Trading Plan is approved, you must wait for a cooling-off period before the first trade is made under the Trading Plan, the length of which will be determined by the

Policy Administrator. Once the Trading Plan is adopted, you must not exercise any influence over the amount of securities to be traded, the price at which they are to be traded or the dates of the trades. The Trading Plan must either specify the amount, pricing and timing of transactions in advance or delegate discretion on these matters to an independent third party. Any modification of a Trading Plan is the equivalent of entering into a new Trading Plan and cancelling the old Trading Plan. Company personnel seeking to establish, modify or cancel a Trading Plan should contact the Policy Administrator.

Pre-Clearance of All Acquisitions, Sales and Other Transfers by Certain Company Personnel. In order to ensure compliance with this Policy and with any Section 16 reporting requirements, all transactions in the Company's securities (including acquisitions, sales, gifts and other transfers, whether or not for value), including the execution of Trading Plans (as defined below), by members of the Company's Board of Directors, Senior Management, Financial Team Members, Designated Employees and Related Persons, must be pre-cleared by the Policy Administrator. If you are a member of one of the groups listed above and you contemplate a transaction in the Company's securities, you must contact the Policy Administrator or other designated individual prior to executing the transaction. The Policy Administrator will use his or her reasonable best efforts to provide approval or disapproval within two business days. You must wait until receiving pre-clearance to execute the transaction. Neither the Company nor the Policy Administrator shall be liable for any delays that may occur due to the pre-clearance process. If the transaction is pre-cleared by the Policy Administrator, it must be executed by the end of the second business day after receipt of pre-clearance. Notwithstanding receipt of pre-clearance of a transaction, if you become aware of material non-public information about the Company after receiving the pre-clearance but prior to the execution of the transaction, you may not execute the transaction. The responsibility for determining whether you are in possession of material non-public information rests with you, as discussed below in Section V. If you are a Section 16 reporting person, promptly following execution of the transaction, but in no event later than the end of the first business day after the execution of the transaction, you must notify the Policy Administrator and provide details regarding the transaction sufficient to complete the required Section 16 filing.

Employees of the Company who are not Directors, members of Senior Management, Financial Team Members or Designated Employees may, but are not required to, pre-clear transactions in the Company's securities in the same manner as set forth above. Such employees are not required to notify the Policy Administrator following execution of the transaction.

Please note that pre-clearance does not provide Company personnel with immunity from investigation or suit, for which it is the responsibility of the individual to comply with the federal securities regulations.

V. Individual Responsibility

Persons subject to this Policy have ethical and legal obligations to maintain the confidentiality of information about the Company and to not engage in transactions in the Company's securities while in possession of material non-public information. Each individual is

responsible for making sure that he or she complies with this Policy, and that any Related Person, whose transactions are subject to this Policy, also comply with this Policy. In all cases, the responsibility for determining whether an individual is in possession of material non-public information rests with that individual, and any action on the part of the Company, the Policy Administrator or any other employee or director pursuant to this Policy (or otherwise) does not in any way constitute legal advice or insulate an individual from liability under applicable securities laws. You may be subject to legal penalties and disciplinary action by law enforcement officials and/or the Company for any conduct prohibited by this Policy or applicable securities laws, as described in Section III above.

Tippling Information to Others. Company personnel must not disclose non-public information about the Company to others outside the Company who do not have an obligation to maintain the confidentiality of such information. If the outsider trades on such information, penalties for insider trading may apply in these situations whether or not you derive any monetary benefit from the other person's trading activities. Material non-public information is often inadvertently disclosed or overheard in casual, social conversations. Please take care to avoid such disclosures.

Prevention of Insider Trading by Others. If you become aware of a potential insider trading violation, you must immediately advise our Policy Administrator and/or report the matter using the Company's anonymous whistleblower reporting procedures. You should also take steps, where appropriate, to prevent persons under your supervision and/or control from using material non-public information for trading purposes. Moreover, Company-imposed sanctions, including termination for cause, could result if an employee fails to comply with this Policy.

Confidentiality. Serious problems could be caused for the Company by the unauthorized disclosure of internal information about the Company, whether or not for the purpose of facilitating improper trading in the Company's securities. Company personnel should not discuss internal Company matters or developments (whether or not you think such information is material) with anyone outside of the Company (including, but not limited to, family, friends, business associates, investors and expert consulting firms), except as required in the performance of regular corporate duties. This prohibition applies specifically (but not exclusively) to inquiries about the Company that may be made by the financial press, investment analysts or others in the financial community and also includes posting material non-public information on any social media outlets such as Facebook, Twitter, etc. It is important that all such communications on behalf of the Company be made only through an authorized officer under carefully controlled circumstances. Unless you are expressly authorized to the contrary, if you receive any inquiries of this nature, you should decline comment and refer the inquirer to the Company's Chief Financial Officer or General Counsel.

VI. Additional Prohibited Transactions

Because we believe it is generally improper and inappropriate for Company personnel to engage in short-term or speculative transactions involving the Company's securities, it is our policy that Company personnel and Related Persons not engage in any of the following activities, except in each case in limited circumstances with prior approval of the Policy Administrator:

- trading in the Company's securities on a short-term basis. Any shares of Company common stock purchased in the open market must be held for a minimum of six months and ideally longer;
- short sales of the Company's securities;
- use of the Company's securities to secure a margin or other loan;
- transactions in straddles, collars or other similar risk reduction or hedging devices; and
- transactions in publicly-traded options relating to the Company's securities (i.e., options that are not granted by the Company).

VII. Post-Termination Transactions

This Policy will no longer apply after termination of service to the Company. However, if an individual is in possession of material non-public information when his or her service terminates, that individual may not trade in the Company's securities until that information has become public or is no longer material, and it would be prudent for the individual, if he or she is subject to a black-out period upon termination of service, to refrain from trading until those restrictions no longer apply to Company personnel.

VIII. Company Assistance

Any person who has any questions about specific transactions or this Policy in general may obtain additional guidance from the Policy Administrator. Remember, however, the ultimate responsibility for adhering to this Policy and avoiding improper transactions rests with you. In this regard, please use your best judgment when considering a transaction in the Company's securities.

IX. Certifications

As a condition to employment, all employees will be required to certify their understanding of and intent to comply with this Policy. Members of the Board of Directors, Senior Management and other personnel may be required to certify compliance on an annual basis.

Exhibit A

Designated Employees

Legal Department

IT Department

Executive Team Administrative Assistants

Certification Under Insider Trading Policy

The undersigned hereby certifies that he/she has read and understands, and agrees to comply with, the Company's Insider Trading Policy, a copy of which was distributed with this Certification.

Date: _____ Signature _____

Name: _____
(Please Print)

Title: _____

SUBSIDIARIES

Company Name	Jurisdiction of Incorporation
Akoya Biosciences, Inc.	Delaware
Akoya Biosciences UK Ltd.	United Kingdom
Aushon Biosystems, Inc.	Delaware
Emission Inc.	Delaware
Quanterix Bio-tech (Shanghai) Co., Ltd	China
Quanterix Holdings (Hong Kong) Limited	Hong Kong
Quanterix Netherlands B.V.	The Netherlands
Quanterix Security Corporation	Massachusetts
UmanDiagnostics AB	Sweden



KPMG LLP
Two Financial Center
60 South Street
Boston, MA 02111

Consent of Independent Registered Public Accounting Firm

We consent to the incorporation by reference in the registration statements (Nos. 333-223771, 333-231373, 333-240420, 333-256032, 333-264836, 333-270295, 333-277643, 333-285924, 333-288648, 333-292362, 333-292759, 333-284932) on Form S-8 of our report dated March 2, 2026, with respect to the consolidated financial statements of Quanterix Corporation and subsidiaries and the effectiveness of internal control over financial reporting.

/s/ KPMG LLP

Boston, Massachusetts
March 2, 2026

CONSENT OF INDEPENDENT REGISTERED PUBLIC ACCOUNTING FIRM

We consent to the incorporation by reference in the following Registration Statements:

- (1) Registration Statement (Form S-8 No. 333-223771) pertaining to the 2007 Stock Option and Grant Plan, as amended, the 2017 Employee Stock Purchase Plan, and 2017 Employee, Director, and Consultant Equity Incentive Plan,
 - (2) Registration Statement (Form S-8 No. 333-231373) pertaining to the 2017 Employee Stock Purchase Plan, and 2017 Employee, Director and Consultant Equity Incentive Plan,
 - (3) Registration Statement (Form S-8 No. 333-240420) pertaining to the 2017 Employee Stock Purchase Plan, and 2017 Employee, Director and Consultant Equity Incentive Plan,
 - (4) Registration Statement (Form S-8 No. 333-256032) pertaining to the 2017 Employee Stock Purchase Plan, and 2017 Employee, Director and Consultant Equity Incentive Plan,
 - (5) Registration Statement (Form S-8 No. 333-264836) pertaining to the 2017 Employee Stock Purchase Plan, and 2017 Employee, Director and Consultant Equity Incentive Plan,
 - (6) Registration Statement (Form S-8 No. 333-270295) pertaining to the 2017 Employee Stock Purchase Plan, and 2017 Employee, Director and Consultant Equity Incentive Plan,
 - (7) Registration Statement (Form S-8 No. 333-277643) pertaining to the 2017 Employee Stock Purchase Plan, and 2017 Employee, Director and Consultant Equity Incentive Plan,
 - (8) Registration Statement (Form S-8 No. 333-285924) pertaining to the 2017 Employee Stock Purchase Plan, and 2017 Employee, Director and Consultant Equity Incentive Plan,
 - (9) Registration Statement (Form S-8 No. 333-288648) pertaining to the Akoya Biosciences, Inc. 2021 Equity Incentive Plan, and Quanterix Corporation Restricted Stock Unit Inducement Awards,
 - (10) Post-effective Amendment No. 3 on Form S-8 to Form S-4 Registration Statement (Form S-8 POS No. 333-284932) pertaining to the Akoya Biosciences, Inc. 2021 Equity Incentive Plan, and Akoya Biosciences, Inc. 2015 Equity Incentive Plan,
 - (11) Registration Statement (Form S-8 No. 333-292362) pertaining to the 2025 Inducement Plan, and
 - (12) Registration Statement (Form S-8 No. 333-292759) pertaining to the Amended and Restated 2025 Inducement Plan;
- of our report dated March 17, 2025, with respect to the consolidated financial statements of Quanterix Corporation included in this Annual Report (Form 10-K) of Quanterix Corporation for the year ended December 31, 2024.

/s/ Ernst & Young LLP

Boston, Massachusetts
March 2, 2026

CERTIFICATIONS UNDER SECTION 302

I, Everett Cunningham, certify that:

1. I have reviewed this annual report on Form 10-K of Quanterix Corporation;
2. Based on my knowledge, this report does not contain any untrue statement of a material fact or omit to state a material fact necessary to make the statements made, in light of the circumstances under which such statements were made, not misleading with respect to the period covered by this report;
3. Based on my knowledge, the financial statements, and other financial information included in this report, fairly present in all material respects the financial condition, results of operations and cash flows of the registrant as of, and for, the periods presented in this report;
4. The registrant's other certifying officer(s) and I are responsible for establishing and maintaining disclosure controls and procedures (as defined in Exchange Act Rules 13a-15(e) and 15d-15(e)) and internal control over financial reporting (as defined in Exchange Act Rules 13a-15(f) and 15d-15(f)) for the registrant and have:
 - a) designed such disclosure controls and procedures, or caused such disclosure controls and procedures to be designed under our supervision, to ensure that material information relating to the registrant, including its consolidated subsidiaries, is made known to us by others within those entities, particularly during the period in which this report is being prepared;
 - b) designed such internal control over financial reporting, or caused such internal control over financial reporting to be designed under our supervision, to provide reasonable assurance regarding the reliability of financial reporting and the preparation of financial statements for external purposes in accordance with generally accepted accounting principles;
 - c) evaluated the effectiveness of the registrant's disclosure controls and procedures and presented in this report our conclusions about the effectiveness of the disclosure controls and procedures, as of the end of the period covered by this report based on such evaluation; and
 - d) disclosed in this report any change in the registrant's internal control over financial reporting that occurred during the registrant's most recent fiscal quarter (the registrant's fourth fiscal quarter in the case of an annual report) that has materially affected, or is reasonably likely to materially affect, the registrant's internal control over financial reporting; and
5. The registrant's other certifying officer(s) and I have disclosed, based on our most recent evaluation of internal control over financial reporting, to the registrant's auditors and the audit committee of the registrant's board of directors (or persons performing the equivalent functions):
 - a) all significant deficiencies and material weaknesses in the design or operation of internal control over financial reporting which are reasonably likely to adversely affect the registrant's ability to record, process, summarize and report financial information; and
 - b) any fraud, whether or not material, that involves management or other employees who have a significant role in the registrant's internal control over financial reporting.

Date: March 2, 2026

/s/ EVERETT CUNNINGHAM

Everett Cunningham

President and Chief Executive Officer and Director (Principal Executive Officer)

CERTIFICATIONS UNDER SECTION 302

I, Vandana Sriram, certify that:

1. I have reviewed this annual report on Form 10-K of Quanterix Corporation;
2. Based on my knowledge, this report does not contain any untrue statement of a material fact or omit to state a material fact necessary to make the statements made, in light of the circumstances under which such statements were made, not misleading with respect to the period covered by this report;
3. Based on my knowledge, the financial statements, and other financial information included in this report, fairly present in all material respects the financial condition, results of operations and cash flows of the registrant as of, and for, the periods presented in this report;
4. The registrant's other certifying officer(s) and I are responsible for establishing and maintaining disclosure controls and procedures (as defined in Exchange Act Rules 13a-15(e) and 15d-15(e)) and internal control over financial reporting (as defined in Exchange Act Rules 13a-15(f) and 15d-15(f)) for the registrant and have:
 - a) designed such disclosure controls and procedures, or caused such disclosure controls and procedures to be designed under our supervision, to ensure that material information relating to the registrant, including its consolidated subsidiaries, is made known to us by others within those entities, particularly during the period in which this report is being prepared;
 - b) designed such internal control over financial reporting, or caused such internal control over financial reporting to be designed under our supervision, to provide reasonable assurance regarding the reliability of financial reporting and the preparation of financial statements for external purposes in accordance with generally accepted accounting principles;
 - c) evaluated the effectiveness of the registrant's disclosure controls and procedures and presented in this report our conclusions about the effectiveness of the disclosure controls and procedures, as of the end of the period covered by this report based on such evaluation; and
 - d) disclosed in this report any change in the registrant's internal control over financial reporting that occurred during the registrant's most recent fiscal quarter (the registrant's fourth fiscal quarter in the case of an annual report) that has materially affected, or is reasonably likely to materially affect, the registrant's internal control over financial reporting; and
5. The registrant's other certifying officer(s) and I have disclosed, based on our most recent evaluation of internal control over financial reporting, to the registrant's auditors and the audit committee of the registrant's board of directors (or persons performing the equivalent functions):
 - a) all significant deficiencies and material weaknesses in the design or operation of internal control over financial reporting which are reasonably likely to adversely affect the registrant's ability to record, process, summarize and report financial information; and
 - b) any fraud, whether or not material, that involves management or other employees who have a significant role in the registrant's internal control over financial reporting.

Date: March 2, 2026

/s/ VANDANA SRIRAM

Vandana Sriram

Chief Financial Officer

(principal financial officer and principal accounting officer)

CERTIFICATIONS UNDER SECTION 906

Pursuant to section 906 of the Sarbanes-Oxley Act of 2002 (subsections (a) and (b) of section 1350, chapter 63 of title 18, United States Code), each of the undersigned officers of Quanterix Corporation, a Delaware corporation (the "Company"), does hereby certify, to such officer's knowledge, that:

The Annual Report for the year ended December 31, 2025 (the "Form 10-K") of the Company fully complies with the requirements of Section 13(a) or 15(d) of the Securities Exchange Act of 1934, and the information contained in the Form 10-K fairly presents, in all material respects, the financial condition and results of operations of the Company.

Dated: March 2, 2026

/s/ EVERETT CUNNINGHAM

Everett Cunningham
President and Chief Executive Officer and Director (Principal Executive Officer)

Dated: March 2, 2026

/s/ VANDANA SRIRAM

Vandana Sriram
Chief Financial Officer (Principal Financial Officer and Principal Accounting Officer)