

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

FORM 8-K

CURRENT REPORT
Pursuant to Section 13 or 15(d)
of the Securities Exchange Act of 1934

Date of Report (Date of earliest event reported): November 10, 2025

QUANTERIX CORPORATION

(Exact name of registrant as specified in its charter)

Delaware
(State or other jurisdiction
of incorporation)

900 Middlesex Turnpike
Billerica, MA
(Address of principal executive offices)

001-38319
(Commission File Number)

20-8957988
(IRS Employer
Identification No.)

01821
(Zip Code)

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

Check the appropriate box below if the Form 8-K filing is intended to simultaneously satisfy the filing obligation of the registrant under any of the following provisions:

- Written communications pursuant to Rule 425 under the Securities Act (17 CFR 230.425)
- Soliciting material pursuant to Rule 14a-12 under the Exchange Act (17 CFR 240.14a-12)
- Pre-commencement communications pursuant to Rule 14d-2(b) under the Exchange Act (17 CFR 240.14d-2(b))
- Pre-commencement communications pursuant to Rule 13e-4(c) under the Exchange Act (17 CFR 240.13e-4(c))

Securities registered pursuant to Section 12(b) of the 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

Indicate by check mark whether the registrant is an emerging growth company as defined in Rule 405 of the Securities Act of 1933 (17 CFR §230.405) or Rule 12b-2 of the Securities Exchange Act of 1934 (17 CFR §240.12b-2).

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.

Item 2.02 Results of Operations and Financial Condition.

On November 10, 2025, Quanterix Corporation ("Quanterix") issued a press release announcing its financial results for the third quarter ended September 30, 2025 (the "Earnings Release"). A copy of the Earnings Release is furnished as Exhibit 99.1 and is incorporated herein by reference.

The information in this Item 2.02 of this Form 8-K (including Exhibits 99.1 and 99.2) shall not be deemed "filed" for purposes of Section 18 of the Securities Exchange Act of 1934 (the "Exchange Act") or otherwise subject to the liabilities of that section, nor shall it be deemed incorporated by reference in any filing under the Securities Act of 1933 or the Exchange Act, except as expressly set forth by specific reference in such a filing.

Item 9.01 Financial Statements and Exhibits.

(d) Exhibits

Exhibit No.	Description
99.1	Earnings Release dated November 10, 2025
99.2	Slides from November 10, 2025 Earnings Call
104	Cover Page Interactive Data File (embedded within the inline XBRL document)

SIGNATURES

Pursuant to the requirements of the Securities Exchange Act of 1934, the Registrant has duly caused this report to be signed on its behalf by the undersigned thereunto duly authorized.

QUANTERIX CORPORATION

By: /s/ Vandana Sriram
Vandana Sriram
Chief Financial Officer

Date: November 10, 2025

Quanterix Releases Financial Results for the Third Quarter of 2025

BILLERICA, Mass. – November 10, 2025 - Quanterix Corporation (NASDAQ: QTRX), a company transforming healthcare by accelerating biomarker breakthroughs from discovery to diagnostics, today announced financial results for the third quarter ended September 30, 2025.

“During the third quarter, we delivered on our revenue expectations despite challenging market conditions,” said Masoud Toloue, Chief Executive Officer of Quanterix. “Equally important, we achieved key integration milestones from our Akoya transaction — keeping us firmly on track to realize \$85 million in annualized synergies and reach cash flow break-even in 2026. Together, we have built a stronger, more scalable company with the foundation to drive sustained growth, profitability, and cash flow in the years ahead.”

“Our Alzheimer’s diagnostics business also continues to accelerate. With a positive Medicare pricing recommendation and expanding partnerships across Asia, we are advancing toward a future where blood-based testing for neurodegenerative disease becomes standard clinical practice — and Quanterix is at the center of that transformation.”

Third Quarter Financial Highlights

- Revenue of \$40.2 million, an increase of 12.3% compared to \$35.8 million in the prior year.
- GAAP gross margin of 42.8%, as compared to 56.3% in the prior year. Adjusted gross margin (non-GAAP) of 45.9% as compared to 50.7% in the prior year.
- Adjusted EBITDA (non-GAAP) loss of \$11.9 million, compared to a loss of \$5.5 million in the prior year. The increase is primarily related to higher acquisition and restructuring charges.
- The Company ended the third quarter with \$138.1 million of cash, cash equivalents, marketable securities, and restricted cash. Adjusted cash usage, after accounting for one-time deal and restructuring costs, was \$16.1 million. Adjusted cash usage is expected to decline in Q4 2025 due to improved working capital performance and a full quarter effect of cost synergies. The company continues to expect it will end 2025 with approximately \$120 million in cash, cash equivalents, marketable securities, and restricted cash.

Operational and Business Highlights

- Executed key integration activities in the quarter, capturing \$67 million of cost synergies and savings. Key integration milestones include:
 - *Commercial*: Combined the Quanterix and Akoya sales organizations, under one sales leadership, selling one combined portfolio of products.
 - *Operations*: Consolidated four manufacturing and lab facilities into two main sites in Billerica and Burlington Massachusetts.
 - *Back Office/G&A*: Eliminated nearly all redundant public company costs and initiated a series of IT systems integrations.
- Established new partnerships in the Asia-Pacific region, including All-Eight (Singapore); Union Clinical Laboratory (Taiwan); and NSW Health Pathology (Australia).
- Launched eight new Simoa assays through the first nine months of the year.
- Delivered multiple recent spatial and immuno-oncology presentations, including more than a dozen poster presentations, at the 40th annual Society for Immunotherapy of Cancer (SITC), demonstrating growing adoption of Quanterix’s high-plex spatial proteomics and AI-driven analysis platforms, underscoring their translational impact in uncovering immune resistance mechanisms and predictive biomarkers across multiple cancer indications. The Company also launched a new Metabolism Spike-in Module, designed for the PhenoCycler-Fusion and the PhenoCode™ Discovery IO60 Human Protein Panel.

2025 Full Year Business Outlook

The Company is reiterating its previous guidance for 2025 revenues to be in a range of \$130 to \$135 million, which includes approximately two quarters of performance from Akoya. On a pro forma basis, assuming the two companies were combined for the full year, the Company expects to generate 2025 revenue of between \$165 and \$170 million.

Quanterix expects GAAP and adjusted gross margin (non-GAAP) to be in the range of 45% to 47%.

Finally, the Company anticipates 2025 adjusted cash usage to be approximately \$34 to \$38 million. With the previously announced expense reductions, Quanterix expects to exit 2025 with approximately \$120 million in cash, cash equivalents, marketable securities, and restricted cash and expects to achieve cash flow breakeven performance in 2026.

Conference Call

In conjunction with this announcement, the Company will host a conference call on November 10, 2025, at 4:30 PM E.T. The dial-in number for USA & Canada is Toll-Free (800) 715-9871 or (646) 307-1963 and the conference ID is 5591565.

Interested investors can also listen to the live webcast from the Event Details page in the Investors section of the Quanterix website at <https://ir.quanterix.com>. An archived webcast replay will be available on the Company's website for one year.

About Quanterix

Quanterix is a global leader in ultra-sensitive biomarker detection, enabling breakthroughs in disease research, diagnostics, and drug development. Its proprietary Simoa® technology delivers industry-leading sensitivity, allowing researchers to detect and quantify biomarkers in blood and other fluids at concentrations far below traditional limits. With approximately 6,000 peer-reviewed publications, Quanterix has been a trusted partner to the scientific community for nearly two decades. In 2025, Quanterix acquired Akoya Biosciences, The Spatial Biology Company®, adding multiplexed tissue imaging with single-cell resolution to its portfolio and 1,396 installed instruments. Together, the combined company offers a uniquely integrated platform that connects biology across blood and tissue—advancing precision medicine from discovery to diagnostics. Learn more at www.quanterix.com.

Financial Highlights

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

	Three Months Ended September 30,		Nine Months Ended September 30,	
	2025	2024	2025	2024
Revenues:				
Product revenue	\$ 26,151	\$ 19,694	\$ 63,723	\$ 59,251
Service and other revenue	13,953	13,845	29,806	39,323
Collaboration and license revenue	46	1,872	1,348	2,756
Grant revenue	83	402	165	930
Total revenues	40,233	35,813	95,042	102,260
Costs of goods sold and services:				
Cost of product revenue	15,379	10,554	34,438	25,461
Cost of service and other revenue	7,648	5,106	15,683	15,864
Total costs of goods sold and services	23,027	15,660	50,121	41,325
Gross profit	17,206	20,153	44,921	60,935
Operating expenses:				
Research and development	8,009	8,104	27,126	23,015
Selling, general and administrative	39,062	22,908	102,872	73,027
Other lease costs	286	889	870	2,740
Impairment and restructuring costs	7,174	—	14,844	—
Total operating expenses	54,531	31,901	145,712	98,782
Loss from operations	(37,325)	(11,748)	(100,791)	(37,847)
Other income (expense):				
Interest income	1,448	3,535	7,408	11,165
Change in fair value of contingent liabilities	58	—	3,952	—
Other income (expense), net	(178)	5	(68)	221
Loss before income taxes	(35,997)	(8,208)	(89,499)	(26,461)
Income tax benefit (expense)	2,480	(145)	5,466	(442)
Net loss	\$ (33,517)	\$ (8,353)	\$ (84,033)	\$ (26,903)
Net loss per common share, basic and diluted	\$ (0.73)	\$ (0.22)	\$ (2.04)	\$ (0.70)
Weighted-average common shares outstanding, basic and diluted	46,060	38,449	41,243	38,305

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

	September 30, 2025	December 31, 2024
ASSETS		
Current assets:		
Cash and cash equivalents	\$ 38,298	\$ 56,709
Marketable securities	96,511	232,413
Accounts receivable, net of allowance for expected credit losses	33,029	32,141
Inventory	54,957	32,775
Prepaid expenses and other current assets	10,483	9,556
Total current assets	233,278	363,594
Restricted cash	3,336	2,610
Property and equipment, net	25,765	17,150
Intangible assets, net	135,148	4,031
Goodwill	23,460	—
Operating lease right-of-use assets	18,184	16,339
Financing lease right-of-use assets	917	—
Other non-current assets	4,052	2,809
Total assets	\$ 444,140	\$ 406,533
LIABILITIES AND STOCKHOLDERS' EQUITY		
Current liabilities:		
Accounts payable	\$ 10,633	6,953
Accrued compensation and benefits	12,582	12,620
Accrued expenses and other current liabilities	24,812	8,851
Deferred revenue	20,913	8,827
Operating lease liabilities	7,984	4,756
Financing lease liabilities	428	—
Total current liabilities	77,352	42,007
Deferred revenue, net of current portion	7,341	1,073
Operating lease liabilities, net of current portion	31,212	32,615
Financing lease liabilities, net of current portion	496	—
Non-current portion of contingent liabilities	5,904	—
Other non-current liabilities	7,301	800
Total liabilities	129,606	76,495
Total stockholders' equity	314,534	330,038
Total liabilities and stockholders' equity	\$ 444,140	\$ 406,533

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

	Nine Months Ended September 30,	
	2025	2024
Cash flows from operating activities:		
Net loss	\$ (84,033)	\$ (26,903)
Adjustments to reconcile net loss to net cash used in operating activities:		
Depreciation and amortization expense	9,618	4,740
Credit losses on accounts receivable	284	744
Accretion of marketable securities	(1,851)	(5,317)
Impairment	7,269	—
Operating lease right-of-use asset amortization	2,058	1,371
Stock-based compensation expense	16,303	15,150
Change in fair value of contingent liabilities	(3,952)	—
Deferred taxes	(5,678)	—
Other operating activity	(76)	(389)
Changes in assets and liabilities:		
Accounts receivable	7,901	(6,402)
Inventory	4,204	(6,845)
Prepaid expenses and other current assets	2,198	(168)
Other non-current assets	110	(479)
Accounts payable	(3,537)	588
Accrued compensation and benefits, accrued expenses, and other current liabilities	(5,847)	(3,166)
Deferred revenue	(526)	(826)
Operating lease liabilities	(4,003)	(2,971)
Other non-current liabilities	(1,234)	11
Net cash used in operating activities	(60,792)	(30,862)
Cash flows from investing activities:		
Purchases of marketable securities	(45,658)	(270,972)
Proceeds from sales and maturities of marketable securities	183,389	159,279
Purchases of property and equipment	(2,710)	(2,956)
Acquisitions, net of cash acquired	(93,229)	—
Net cash provided by (used in) investing activities	41,792	(114,649)
Cash flows from financing activities:		
Principal payments on financing leases	(106)	—
Proceeds from common stock issued under stock plans	930	2,999
Payments for employee taxes withheld on stock-based compensation awards	(1,010)	(2,584)
Net cash provided by (used in) financing activities	(186)	415
Net increase (decrease) in cash, cash equivalents, and restricted cash	(19,186)	(145,096)
Effect of exchange rate changes on cash, cash equivalents, and restricted cash	1,501	18
Cash, cash equivalents, and restricted cash at beginning of period	59,319	177,026
Cash, cash equivalents, and restricted cash at end of period	\$ 41,634	\$ 31,948

Use of 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 EBITDA and adjusted EBITDA margin:** We define adjusted EBITDA as net income (loss) adjusted to exclude interest income, income tax (expense) benefit, depreciation and amortization expense, stock-based compensation expense, acquisition and integration related costs, impairment and restructuring, 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. These items are discussed in more detail below the tables reconciling the GAAP to non-GAAP measures. Adjusted EBITDA margin is calculated as adjusted EBITDA divided by total revenues.
- **Adjusted cash usage:** We calculate cash usage as the total change in cash, cash equivalents, and restricted cash adjusted to include the net change from purchases, sales, and maturities of marketable securities (excluding any interest receivable). Adjusted cash usage is calculated as cash usage further adjusted to exclude cash payments related to 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 profit, adjusted gross margin, adjusted total operating expenses, and adjusted loss from operations:** We calculate these non-GAAP financial measures by including shipping and handling costs for product sales within cost of product revenue instead of within selling, general and administrative expenses. Additionally, we exclude 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 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. For example, adjusted EBITDA excludes a number of expense items that are included in net loss and adjusted cash usage excludes certain actual cash payments. As a result, positive adjusted EBITDA or positive adjusted cash usage may be achieved even where we record a significant net loss or reduction in our cash and marketable securities balances in accordance with U.S. GAAP.

Investors are encouraged to review the reconciliation of these non-GAAP financial measures to their most directly comparable GAAP financial measures set forth in the tables captioned "Reconciliation of GAAP to Non-GAAP Financial Measures" in the section below.

Additionally, we make certain forward-looking statements about our future financial performance that include non-GAAP financial measures, which are difficult to predict for future periods because the nature of the adjustments pertains to events that have not yet occurred. We do not forecast many of the excluded items for internal use and therefore information reconciling forward-looking non-GAAP financial measures to U.S. GAAP financial measures is not available without unreasonable effort and is not provided. The occurrence, timing, and amount of any of the items excluded from U.S. GAAP to calculate non-GAAP financial measures could significantly impact our U.S. GAAP results.

QUANTERIX CORPORATION
RECONCILIATIONS OF GAAP TO NON-GAAP FINANCIAL MEASURES

Reconciliation of Net Loss to Adjusted EBITDA (non-GAAP) and Adjusted EBITDA Margin (non-GAAP)
(Unaudited, amounts in thousands except percentages)

	Three Months Ended September 30,		Nine Months Ended September 30,	
	2025	2024	2025	2024
Net loss	\$ (33,517)	\$ (8,353)	\$ (84,033)	\$ (26,903)
Interest income	(1,448)	(3,535)	(7,408)	(11,165)
Income tax expense (benefit)	(2,480)	145	(5,466)	442
Depreciation and amortization	5,431	1,616	9,618	4,740
Stock-based compensation expense	5,469	4,657	16,303	15,150
Acquisition and integration related costs (1)	7,315	—	15,032	—
Earnout recorded as compensation expense (2)	229	—	8,129	—
Changes in contingent liabilities (3)	(58)	—	(3,952)	—
Impairment and restructuring costs (4)	7,174	—	14,844	—
Adjusted EBITDA (non-GAAP)	<u>\$ (11,885)</u>	<u>\$ (5,470)</u>	<u>\$ (36,933)</u>	<u>\$ (17,736)</u>
Total revenues	\$ 40,233	\$ 35,813	\$ 95,042	\$ 102,260
Adjusted EBITDA margin (non-GAAP) (adjusted EBITDA as a % of revenue)	(29.5)%	(15.3)%	(38.9)%	(17.3)%

(1) 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.

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

(3) Consists of fair value adjustments for contingent consideration liabilities related to acquisitions.

(4) Impairment charges for goodwill and an acquired lease facility not in use and severance and related benefit costs from the restructuring plan announced in 2025.

Reconciliation of Net Increase (Decrease) in Cash, Cash Equivalents, and Restricted Cash to Adjusted Cash Usage (non-GAAP)
(Unaudited, amounts in thousands)

	Three Months Ended September 30,		Nine Months Ended September 30,	
	2025	2024	2025	2024
Net increase (decrease) in cash, cash equivalents, and restricted cash	\$ (93,949)	\$ (18,032)	\$ (19,186)	\$ (145,096)
Effect of exchange rate changes on cash, cash equivalents, and restricted cash	45	371	1,501	18
Net change in marketable securities	(31,765)	14,331	(135,902)	117,282
Cash usage	(125,669)	(3,330)	(153,587)	(27,796)
Adjustments:				
Cash acquired from acquisitions	(16,799)	—	(16,822)	—
Acquisition and integration related payments (1)	120,310	—	134,387	—
Payments of restructuring costs (2)	6,095	—	7,075	—
Payments related to restatement costs (3)	—	—	1,102	—
Adjusted cash usage (non-GAAP)	\$ (16,063)	\$ (3,330)	\$ (27,845)	\$ (27,796)

(1) Represents cash payments towards acquisition and integration related activities, including the cash purchase price of an acquired business.

(2) Represents cash payments for severance and related benefits from the restructuring plan announced in 2025.

(3) Payment of costs associated with the restatement of previously issued financial statements that was completed at the end of 2024.

**Reconciliation of Gross Profit, Gross Margin, Total Operating Expenses and Loss from Operations to
Non-GAAP Financial Measures**
(Unaudited, amounts in thousands except percentages)

	Three Months Ended September 30,				Nine Months Ended September 30,			
	2025		2024		2025		2024	
Gross profit	\$	17,206	\$	20,153	\$	44,921	\$	60,935
Shipping and handling costs		(1,305)		(2,011)		(4,181)		(6,228)
Purchase accounting impact on inventory and property and equipment (1)		35		—		35		—
Amortization of acquired intangible assets (2)		2,532		—		2,993		—
Adjusted gross profit (non-GAAP)		18,468		18,142		43,768		54,707
Total revenues	\$	40,233	\$	35,813	\$	95,042	\$	102,260
Gross margin (gross profit as % of total revenues)		42.8%		56.3%		47.3%		59.6%
Adjusted gross margin (non-GAAP) (adjusted gross profit as % of total revenues)		45.9%		50.7%		46.1%		53.5%
Total operating expenses	\$	54,531	\$	31,901	\$	145,712	\$	98,782
Shipping and handling costs		(1,305)		(2,011)		(4,181)		(6,228)
Purchase accounting impact on property and equipment (1)		(212)		—		(212)		—
Amortization of acquired intangible assets (2)		(73)		—		(73)		—
Acquisition and integration related costs (3)		(7,315)		—		(15,032)		—
Earnout recorded as compensation expense (4)		(229)		—		(8,129)		—
Impairment and restructuring costs (5)		(7,174)		—		(14,844)		—
Adjusted total operating expenses (non-GAAP)	\$	38,223	\$	29,890	\$	103,241	\$	92,554
Loss from operations	\$	(37,325)	\$	(11,748)	\$	(100,791)	\$	(37,847)
Purchase accounting impact on inventory and property and equipment (1)		247		—		247		—
Amortization of acquired intangible assets (2)		2,605		—		3,066		—
Acquisition and integration related costs (3)		7,315		—		15,032		—
Earnout recorded as compensation expense (4)		229		—		8,129		—
Impairment and restructuring costs (5)		7,174		—		14,844		—
Adjusted loss from operations (non-GAAP)	\$	(19,755)	\$	(11,748)	\$	(59,473)	\$	(37,847)

(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 an acquired lease facility not in use and severance and related benefit costs from the restructuring plan announced in 2025.

CAUTIONARY STATEMENT REGARDING FORWARD-LOOKING STATEMENTS

Statements included in this press release that are not historical in nature or do not relate to current facts are intended to be, and are hereby identified as, forward-looking statements for purposes of the safe harbor provided by Section 27A of the Securities Act of 1933, as amended, and Section 21E of the Securities Exchange Act of 1934, as amended. Forward-looking statements include, among other things, statements about Quanterix's future business outlook, operations, strategy and financial performance, including statements related to our expectations about the development and commercialization of our products, about the benefits we may realize from the acquisition of Akoya Biosciences Inc., and under the header "2025 Full Year Business Outlook." Words and phrases such as "may," "approximately," "continue," "should," "expects," "projects," "anticipates," "is likely," "look ahead," "look forward," "believes," "will," "intends," "estimates," "strategy," "plan," "could," "potential," "possible" and variations of such words and similar expressions are intended to identify such forward-looking statements. Forward-looking statements are subject to certain risks and uncertainties that are difficult to predict with regard to, among other things, timing, extent, likelihood and degree of occurrence, which could cause actual results to differ materially from anticipated results. Such risks and uncertainties include, among others, the following possibilities with respect to Quanterix's future business, operations, strategy and financial performance: risks related to the impact of recent U.S. government policies, including reductions in federal research funding and increased tariffs; risks that we may not realize the expected benefits of our cost reduction actions; risks associated with the anticipated timing for launch of, and features of, Quanterix's next-generation instrument, Simoa One; risks that Quanterix may fail to realize the anticipated benefits and synergies of its recent acquisitions of Emission, Inc. and Akoya Biosciences Inc.; risk that integrating Quanterix's business with that of Akoya could be more difficult, costly or time-consuming than expected; Quanterix risks that Quanterix's estimates regarding expenses, future revenues, capital requirements, and needs for additional financing could be incorrect; risks related to the restatement of Quanterix's consolidated financial statements, including risks of increased costs and the increased possibility of legal proceedings and regulatory inquiries, sanctions, or investigation; risks related to Quanterix's ability to maintain effective internal control over financial reporting and disclosure controls and procedures, including its ability to remediate existing material weaknesses in its internal control over financial reporting and the timing of any such remediation; risks related to defects or other quality issues in Quanterix's products that could lead to unforeseen costs, product recalls, adverse regulatory actions, negative publicity and litigation; risks related to Quanterix's ability to retain and expand its customer base and achieve sufficient market acceptance of its products; and other factors that may affect future results of Quanterix. Additional factors that could cause results to differ materially from those described above can be found in the periodic reports filed by Quanterix with the SEC, including the "Risk Factors" sections contained therein, which are available on the SEC's website at www.sec.gov.

All forward-looking statements, expressed or implied, included in this press release are expressly qualified in their entirety by the cautionary statements contained or referred to herein. If one or more events related to these or other risks or uncertainties materialize, or if Quanterix's underlying assumptions prove to be incorrect, actual results may differ materially from what Quanterix anticipates. Quanterix cautions readers not to place undue reliance on any such forward-looking statements, which speak only as of the date they are made and are based on information available at that time. Quanterix does not assume any obligation to update or otherwise revise any forward-looking statements to reflect circumstances or events that occur after the date the forward-looking statements were made or to reflect the occurrence of unanticipated events except as required by federal securities laws.

Contact:

Joshua Young
(508) 846-3327
ir@quanterix.com

Third Quarter 2025 Earnings

November 10, 2025



Legal Information

CAUTIONARY STATEMENT REGARDING FORWARD-LOOKING STATEMENTS

Statements included in this press release that are not historical in nature or do not relate to current facts are intended to be, and are hereby identified as, forward-looking statements for purposes of the safe harbor provided by Section 27A of the Securities Act of 1933, as amended, and Section 21E of the Securities Exchange Act of 1934, as amended. Forward-looking statements include, among other things, statements about Quanterix's future business outlook, operations, strategy and financial performance, including statements related to our expectations about the development and commercialization of our products, about the benefits we may realize from the acquisition of Akoya Biosciences Inc., and under the header "2025 Full Year Business Outlook.". Words and phrases such as "may," "approximately," "continue," "should," "expects," "projects," "anticipates," "is likely," "look ahead," "look forward," "believes," "will," "intends," "estimates," "strategy," "plan," "could," "potential," "possible" and variations of such words and similar expressions are intended to identify such forward-looking statements. Forward-looking statements are subject to certain risks and uncertainties that are difficult to predict with regard to, among other things, timing, extent, likelihood and degree of occurrence, which could cause actual results to differ materially from anticipated results. Such risks and uncertainties include, among others, the following possibilities with respect to Quanterix's future business, operations, strategy and financial performance: risks related to the impact of recent U.S. government policies, including reductions in federal research funding and increased tariffs; risks that we may not realize the expected benefits of our cost reduction actions; risks associated with the anticipated timing for launch of, and features of, Quanterix's next-generation instrument, Simoa One; risks that Quanterix may fail to realize the anticipated benefits and synergies of its recent acquisitions of Emission, Inc. and Akoya Biosciences Inc.; risk that integrating Quanterix's business with that of Akoya could be more difficult, costly or time-consuming than expected; Quanterix risks that Quanterix's estimates regarding expenses, future revenues, capital requirements, and needs for additional financing could be incorrect; risks related to the restatement of Quanterix's consolidated financial statements, including risks of increased costs and the increased possibility of legal proceedings and regulatory inquiries, sanctions, or investigation; risks related to Quanterix's ability to maintain effective internal control over financial reporting and disclosure controls and procedures, including its ability to remediate existing material weaknesses in its internal control over financial reporting and the timing of any such remediation; risks related to defects or other quality issues in Quanterix's products that could lead to unforeseen costs, product recalls, adverse regulatory actions, negative publicity and litigation; risks related to Quanterix's ability to retain and expand its customer base and achieve sufficient market acceptance of its products; and other factors that may affect future results of Quanterix. Additional factors that could cause results to differ materially from those described above can be found in the periodic reports filed by Quanterix with the SEC, including the "Risk Factors" sections contained therein, which are available on the SEC's website at www.sec.gov.

All forward-looking statements, expressed or implied, included in this press release are expressly qualified in their entirety by the cautionary statements contained or referred to herein. If one or more events related to these or other risks or uncertainties materialize, or if Quanterix's underlying assumptions prove to be incorrect, actual results may differ materially from what Quanterix anticipates. Quanterix cautions readers not to place undue reliance on any such forward-looking statements, which speak only as of the date they are made and are based on information available at that time. Quanterix does not assume any obligation to update or otherwise revise any forward-looking statements to reflect circumstances or events that occur after the date the forward-looking statements were made or to reflect the occurrence of unanticipated events except as required by federal securities laws.

USE OF NON-GAAP FINANCIAL MEASURES

To supplement Quanterix's preliminary financial information presented on a GAAP basis, Quanterix has provided certain non-GAAP financial measures, including adjusted EBITDA, adjusted EBITDA margin, adjusted cash usage, adjusted gross profit, adjusted gross margin, adjusted total operating expenses, and adjusted loss from operations. Management uses 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. Management believes that presentation of these non-GAAP financial measures provides useful information to investors in assessing our operating performance within our industry and in order to allow comparability to the presentation of other companies in our industry. The non-GAAP financial measures presented herein should be considered in conjunction with, and not as a substitute for, the financial information presented in accordance with GAAP. For example, adjusted EBITDA excludes a number of expense items that are included in net loss and adjusted cash usage excludes certain actual cash payments. As a result, positive adjusted EBITDA or positive adjusted cash usage may be achieved even where we record a significant net loss or reduction in our cash and marketable securities balances in accordance with U.S. GAAP.

Investors are encouraged to review the reconciliation of these non-GAAP financial measures to their most directly comparable GAAP financial measures set forth herein. Quanterix does not forecast many of the excluded items for internal use and therefore information reconciling forward-looking non-GAAP financial measures to U.S. GAAP financial measures is not available without unreasonable effort and is not provided. The occurrence, timing, and amount of any of the items excluded from U.S. GAAP to calculate non-GAAP financial measures could significantly impact our U.S. GAAP results.

Please refer to our third quarter 2025 earnings release for additional discussion of non-GAAP financial measures.

Unless otherwise specified, all information contained herein is provided as of Sep 30, 2025.

Q3-25: Updates from Last Quarter

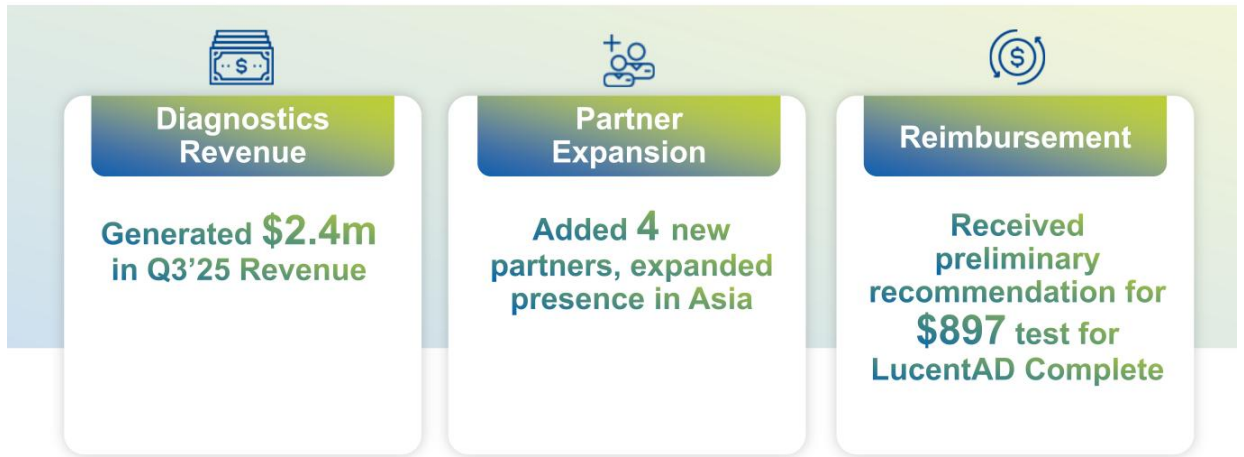
- ✓ **Delivered on revenue expectations in a difficult environment**
Reported third quarter 2025 revenue of \$40.2M
- ✓ **Moving fast and capturing significant synergies through Akoya integration**
Greater scale establishing foundation for future growth, profitability and cash flow
- ✓ **Making key investments for future growth**
YTD R&D spending is just under 30% of revenue
- ✓ **Carefully managing cash balance**
Expect to end 2025 with approximately \$120M of cash and no debt
- ✓ **Impact of synergies expected result in cash flow breakeven performance in 2026**
Synergy and cost actions of \$85M in 2026; ~80% of expense reductions already implemented

Synergies Leading to Cash Flow Breakeven in 2026

	Q2 2025	Q3 2025	Q4 2025	Q1 2026
Major Milestones	<ul style="list-style-type: none"> ✓ Pre-close cost actions in commercial and operations 	<ul style="list-style-type: none"> ✓ Complete physical consolidation ✓ Implement one commercial team ✓ Eliminate duplicate G&A 	<ul style="list-style-type: none"> ✓ Implement one manufacturing team • Combine Lab Services 	<ul style="list-style-type: none"> • Complete all Systems and Financial integration
Cost Reduction Implemented (Annualized)	\$29M	\$64M	\$67M	\$85M
Cost Reduction Realized (in the quarter)	\$3M	\$12M	\$15M	\$21M

\$67M of the total \$85M annualized cost reduction already implemented

We continue to make solid progress in advancing our Alzheimer's Diagnostics Business



Scientific Validation Driving Adoption (Pro Forma)

■ Spatial ■ Simoa
—●— Number of Drug Trial Projects (Simoa Only)



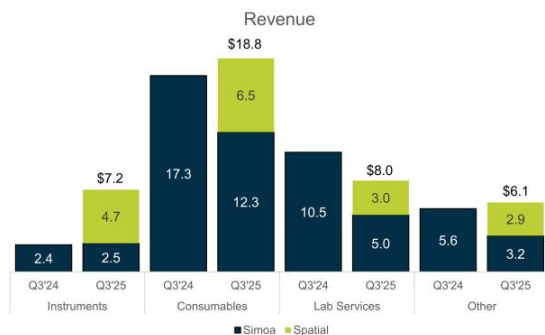
Q3'25 Results vs PYQ3

(in \$M)

	Q3 GAAP		Q3 Non-GAAP		
	2024	2025	2024	2025	Var %
Revenue	35.8	40.2	35.8	40.2	12%
Gross Margin \$	20.2	17.2	18.1	18.5	2%
Gross Margin %	56.3%	42.8%	50.7%	45.9%	-475 bps
Operating Expense	31.9	54.5	29.9	38.2	-22%
Operating Loss	-11.7	-37.3	-11.7	-19.8	-41%
Cash Usage Net of Cash Acquired	-3.3	-125.7	-3.3	-16.1	-79%
Adj EBITDA %			-15%	-29%	-1410 bps

Quanterix
Discovery Fueled by Ultra Sensitivity

	INSTRUMENTS	CONSUMABLES	ACCELERATOR	OTHER
Revenue Mix	18%	47%	20%	15%



2025 Guidance

Full Year Revenue: \$130 to \$135 million

~\$30 million contribution from Spatial.

Gross Margin

GAAP and Adjusted Gross Margin (Non-GAAP) between 45% to 47%.

\$34 to \$38 million of adjusted cash usage

\$136 million payment for Akoya and Emission acquisitions and restructuring costs, net of cash acquired

\$170 to \$174 million of total cash usage

'25 Exit Cash ~\$120 million; '26 Cash >\$100 million

Adjusted EBITDA (non-GAAP)

QUANTERIX CORPORATION RECONCILIATIONS OF GAAP TO NON-GAAP FINANCIAL MEASURES

Reconciliation of Net Loss to Adjusted EBITDA (non-GAAP) and Adjusted EBITDA Margin (non-GAAP) (Unaudited, amounts in thousands except percentages)

	Three Months Ended September 30,		Nine Months Ended September 30,	
	2025	2024	2025	2024
Net loss	\$ (33,517)	\$ (8,353)	\$ (84,033)	\$ (26,903)
Interest income	(1,448)	(3,535)	(7,408)	(11,165)
Income tax expense (benefit)	(2,480)	145	(5,466)	442
Depreciation and amortization	5,431	1,616	9,618	4,740
Stock-based compensation expense	5,469	4,657	16,303	15,150
Acquisition and integration related costs (1)	7,315	—	15,032	—
Earnout recorded as compensation expense (2)	229	—	8,129	—
Changes in contingent liabilities (3)	(58)	—	(3,952)	—
Impairment and restructuring costs (4)	7,174	—	14,844	—
Adjusted EBITDA (non-GAAP)	\$ (11,885)	\$ (5,470)	\$ (36,933)	\$ (17,736)
Total revenues	\$ 40,233	\$ 35,813	\$ 95,042	\$ 102,260
Adjusted EBITDA margin (non-GAAP) (adjusted EBITDA as a % of revenue)	(29.5)%	(15.3)%	(38.9)%	(17.3)%

(1) 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.

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

(3) Consists of fair value adjustments for contingent consideration liabilities related to acquisitions.

(4) Impairment charges for goodwill and an acquired lease facility not in use and severance and related benefit costs from the restructuring plan announced in 2025.

Adjusted Cash Usage (non-GAAP)

Reconciliation of Net Increase (Decrease) in Cash, Cash Equivalents, and Restricted Cash to Adjusted Cash Usage (non-GAAP)

(Unaudited, amounts in thousands)

	Three Months Ended September 30,		Nine Months Ended September 30,	
	2025	2024	2025	2024
Net increase (decrease) in cash, cash equivalents, and restricted cash	\$ (93,949)	\$ (18,032)	\$ (19,186)	\$ (145,096)
Effect of exchange rate changes on cash, cash equivalents, and restricted cash	45	371	1,501	18
Net change in marketable securities	(31,765)	14,331	(135,902)	117,282
Cash usage	(125,669)	(3,330)	(153,587)	(27,796)
Adjustments:				
Cash acquired from acquisitions	(16,799)	—	(16,822)	—
Acquisition and integration related payments (1)	120,310	—	134,387	—
Payments of restructuring costs (2)	6,095	—	7,075	—
Payments related to restatement costs (3)	—	—	1,102	—
Adjusted cash usage (non-GAAP)	\$ (16,063)	\$ (3,330)	\$ (27,845)	\$ (27,796)

(1) Represents cash payments towards acquisition and integration related activities, including the cash purchase price of an acquired business.

(2) Represents cash payments for severance and related benefits from the restructuring plan announced in 2025.

(3) Payment of costs associated with the restatement of previously issued financial statements that was completed at the end of 2024.

Additional Non-GAAP Financial Measures

Reconciliation of Gross Profit, Gross Margin, Total Operating Expenses and Loss from Operations to Non-GAAP Financial Measures

(Unaudited, amounts in thousands except percentages)

	Three Months Ended September 30,		Nine Months Ended September 30,	
	2025	2024	2025	2024
Gross profit	\$ 17,206	\$ 20,153	\$ 44,921	\$ 60,935
Shipping and handling costs	(1,305)	(2,011)	(4,181)	(6,228)
Purchase accounting impact on inventory and property and equipment (1)	35	—	35	—
Amortization of acquired intangible assets (2)	2,532	—	2,993	—
Adjusted gross profit (non-GAAP)	\$ 18,468	\$ 18,142	\$ 43,768	\$ 54,707
Total revenues	\$ 40,233	\$ 35,813	\$ 95,042	\$ 102,260
Gross margin (gross profit as % of total revenues)	42.8%	56.3%	47.3%	59.6%
Adjusted gross margin (non-GAAP) (adjusted gross profit as % of total revenues)	45.9%	50.7%	46.1%	53.5%
Total operating expenses	\$ 54,531	\$ 31,901	\$ 145,712	\$ 98,782
Shipping and handling costs	(1,305)	(2,011)	(4,181)	(6,228)
Purchase accounting impact on property and equipment (1)	(212)	—	(212)	—
Amortization of acquired intangible assets (2)	(73)	—	(73)	—
Acquisition and integration related costs (3)	(7,315)	—	(15,032)	—
Earnout recorded as compensation expense (4)	(229)	—	(8,129)	—
Impairment and restructuring costs (5)	(7,174)	—	(14,844)	—
Adjusted total operating expenses (non-GAAP)	\$ 38,223	\$ 29,890	\$ 103,241	\$ 92,554
Loss from operations	\$ (37,325)	\$ (11,748)	\$ (100,791)	\$ (37,847)
Purchase accounting impact on inventory and property and equipment (1)	247	—	247	—
Amortization of acquired intangible assets (2)	2,605	—	3,066	—
Acquisition and integration related costs (3)	7,315	—	15,032	—
Earnout recorded as compensation expense (4)	229	—	8,129	—
Impairment and restructuring costs (5)	7,174	—	14,844	—
Adjusted loss from operations (non-GAAP)	\$ (19,755)	\$ (11,748)	\$ (59,473)	\$ (37,847)

(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 an acquired lease facility not in use and severance and related benefit costs from the restructuring plan announced in 2025.

