



# First Integrated Solution for Ultra-Sensitive Detection of Blood and Tissue Based Protein Biomarkers

Uniquely positioned to speed up market development of new diagnostic testing solutions

January 10, 2025

# Legal Information

#### **IMPORTANT ADDITIONAL INFORMATION**

In connection with the proposed transaction, Quanterix Corporation ("Quanterix" or "QTRX") will file with the U.S. Securities and Exchange Commission (the "SEC") a registration statement on Form S-4 (the "registration statement"), which will contain a joint proxy statement of Quanterix and Akoya Biosciences, Inc. ("Akoya" or "AKYA") and a prospectus of Quanterix (the "joint proxy statement/prospectus"), and each of Quanterix and Akoya may file with the SEC other relevant documents regarding the proposed transaction. INVESTORS AND SECURITY HOLDERS ARE URGED TO READ THE REGISTRATION STATEMENT AND THE JOINT PROXY STATEMENT/PROSPECTUS CAREFULLY AND IN THEIR ENTIRETY AND ANY OTHER RELEVANT DOCUMENTS FILED WITH THE SEC BY QUANTERIX AND AKOYA, AS WELL AS ANY AMENDMENTS OR SUPPLEMENTS TO THOSE DOCUMENTS WHEN THEY BECOME AVAILABLE BECAUSE THEY WILL CONTAIN IMPORTANT INFORMATION ABOUT QUANTERIX, AKOYA AND THE PROPOSED TRANSACTION. A definitive copy of the joint proxy statement/prospectus will be mailed to Quanterix and Akoya or from the SEC's website when they are filed. The documents filed by Quanterix with the SEC may be obtained free of charge at Quanterix's website, at www.quanterix.com, or by requesting them by mail at Akoya Investor Relations, 100 Campus Drive, 6th floor, Marlboro, MA 01752.

#### PARTICIPANTS IN THE SOLICITATION

Quanterix and Akoya and certain of their respective directors and executive officers may be deemed to be participants in the solicitation of proxies from the stockholders of Quanterix or Akoya in respect of the proposed transaction. Information about Quanterix's directors and executive officers is available in Quanterix's proxy statement dated April 15, 2024, for its 2024 Annual Meeting of Stockholders, and other documents filed by Quanterix with the SEC. Information about Akoya's directors and executive officers is available in Akoya's proxy statement dated April 23, 2024, for its 2024 Annual Meeting of Stockholders, and other documents filed by Akoya with the SEC. Other information regarding the persons who may, under the rules of the SEC, be deemed participants in the proxy solicitation and a description of their direct and indirect interests, by security holdings or otherwise, will be contained in the joint proxy statement/prospectus and other relevant materials to be filed with the SEC regarding the proposed transaction when they become available. Investors should read the joint proxy statement/prospectus carefully when it becomes available before making any voting or investment decisions. You may obtain free copies of these documents from Quanterix or Akoya as indicated above.

#### **NO OFFER OR SOLICITATION**

This presentation shall not constitute an offer to sell or the solicitation of an offer to buy any securities or a solicitation of any vote or approval with respect to the proposed merger of Quanterix and Akoya, nor shall there be any sale of securities in any jurisdiction in which such offer, solicitation or sale would be unlawful prior to registration or qualification under the securities laws of any such jurisdiction. No offering of securities shall be made except by means of a prospectus meeting the requirements of Section 10 of the U.S. Securities Act of 1933, as amended.

#### CAUTIONARY STATEMENT REGARDING FORWARD-LOOKING STATEMENTS

Statements included in this presentation which 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 are based on, among other things, projections as to the anticipated benefits of the proposed transaction. Mords 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. Quanterix and Akoya caution readers that 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 statements are based transaction of a such adversely affect the combined company or the expected benefits of the proposed transaction as well as a result of the proposed transaction and stockholder approvals or to satisfy any of the other conditions to the proposed transaction on a timely basis or at all, the possibility that the anticipated benefits of the proposed transaction may be more expensive to complete than anticipated, diversion of the strategy of the strategy and competitive factors in the areas where Quanterix and Akoya do business; the possibility that the anticipated benefits and synergies of the association are not realized benefits of the proposed transaction and the closing of the proposed transaction and the closing of the proposed transactin the definitive merger agreement entered into be

### **Quanterix**<sup>®</sup> Discovery Fueled by Ultra-Sensitivity

# Legal Information

### CAUTIONARY STATEMENT REGARDING FORWARD-LOOKING STATEMENTS (CONTD.)

All forward-looking statements, expressed or implied, included in this presentation 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 or Akoya's underlying assumptions prove to be incorrect, actual results may differ materially from what Quanterix and Akoya anticipate. Quanterix and Akoya caution 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. Neither Quanterix nor Akoya assumes 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 FINANCIAL PROJECTIONS

This presentation contains projected financial information with respect to the combined company that will result from the proposed transaction, if consummated. Such projected financial information constitutes forward-looking information, is for illustrative purposes only and should not be relied upon as necessarily being indicative of future results. The assumptions and estimates underlying such projected financial information are inherently uncertain and are subject to a wide variety of significant business, economic, competitive and other risks and uncertainties that could cause actual results to differ materially from those contained in the projected financial information. Actual results may differ materially from the results contemplated by the projected financial information contained in this presentation, and the inclusion of such information in this presentation should not be regarded as a representation by any person that the results reflected in such forecasts will be achieved. Neither Quanterix's nor Akoya's independent auditors have audited, reviewed, compiled or performed any procedures with respect to the projections for purposes of their inclusion in this presentation, and accordingly, neither of them expressed an opinion or provided any other form of assurance with respect thereto for purposes of this presentation.

### 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 non-GAAP gross margin. Management uses these non-GAAP measures to evaluate our operating performance in 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 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 information presented herein should be considered in conjunction with, and not as a substitute for, the financial information presented in accordance with GAAP.

Unless otherwise specified, all information contained herein is provided as of September 30, 2024.



# **Transaction Overview**

Consideration	All-stock transaction AKYA shareholders to receive <b>0.318x</b> QTRX shares per share of AKYA
Ownership	QTRX shareholders to own <b>~70%</b> of combined entity AKYA shareholders to own <b>~30%</b> of combined entity
Governance	AKYA will nominate 2 candidates to join QTRX's Board of Directors, with the total Board size at 9 members Masoud Toloue will serve as Chief Executive Officer and Vandana Sriram will serve as Chief Financial Officer; combined company will continue to <b>operate under the Quanterix name</b>
Timing	Expected transaction close in Q2 2025 Subject to shareholder approvals and regulatory approvals and other customary closing conditions Shareholders owning more than <b>50%</b> of AKYA common stock have entered into a voting agreement in support of the transaction



### Accelerated Path to Scale and Profitability with Focus on High Growth Markets

01	Growth	Attractive growth expected in a \$5B <sup>1</sup> addressable market
	Scale	Combined revenue of ~\$220M <sup>2</sup> Expected strong double-digit organic revenue growth by 2026
	Efficiency	Significant annual cost synergies of <b>~\$40M</b> by end of 2026 ( <b>~\$20M</b> by end of year 1 post close)
	Acceleration	Path to cash flow positive in 2026
	Strong Runway	Significant combined balance sheet with ~\$ 175M <sup>3</sup> in cash expected at time of closing
		<ol> <li>Management's estimate; Excludes Dx and CDx</li> <li>TTM ended Q3 2024</li> <li>Assumes that the deal closes in Q2 2025.</li> </ol>



### Quanterix Today Leader in Ultra-Sensitive Biomarker Detection

Proven track record of

operating discipline



Unmatched Technology



## Quanterix

Early penetration in multi-billion diagnostics



CLIA lab running validated neuro LDTs

**\$134-138M** 2024 Revenue Guidance<sup>1</sup>

**3,100+** 

>1,000 Installed Base

**51-55%** 2024 Non-GAAP Gross Margin<sup>1</sup> Guidance

North America: 65% | EMEA: 26% | APAC: 9% Revenue by Geography

1. Guidance as of November 12, 2024

2. GAAP gross margin guidance range of 57-61%. Projected GAAP gross margin does not include shipping and handling costs, which include freight and other activities costs associated with product shipments. Non-GAAP gross margin includes these shipping and handling costs.



### Akoya Biosciences Leading the Spatial Biology Revolution



Established leader with largest install base in market



Fastest and most robust platform for tissue biomarker detection

Recurring revenue through reagents, software and services



Growing clinical business via Clinical Services and future CDx **\$80-85M** 2024 Revenue Guidance

1,600+

~1,300 Installed Base

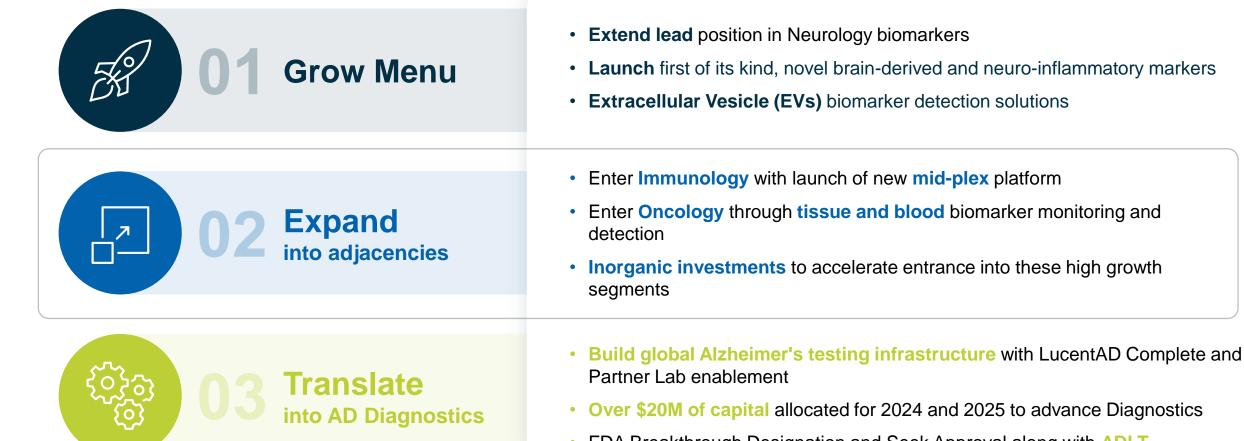
**59%** Q1-Q3'24 Non-GAAP Gross Margin<sup>1</sup>

North America: 58% | EMEA: 26% | APAC: 16% Revenue by Geography

1. Non-GAAP gross margin does not include write-offs from discontinued legacy products in the first quarter of 2024. Q1-Q3'24 GAAP gross margin was 56%.



# Continued Execution of Quanterix's Growth Plan



 FDA Breakthrough Designation and Seek Approval along with ADLT reimbursement



# Immediate Opportunity to Increase Serviceable Addressable Market



Combined Instrument

Installed-Base

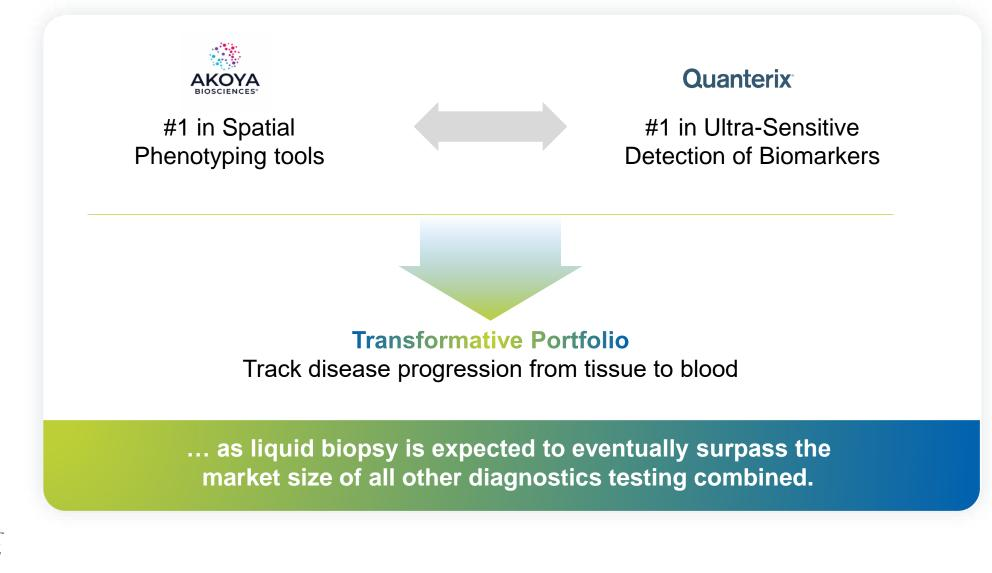
 Combined commercial platform and cross-selling opportunities enable further market development

\* Management's estimate; Excludes Dx and CDx

~2,300

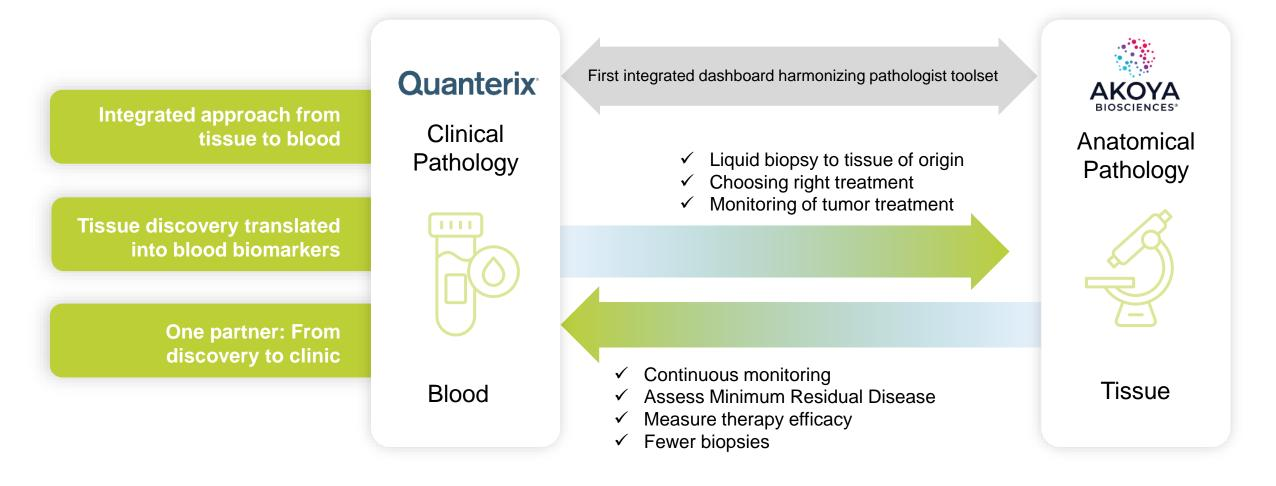


# Uniquely positioned to accelerate market development of new liquid biopsy tests ...





# Uniting Clinical and Anatomical Pathology





### Complementary Technology and Service Offerings Expand Commercial Reach



Discovery Fueled by Ultra-Sensitivit

12

# Combined Company: Strong Double-Digit Organic Growth by 2026

Uniquely positioned to speed market development of new liquid biopsy tests

### **Near-term Growth Drivers**

Menu

**Services** 

Apply consumables product development engine to accelerate growth of new biomarker pairs across tissue and blood. Suite of synergistic biomarkers already in development. **Over past 2 years, Quanterix grew reagents by 24%.** 

Replicate Quanterix's Accelerator model to expand Akoya's Tissue testing services portfolio. Over past 2 years, Quanterix's Accelerator services grew 33%

**Doubling**<br/>of footprintInstall base expansion creates new cross-selling opportunities.<br/>Combined install base of ~2,300 instruments

With upside opportunity in a potential multi-billion Alzheimer's and Companion Diagnostics segment



# Significant Cost Synergies

COSTS			
Cost of Goods	$\oslash$	Combined reagent manufacturing capabilities / infrastructure to improve overall margins	~
Commercial	$\oslash$	Integrated commercial organization capable of cross selling products / solutions into a large common Biomarker lab customer base	In Qu
General and Admin	$\oslash$	Overlap of Corporate, Supply chain and other Support functions offer significant optimization opportunity	op tra
FOOTPRINT			•
Infrastructure	$\oslash$	Geographical proximity of Operations and CLIA lab of both entities (in MA) enables efficient consolidation and integration	lm co

# \$20M

n first 12 months

Annual cost synergies by end of 2026

~\$40M

Quanterix has a proven track record of operating discipline; 2022 Corporate transformation delivered<sup>1</sup>:

- Margin expansion of 1,850 bps
- /2/3<sup>rd</sup> reduction in cash burn

Improvements have resulted in 7 consecutive quarters of double-digit growth

1. Improvement from FY22 to FY23



# Uniquely Positioned to Deliver Value

# Transformative **Portfolio**

Creates first integrated platform of liquid and tissue biomarkers

High Growth Expanded customer relationships & continuum of solution offerings maximize cross-selling opportunities Unlocks a high-growth \$5B<sup>1</sup> addressable market across Neurology, Immunology and Oncology

Combined revenue of ~\$220M<sup>2</sup>

Expected strong **double-digit** organic revenue growth by 2026

### Accelerated Returns

Significant synergy opportunities across Operations and SG&A Significant annual cost synergies of ~\$40M by end of 2026

Path to free cash flow positive in 2026

1. Management's estimate; Excludes Dx and CDx

2. TTM ended Q3 2024

