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

SCHEDULE 14A
(Rule 14a-101)

SCHEDULE 14A INFORMATION
Proxy Statement Pursuant to Section 14(a) of
the Securities Exchange Act of 1934

Filed by the Registrant

Filed by a party other than the Registrant

Check the appropriate box:

- Preliminary Proxy Statement
- Confidential, for Use of the Commission Only (as permitted by Rule 14a-6(e)(2))
- Definitive Proxy Statement
- Definitive Additional Materials
- Soliciting Material Pursuant to §240.14a-12

QUANTERIX CORPORATION

(Name of Registrant as Specified In Its Charter)

KENT LAKE PARTNERS LP
KENT LAKE PR LLC
BENJAMIN NATTER
ALEXANDER G. DICKINSON
BRUCE FELT
HAKAN SAKUL

(Name of Person(s) Filing Proxy Statement, if other than the Registrant)

Payment of Filing Fee (Check the appropriate box):

- No fee required.
- Fee paid previously with preliminary materials.
- Fee computed on table in exhibit required by Item 25(b) per Exchange Act Rules 14a-6(i)(1) and 0-11.
-

Kent Lake Partners LP (“Kent Lake Partners”), together with the other Participants (as defined below), intends to file a preliminary proxy statement and an accompanying GOLD universal proxy card with the Securities and Exchange Commission (“SEC”) to be used to solicit votes for, among other matters, the election of its slate of highly-qualified director nominees at the 2025 annual meeting of stockholders of Quanterix Corporation, a Delaware corporation (the “Company”).

On March 11, 2025, Kent Lake Partners issued the following press release and Investor Presentation, a copy of which is attached hereto as Exhibit 1 and is incorporated herein by reference.

Kent Lake Issues Investor Presentation Detailing Opposition to Quanterix’s Proposed Merger with Akoya Biosciences

Believes Transaction Significantly Undervalues Quanterix

Highlights that Quanterix’s Board is Essentially Rescuing Akoya at the Direct Expense of Quanterix Shareholders

Transaction Would Introduce Unnecessary Risk and Divert Cash and Management Focus Away from Quanterix’s Alzheimer’s Diagnostics Program, a Core Growth Opportunity

Calls on Quanterix Board to Answer Pressing Questions around Valuation, Process and Governance Issues in Connection with Transaction

Urges Shareholders to Vote Against Merger with Akoya to Protect the Future Value of Their Investment

RINCON, Puerto Rico – March 11, 2025 – Kent Lake PR LLC (“Kent Lake”) a holder of 7.25% of the outstanding common stock of Quanterix Corporation (“Quanterix” or the “Company”) (NASDAQ: QTRX), today issued a presentation to Quanterix shareholders in connection to Kent Lake’s opposition to the Company’s proposed transaction (the “Transaction”) with Akoya Biosciences (“Akoya”) (NASDAQ: AKYA).

Kent Lake’s full presentation is available at https://kentlake.s3.us-west-1.amazonaws.com/QTRX-AKYA_KLPRVoteNoDeck.pdf.

Highlights of the presentation include:

- **The Transaction is a bad deal for Quanterix shareholders** – The Transaction inexplicably values Akoya nearly 5x higher than Quanterix, despite Quanterix’s superior mix, growth and market opportunity. It is essentially a bailout of Akoya. Quanterix’s shares have declined nearly 40% since the deal announcement.
 - **Quanterix ended up bidding against itself for Akoya** – As revealed in the S-4 related to the Transaction, other motivated, strategic potential buyers of Akoya walked away due to concerns around profitability, cash burn and leverage, leaving Quanterix as the only bidder. Reading the S-4 raises serious issues about the process that led to this Transaction.
 - **The merger introduces unnecessary risk for Quanterix** – Expected National Institute of Health funding cuts will further strain Akoya’s business, making Akoya’s balance sheet situation untenable and reinforcing Kent Lake’s belief that Akoya currently has little or no positive equity value.
 - **Management is entrenched and conflicted** – Certain directors remain on the Quanterix Board of Directors (the “Board”) despite failing to obtain majority voting from shareholders, while others appear to have undisclosed conflicts of interest that compromise their fiduciary duty to Quanterix shareholders, such as Dr. Ivana Magovčević-Liebisch’s director position at a Company with a material interest in Akoya.
-

- **The Transaction includes a provision where Quanterix would provide Akoya with \$30 million in bridge financing absent any clear shareholder mandate** – This is an egregious breach of the Board’s fiduciary duty to shareholders.
- **Quanterix has significant potential as a standalone company** – The Transaction will divert focus away from Quanterix’s transformational Alzheimer’s testing opportunity, which, if pursued, could unlock substantial value.

In Kent Lake’s view, the incumbent Board can no longer be trusted to represent shareholders’ best interests. That is why Kent Lake recently nominated three highly qualified, independent candidates for election to Quanterix’s Board at the 2025 Annual Meeting of Shareholders. There are clear, superior alternatives for Quanterix on a standalone basis, and Kent Lake believes that abandoning the ill-fated Transaction would allow shareholders to realize greater value creation than combining with Akoya.

About Kent Lake

Kent Lake Partners LP is an investment fund founded by Ben Natter in 2019 with a focus on small and mid-capitalization public equities, particularly in the healthcare space. Mr. Natter has over a decade of successful public healthcare equity investing experience.

Certain Information Concerning the Participants

Kent Lake Partners LP (“Kent Lake Partners”), together with the other Participants (as defined below), intends to file a preliminary proxy statement and an accompanying GOLD universal proxy card with the Securities and Exchange Commission (“SEC”) to be used to solicit votes for, among other matters, the election of its slate of highly-qualified director nominees at the 2025 annual meeting of stockholders of Quanterix Corporation, a Delaware corporation (the “Company”).

KENT LAKE STRONGLY ADVISES ALL SHAREHOLDERS OF THE COMPANY TO READ THE PROXY STATEMENT AND OTHER PROXY MATERIALS, INCLUDING ITS GOLD PROXY CARD, AS THEY BECOME AVAILABLE BECAUSE THEY WILL CONTAIN IMPORTANT INFORMATION. SUCH PROXY MATERIALS WILL BE AVAILABLE AT NO CHARGE ON THE SEC’S WEB SITE AT [HTTP://WWW.SEC.GOV](http://www.sec.gov). IN ADDITION, THE PARTICIPANTS IN THIS PROXY SOLICITATION WILL PROVIDE COPIES OF THE PROXY STATEMENT WITHOUT CHARGE, WHEN AVAILABLE, UPON REQUEST. REQUESTS FOR COPIES SHOULD BE DIRECTED TO THE PARTICIPANTS’ PROXY SOLICITOR.

The participants in the proxy solicitation are currently anticipated to be Kent Lake Partners, Kent Lake PR LLC (“Kent Lake PR”) and Benjamin Natter (collectively, the “Kent Lake Parties”); and Alexander G. Dickinson, Bruce Felt and Hakan Sakul (the “Kent Lake Nominees” and collectively with the Kent Lake Parties, the “Participants”).

As of the date hereof, Kent Lake Partners directly beneficially owned 2,793,405 shares of the Company’s Common Stock, \$0.001 par value per share (the “Common Stock”). Kent Lake PR, as the investment adviser and as the general partner to Kent Lake Partners, may be deemed to beneficially own the 2,793,405 shares of Common Stock beneficially owned by Kent Lake Partners. Mr. Natter, as the Managing Member of Kent Lake PR, may be deemed to beneficially own the 2,793,405 shares of Common Stock beneficially owned by Kent Lake Partners. None of the Kent Lake Nominees beneficially own any shares of Common Stock. All of the foregoing information is as of the date hereof unless otherwise disclosed.

Investor Contacts

Ben Natter, 415-237-0007
info@kentlakecap.com

Saratoga Proxy Consulting LLC
John Ferguson / Ann Marie Mellone
(212) 257-1311 / (888) 368-0379

info@saratogaproxy.com

Media Contact

Joe Germani / Kendall Heebink
Longacre Square Partners
KentLake@longacresquarepartners.com



KENT LAKE PR LLC

Urges Quanterix Shareholders to Vote **AGAINST**
the Proposed Merger with Akoya Biosciences

THIS IS NOT A SOLICITATION OF AUTHORITY TO VOTE YOUR PROXY. DO NOT SEND US YOUR PROXY CARD. KENT LAKE IS NOT ABLE TO VOTE YOUR PROXY, NOR DOES THIS COMMUNICATION CONTEMPLATE SUCH AN EVENT.

Disclaimer

The materials contained herein (the "Materials") represent the opinions of Kent Lake PR LLC and its affiliates (collectively, "Kent Lake") and are based on publicly available information with respect to the Quanterix Corporation, a Delaware corporation (the "Company"). Kent Lake recognizes that there may be confidential information in the possession of the Company that could lead it or others to disagree with Kent Lake's conclusions. Kent Lake reserves the right to change any of its opinions expressed herein at any time as it deems appropriate and disclaims any obligation to notify the market or any other party of any such changes. Kent Lake disclaims any obligation to update the information or opinions contained herein. Certain financial projections and statements made herein have been derived or obtained from filings made with the Securities and Exchange Commission ("SEC") or other regulatory authorities and from other third party reports. There is no assurance or guarantee with respect to the prices at which any securities of the Company will trade, and such securities may not trade at prices that may be implied herein. The estimates, projections and potential impact of the opportunities identified by Kent Lake herein are based on assumptions that Kent Lake believes to be reasonable as of the date of the Materials, but there can be no assurance or guarantee that actual results or performance of the Company will not differ, and such differences may be material. The Materials are provided merely as information and are not intended to be, nor should they be construed as, an offer to sell or a solicitation of an offer to buy any security.

Certain members of Kent Lake currently beneficially own, and/or have an economic interest in, securities of the Company. It is possible that there will be developments in the future (including changes in price of the Company's securities) that cause one or more members of Kent Lake from time to time to sell all or a portion of their holdings of the Company in open market transactions or otherwise (including via short sales), buy additional securities (in open market or privately negotiated transactions or otherwise), or trade in options, puts, calls or other derivative instruments relating to some or all of such securities. To the extent that Kent Lake discloses information about its position or economic interest in the securities of the Company in the Materials, it is subject to change and Kent Lake expressly disclaims any obligation to update such information.

The Materials contain forward-looking statements. All statements contained herein that are not clearly historical in nature or that necessarily depend on future events are forward-looking, and the words "anticipate," "believe," "expect," "potential," "opportunity," "estimate," "plan," "may," "will," "projects," "targets," "forecasts," "seeks," "could," and similar expressions are generally intended to identify forward-looking statements. The projected results and statements contained herein that are not historical facts are based on current expectations, speak only as of the date of the Materials and involve risks, uncertainties and other factors that may cause actual results, performance or achievements to be materially different from any future results, performance or achievements expressed or implied by such projected results and statements. Assumptions relating to the foregoing involve judgments with respect to, among other things, future economic, competitive and market conditions and future business decisions, all of which are difficult or impossible to predict accurately and many of which are beyond the control of Kent Lake. Although Kent Lake believes that the assumptions underlying the projected results or forward-looking statements are reasonable as of the date of the Materials, any of the assumptions could be inaccurate and therefore, there can be no assurance that the projected results or forward-looking statements included herein will prove to be accurate. In light of the significant uncertainties inherent in the projected results and forward-looking statements included herein, the inclusion of such information should not be regarded as a representation as to future results or that the objectives and strategic initiatives expressed or implied by such projected results and forward-looking statements will be achieved. Kent Lake will not undertake and specifically declines any obligation to disclose the results of any revisions that may be made to any projected results or forward-looking statements herein to reflect events or circumstances after the date of such projected results or statements or to reflect the occurrence of anticipated or unanticipated events.

Unless otherwise indicated herein, Kent Lake has not sought or obtained consent from any third party to use any statements, photos or information indicated herein as having been obtained or derived from statements made or published by third parties. Any such statements or information should not be viewed as indicating the support of such third party for the views expressed herein. No warranty is made as to the accuracy of data or information obtained or derived from filings made with the SEC by the Company or from any third-party source. All trade names, trademarks, service marks, and logos herein are the property of their respective owners who retain all proprietary rights over their use.

Executive Summary: Why Quanterix Shareholders Should Vote Against

Merger is a Bailout of Akoya at Quanterix Shareholders' Expense	Quanterix Board is Entrenched and Conflicted	Merger Introduces Unnecessary Risk for Quanterix	Distracts from Core Growth Opportunity
<ul style="list-style-type: none"> ▪ The deal assigns Quanterix an enterprise value of only ~\$40M, while valuing Akoya at ~\$190M – nearly 5x higher than Quanterix – despite Quanterix's superior mix, growth, and market opportunity¹ ▪ Quanterix bid against itself despite clear issues with Akoya's financials that caused all other bidders to walk away ▪ A standalone Akoya would require dilutive financing to survive – Quanterix's board is rescuing Akoya at the direct expense of Quanterix investors 	<ul style="list-style-type: none"> ▪ Quanterix shareholders have a clear lack of confidence in Board Chairman Dr. Martin Madaus – he failed to obtain support from a majority of voters at the 2024 annual meeting yet did not resign ▪ Quanterix director Dr. Ivana Magovčević-Liebisch's director position at a Company with a material interest in Akoya suggests a significant undisclosed conflict of interest ▪ We have serious reservations about the process that led to this deal 	<ul style="list-style-type: none"> ▪ Post-merger, Quanterix's cash runway would shrink from ~6.5 years to ~2.5 years², greatly increasing the risk of future dilutive capital raises and undermining its previously robust balance sheet ▪ Anticipated NIH funding cuts would likely further worsen Akoya's balance sheet situation ▪ Quanterix shares deserve a higher revenue multiple in a stand-alone scenario than in a combined company scenario 	<ul style="list-style-type: none"> ▪ Quanterix is at a critical juncture with its Alzheimer's diagnostics program, and the merger would divert focus and resources away from FDA approval, reimbursement, and commercialization ▪ Quanterix should be concentrating on its biggest growth opportunity... not absorbing a struggling Akoya ▪ Market reaction to the deal speaks volumes – Quanterix's shares have declined more than 30% since the deal announcement
<p>To protect their investment, Kent Lake urges its fellow Quanterix shareholders to <u>vote against</u> a highly dilutive, risky merger with Akoya Biosciences that devalues and distracts from Quanterix's standalone potential.</p>			

Source: Capital IQ, and public filings

1) Uses 2/10/25 \$7.83 Quanterix share price and consideration allocation as disclosed in the S-4 Filing.

2) Calculated as LTM CFO less CA, divided by Net Cash. Quanterix cash reflects \$296M unrestricted and restricted cash and marketable securities at 9/30/24, adjusted for \$4.4M 4Q24 outflows and \$20M related to the Emission acquisition.

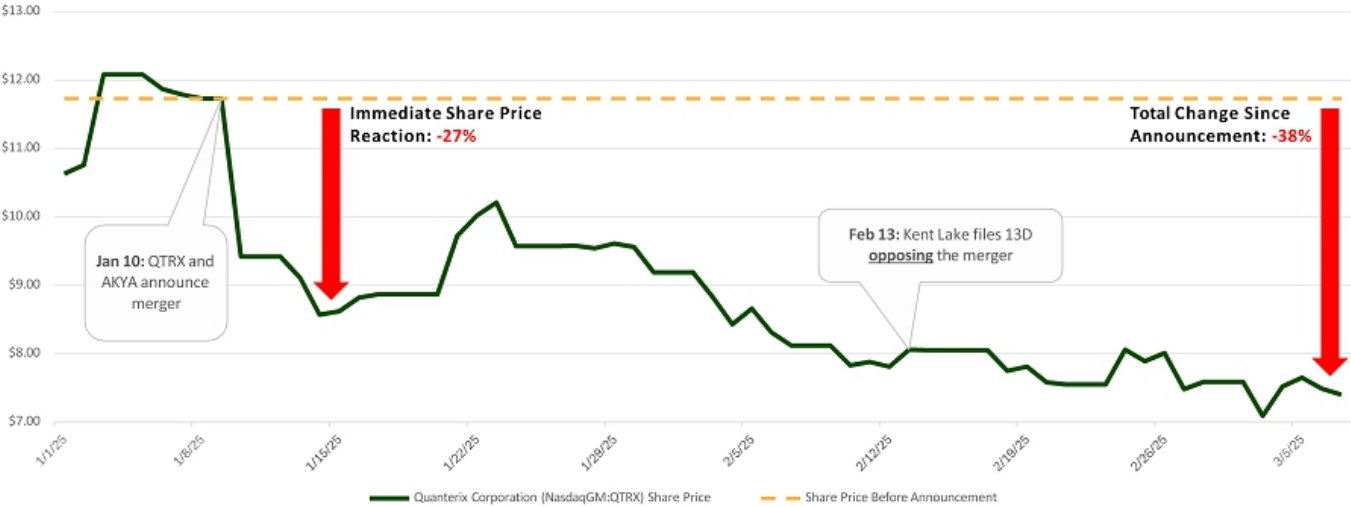
3) Per Akoya CEO on 2Q24 CC: "In aggregate, we implemented a workforce reduction of approximately 35% compared to end of year 2023."

Valuation and Market Reaction



QTRX's Value Declined Significantly Since the Announcement

Since announcing the merger, QTRX's share price has declined **nearly 40%**. Why do management and the Board insist on a transaction that the owners of the company are so clearly against?



Source: Capital IQ, and public filings
Share price calculated as QTRX daily close price dated March 7, 2025

Structurally Unfair Transaction for **QTRX Shareholders Who Receive a Mere 18% of Combined Company Enterprise Value**, While AKYA Shareholders Receive 82%!

Dr. Toloue and the QTRX Board, *which set of shareholders are you really working for?*

AKYA IMPLIED VALUE - S4 DATE (2/10/25)	QTRX	Implied AKYA
Share Price (2/10/25)	\$ 7.83	0.318x
Total Akoya FDSO		51,448
QTRX Equivalent FDSO	39,745	16,361
Market Capitalization	\$ 311,206	\$ 128,105
Replacement of AKYA Options + Frac. Shares		\$ 13,698
Total Value of Consideration		\$ 141,803
Net Cash	\$ 269,123	\$ (46,993)
Enterprise Value	\$ 42,083	\$ 188,796
% of Pro Forma TEV	18%	82%
<i>Forward Revenue (CIQ Est.)</i>		
2024E Revenue	\$ 137,200	\$ 81,790
2025E Revenue	\$ 154,701	\$ 91,571
2026E Revenue	\$ 179,867	\$ 105,306
TEV / 2024E Revenue	0.3x	2.3x
TEV / 2025E Revenue	0.3x	2.1x
TEV / 2026E Revenue	0.2x	1.8x

Source: Capital IQ, and public filings

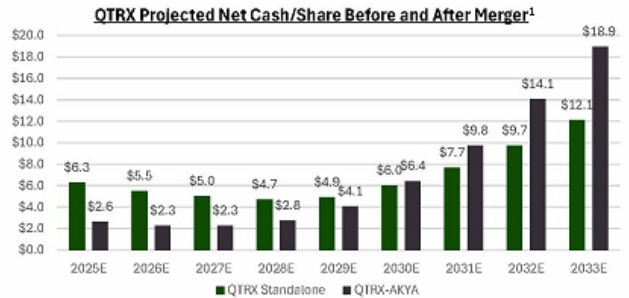
Quanterix's Balance Sheet Is **STRONGER WITHOUT** Akoya

Quanterix would have **>2x the cash/share at year end 2025 in the QTRX stand-alone scenario** than if the merger goes through per mgmt. S-4 projections

- Quanterix currently has ~\$270M in cash (at 12/31/24, net of \$20M related to the Emission acquisition)
 - 2024 cash burn was ~\$40M, implying a **QTRX standalone cash runway of ~6.5 years**
- With \$175M expected cash outstanding following the merger, a combined TTM QTRX-AKYA cash burn of \$91M means that, even after \$20M in optimistic synergies, **the ~\$70M burn rate shrinks cash runway to just 2.5 years**

Fiscal Year	QTRX			AKYA		
	Mgmt. Projections (S-4)	Actual / Consensus ²	Variance %	Mgmt. Projections (S-4)	Actual / Consensus ²	Variance %
FY22A		\$105.5		\$74.9		
FY23A		\$122.4		\$96.6		
FY24E		\$137.2		\$81.8		
FY25E	\$155.0	\$154.7	0%	\$87.0	\$91.6	-5%
FY26E	\$196.0	\$179.9	9%	\$99.0	\$105.3	-6%
FY27E	\$273.0	\$215.2	27%	\$128.0	\$110.9	15%
FY28E	\$450.0	\$290.9	79%	\$192.0	\$120.5	59%
FY29E	\$689.0	\$287.5	140%	\$273.0	\$162.7	68%
FY30E	\$1,022.0	\$327.8	212%	\$370.0	\$182.2	103%

Note that management's projections for the inflection years 2028 through 2030 are significantly above current self-side estimates for both companies. While we are extremely bullish on QTRX's potential, we caution reliance on overly optimistic projections to justify this merger



Source: Capital IQ, and public filings

1) Preliminary 2024 cash balance per 1/13/25 press release. Projected cash represents 2024 cash and securities plus QTRX management's unlevered FCF projections from the S-4. FY25 ending cash/share assumes ~1/3 of FY25 cash usage occurs in 2H, with the starting point the guided \$175M net cash post transaction close. Per share assumes ~1% annual FDSO growth

2) Capital IQ sell-side consensus dated 2/26/25. QTRX FY24E revenue reflects preliminary results from the 1/13/25 press release.

QTRX Shares **GO HIGHER** If Merger Is Voted Down

QTRX price is higher as a stand-alone company compared to QTRX-AKYA at any equal multiple:

If QTRX traded at 1x 2024 revenue there is **33% UPSIDE** for shares, but if the combined company traded at 1x revenue there is **5% DOWNSIDE** in the shares

QTRX shares deserve a higher revenue multiple in a stand-alone scenario than in a combined company scenario due to **higher organic revenue growth (12% organic for QTRX in 2024, versus 0% for the combined company)**, and because **stand-alone QTRX has >2x the cash per share** than the merged company

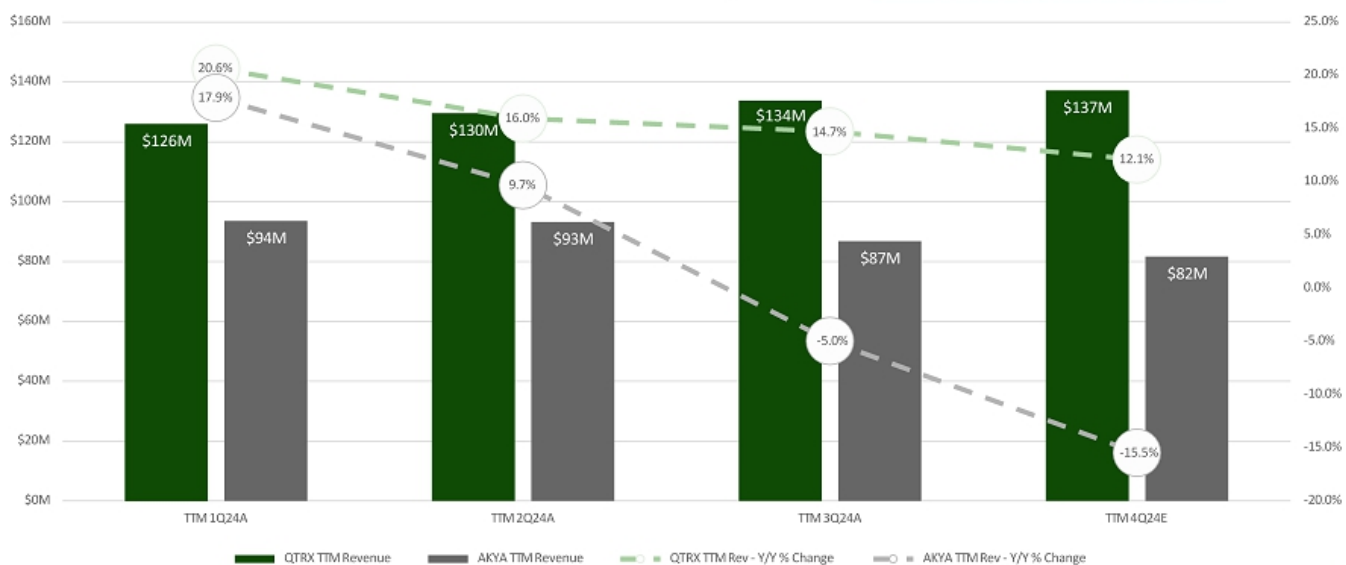
QTRX STANDALONE		
Sensitivity Analysis		
TEV / FY24 Revenue	Share Price	% Upside (Downside)
2.0x	\$13.30	79.7%
1.5x	\$11.57	56.4%
1.0x	\$9.85	33.1%
0.5x	\$8.12	9.7%
0.0x	\$6.39	-13.6%
Capitalization		
QTRX Share Price (3/7/25)		\$7.40
Enterprise Value		\$40.0
FY24 Revenue		\$137.2
Current TEV / FY24 Revenue		0.3x
KLPR Target Multiple		1.5x
KLPR Target Price		\$11.57
% Upside (Downside)		56.4%
<small>KLPR Price Target assumes a combined \$15M in incremental YTD 2025 cash burn and breakup fee, and multiple re-rating in-line with peer group; PT also approximates QTRX price pre-merger announcement. In the stand-alone scenario, QTRX deserves a higher multiple due to higher organic growth rate (12% versus 0% in 2024) as well as significantly more cash per share.</small>		

QTRX-AKYA COMBINED		
Sensitivity Analysis		
TEV / FY24 Revenue	Share Price	% Upside (Downside)
2.0x	\$10.93	47.6%
1.5x	\$8.97	21.3%
1.0x	\$7.02	-5.1%
0.5x	\$5.25	-29.1%
0.0x	\$3.12	-57.9%
Capitalization		
QTRX Share Price (3/7/25)		\$7.40
Enterprise Value		\$253.9
FY24 Revenue		\$219.0
Current TEV / FY24 Revenue		1.2x
KLPR Target Multiple		0.5x
KLPR Target Price		\$5.25
% Upside (Downside)		-29.1%
<small>KLPR Price Target assumes \$175M post-close net cash balance and assigns a 0.5x revenue multiple, which is the average of QTRX stand-alone and Standard BioTools (LAB), given the combined QTRX AKYA has slower proforma growth (0% versus 12% for QTRX stand alone in 2024) and significantly less cash per share.</small>		

Source: Capital IQ, public filings, and QTRX S-4 Prospectus filed 02/13/25

Quanterix Is Clearly the Superior Business

QTRX's growth has shown resilience while AKYA's growth has gone meaningfully negative

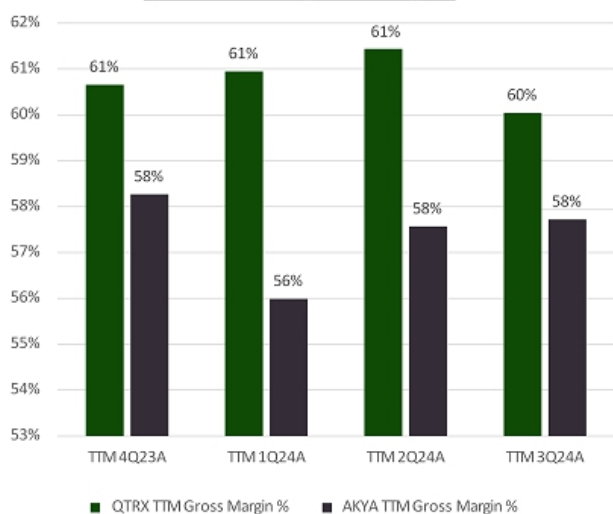


Source: Capital IQ, and public filings

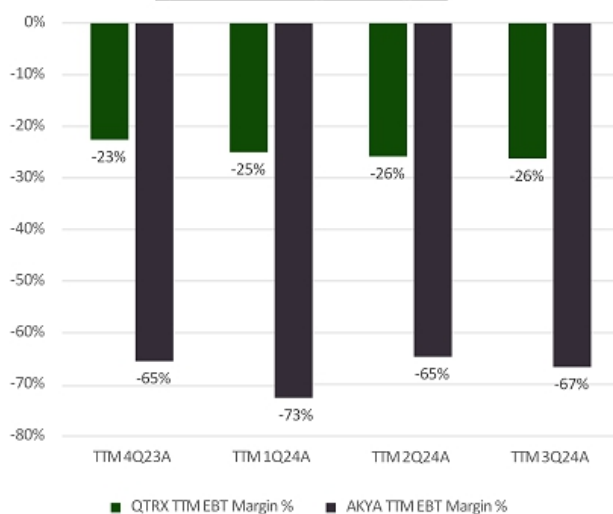
1) QTRX TTM 4Q24E revenue reflects preliminary revenue as reported in the 1/13/25 press release. AKYA 4Q24E reflects Capital IQ consensus estimates as of 2/26/25

Quanterix's Margins Are Superior Compared to Akoya's

QTRX-AKYA Trailing Gross Margin %



QTRX-AKYA Trailing EBT Margin %¹



Source: Capital IQ, and public filings

1) EBT represents earnings before interest and taxes, calculated as GAAP net income plus reported provision for income tax.

Negotiation and Process



Red Flag: Quanterix Ended Up Bidding Against Itself

AKYA contacted 20 parties: 13 signed NDAs, and 3 submitted initial offers – yet QTRX was the only final bidder

- At least **four** bidders walked away due to concerns around AKYA's **profitability, cash burn, and leverage**¹
- Motivated **strategic** buyers were unimpressed by AKYA's CDx pipeline, leading us to question whether this is the high-quality asset the QTRX board purports
- If no serious buyer would touch Akoya over these issues, **why are QTRX shareholders being left holding the bag?**



Source: Capital IQ, and public filings

1) See commentary from Parties E, L, B, and C disclosed within S-4 Prospectus and included on following slides

At Least Two Parties Dropped Explicitly Over Cash Burn

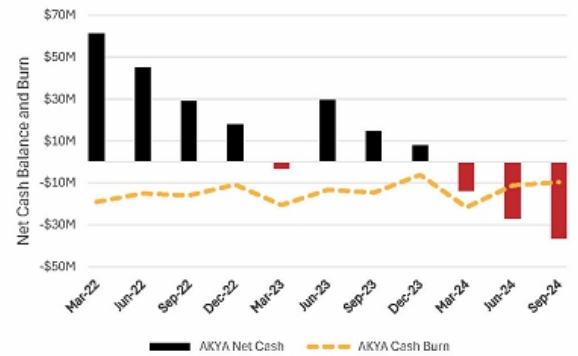
“ Party E indicated that, while they admired the business and product suite that Akoya had developed, **the scale remained small, profitability was uncertain** and Party E believed **Akoya’s financial plan presented significant risks and uncertainties.**

(October 24, 2024) ”

“ Party C advised representatives of PWP that it would **no longer be participating in the process** due to **the extent of Akoya’s projected financial losses and cash burn** and the **dilutive impact** that would represent for Party C. Party C further informed representatives of PWP that Party C had evaluated several strategies to mitigate the dilutive impact of a transaction but was **unable to become comfortable with acquiring Akoya at that time.**

(December 11, 2024) ”

AKYA Cash Burn History
AKYA unrestricted cash and marketable securities net of long-term debt, and CFFO - CX over time



Note that at 9/30/24, AKYA only had \$39.3M in unrestricted cash and marketable securities... meaning **AKYA’s current cash runway is <1 YEAR**

Source: Capital IQ, public filings, and QTRX S-4 Prospectus filed 02/13/25

Two Parties Dropped Due to Akoya’s CDX and R&D Pipeline

- Two bidders – “Party B” and “Party L” – exited after diligence revealed **serious concerns** about Akoya’s CDx path and uncertain R&D pipeline
- Yet QTRX management want to invest **even more** in these projects?

Masoud Toloue
President, CEO & Director

“Akoya has a great companion diagnostics business, and we intend to make sure that, that is moving forward, if not accelerating with investment.

(January 10, 2025) ”

- If major strategic buyers were **unwilling** to bet on Akoya’s uncertain R&D spend, **why should Quanterix shareholders shoulder the burden?**

“Party L indicated that, **after detailed diligence**, Party L’s **conviction around the depth of Akoya’s research and development pipeline** as well as the **time, cost and effort** associated with Akoya’s Companion Diagnostics strategy and cultural compatibility between the companies **had weakened**.

(December 5, 2024) ”

“Party B informed representatives of PWP that it would no longer be participating in the process due to (i) their view that **Akoya’s Companion Diagnostics strategy would take longer to develop and require more investment**, (ii) their perspective on the timing for an overall recovery in the Spatial Biology Instrumentation market in which Akoya operates and (iii) **the amount of balance sheet cash that would need to be allocated to repayment of Akoya debt under a potential transaction**.

(October 31, 2024) ”

Source: Capital IQ, public filings, and QTRX S-4 Prospectus filed 02/13/25

The Merger Is a Bailout for Akoya at QTRX Shareholders' Expense

The S-4 discloses that AKYA's management and advisor knew they would have had to raise dilutive financing to remain solvent

- Absent the merger, **Akoya would have had to raise \$75M** to repay their Nov 2027 MidCap Financial debt
- At 2/10/25, AKYA was trading at ~\$120M market capitalization, or ~\$2.34/share. This means that to raise the \$75M at a 20% discount, AKYA would have had to issue ~**42M new shares**, resulting in **>45% shareholder dilution¹**
- Rather than let AKYA face their own problems, the Quanterix Board is **shifting the burden onto QTRX investors**
- The proposed merger introduces **substantial risk** to Quanterix's business model

“ At the direction of management of Akoya, PWP assumed that, **in order to remain a standalone company, Akoya would need to raise \$75 million in additional capital by issuing Akoya Common Stock at a 20% discount** to the Akoya Current Share Price with a 6% issuance fee, \$25 million the proceeds of which would be used to pay down Akoya's debt at the time of issuance

(QTRX S-4 Prospectus, pg. 165, filed 02/13/25) ”

“ Akoya's management anticipates the Combined Company will experience some **revenue dis-synergies as a result of the Merger**

(QTRX S-4 Prospectus, pg. 153) ”

“ The Combined Company may face **liquidity challenges** during the next few years in light of **significant contingent liabilities and financial obligations and commitments**

(QTRX S-4 Prospectus, pg. 154) ”

Source: Capital IQ, public filings, and QTRX S-4 Prospectus filed 02/13/25

1) Calculated as \$75M plus a 6% issuance fee, divided by the AKYA 2/10/25 share price of \$2.34, discounted by 20%; added to total AKYA common shares outstanding at 1/1/25 of 49.6M per the S4.

RED FLAG: AKYA's Exchange Ratio *Improved* **AFTER** AKYA Missed Q3 and **Shares Fell Over 30%?**

AKYA's pro forma ownership increased from 26% to 30% in the final offer... despite AKYA's Q3 revenue miss and guidance cut, which sent AKYA shares spiraling down over 30%!

- According to the S-4, QTRX's initial November offer gave AKYA shareholders a pro-forma ownership split of 26% (approx. 0.2710x implied exchange ratio)
- On 11/14/24, AKYA reported 3Q24 revenue of \$18.8M (missing FactSet consensus \$25.4M) and cutting FY24 topline guidance by ~17%
- Despite this, on 12/11 – after AKYA's stock fell by **over 20%** – QTRX offered an updated exchange ratio of 0.2970x, or **28% pro forma AKYA ownership**
- Akoya's fundamentals materially worsened throughout the deal process, yet Quanterix settled on an exchange ratio (0.318x) that concedes **>4% of the combined company** to AKYA shareholders
- How did AKYA's worse performance and outlook translate into a higher deal premium? We view this as a sign of poor deal-making by the Quanterix Board and management team, which prioritized an expedited process over shareholder value

Between November 1, 2024, and November 4, 2024, Akoya received first-round proposals from Quanterix, Party A and Party L. Quanterix proposed an all-stock transaction valued at a per share purchase price of \$3.50, which would be translated into a fixed exchange ratio of Quanterix shares for Akoya shares in advance of signing definitive documentation. Based on Quanterix's and Akoya's stock prices at market close on October 31, 2024 the proposal represented a premium of 24% to Akoya's closing stock price and would result in a **pro-forma ownership split of approximately 26% for Akoya stockholders and 74% for Quanterix stockholders on a fully diluted basis**. Quanterix's proposal also anticipated that Quanterix would fully repay Akoya's existing outstanding debt at the closing of a transaction.

On December 11, 2024, Quanterix delivered an updated indication of interest to Akoya (the "December 11 Quanterix Indication of Interest"). The December 11 Quanterix Indication of Interest proposed to acquire 100% of the outstanding shares of capital stock of Akoya and the purchase price would be paid in shares of Quanterix Common Stock based on a 0.2970 exchange ratio, **implying an offer value of \$3.50 per Akoya share**, representing a 49% premium to Akoya's share price based on Akoya's and Quanterix's respective stock closing prices on December 10, 2024 and a **28% pro forma ownership for Akoya stockholders in the combined company**.

Source: Capital IQ, public filings, and QTRX S-4 Prospectus filed 02/13/25

Red Flag: Rescue Loan From QTRX to AKYA

The proposal to offer \$30M bridge financing before the shareholder vote is highly unusual

- Per the S-4, Quanterix must provide Akoya with up to \$30 million in “bridge financing” even before our shareholders approve the merger
- This forces Quanterix to shoulder Akoya’s immediate liquidity crisis, tying up \$30M of our cash in a high-risk convertible note – without any clear shareholder mandate
- If this deal is not consummated, and the convert is issued, QTRX shareholders will be left holding Akoya’s subordinated distressed debt
- Providing backdoor financing before we even vote on the merger is an unacceptable and egregious breach of fiduciary duty
- How did Akoya’s desperate need for this bridge loan factor into Quanterix Board’s valuation of Akoya?
- If Akoya’s really needs this rescue loan, then that calls into question all forecasts utilized by QTRX management to justify this merger

Bridge Financing

Akoya and Quanterix are required to use their respective reasonable best efforts to cooperate in good faith to enter into one or more agreements or instruments pursuant to which Quanterix will provide Akoya with bridge financing in an aggregate principal amount not to exceed \$30,000,000, subject to Akoya having obtained any required consents and satisfied any other conditions under the Akoya Existing Loan Documents, which bridge financing may be drawn by the Company at any time after March 15, 2025 and prior to the Closing. Any such financing would be in the form of subordinated convertible note(s) and would be subject to such other terms agreed between Akoya and Quanterix and in any event as consented to by the lenders or the agent under the Akoya Existing Loan Documents.

Source: Capital IQ, public filings, and QTRX S-4 Prospectus filed 02/13/25

Quanterix Restatement Tainted Merger Negotiations

- QTRX traded at \$15 prior to 11/12/24 announcement that accounting errors necessitated restating financials
- QTRX shares traded from \$15 to \$11 within a week of this disclosure likely due to investor concern over the restatement
- Much of the negotiation for AKYA occurred while QTRX shares were depressed during the restatement process which likely negatively impacted the exchange ratio for QTRX holders

Material weakness over ICFR has been pervasive since 2022

the Treadway Commission (2013 Framework). Based on our evaluation under the framework set forth in Internal Control-Integrated Framework, our management concluded that our internal control over financial reporting was not effective at the reasonable assurance level as of December 31, 2022, due to the material weaknesses described below.

- FY2022 10K (March 2023)

the Treadway Commission (2013 Framework). Based on our evaluation under the framework set forth in Internal Control-Integrated Framework, as well as the actions outlined above, while our management concluded that the Financial Statement Close Process MW and Compensation MW were remediated as of December 31, 2023, management concluded that our internal control over financial reporting was not effective at the reasonable assurance level as of December 31, 2023 due to the material weaknesses described below.

- FY2023 10K (February 2024)

Based on our updated evaluation of the effectiveness of internal control over financial reporting under the Internal Control-Integrated Framework issued by the Committee of Sponsoring Organizations of the Treadway Commission, and in light of the material weaknesses discussed above, our management continued to conclude that our internal control over financial reporting was not effective at the reasonable assurance level as of September 30, 2024.

- 3Q24 10Q (December 2024)

Material weakness led to the restatement

Restatement Background

As previously disclosed by the Company in its Current Report on Form 8-K filed with the SEC on November 12, 2024, in connection with its efforts to remediate a material weakness in its internal control over financial reporting relating to the operating effectiveness of internal controls associated with the accounting for inventory valuation, and while performing closing procedures for the third quarter of 2024, management identified an error related to the capitalization of labor and overhead costs in the Company's inventory balances in prior periods (the "Misstatement"). The error was not caused by any override of controls, misconduct, or fraud. The correction of the Misstatement impacts the previously reported amounts of inventory, cost of product revenue, net loss per common share, and all related financial statement subtotals and totals. In addition to correcting the Misstatement, the Amended Filings also reflect adjustments to correct unrelated errors to other financial statement line items identified by the Company in prior periods which include, but are not limited to, adjustments to property and equipment, accrued compensation and benefits, and operating expenses.

Source: Capital IQ, public filings

A Tainted Process? Half the Transaction Committee Walked Away

Back-to-back director resignations lead us to question deal process further

“ In October of 2023, Akoya and Quanterix had discussions in connection with a potential strategic commercial relationship. From time to time thereafter, Mr. McKelligon and Masoud Toloue, Ph.D., the Chief Executive Officer of Quanterix, **held a number of discussions** ... to explore various ways in which they could collaborate

(QTRX S-4 Prospectus, pg. 132) ”

“ As part of its ongoing evaluation, from time to time, the Quanterix Board has considered various potential strategic transactions and, on October 17, 2023, the Quanterix Board approved the formation of a Transaction Committee (the “Quanterix Transaction Committee”), then comprised of **Brian Blaser, Karen Flynn, Paul Meister and Laurie Olson**, to evaluate certain strategic alternatives

(QTRX S-4 Prospectus, pg. 132) ”

“ On October 2, 2024, the Quanterix Board met and resolved to reconvene the previously formed Quanterix Transaction Committee for the purpose of facilitating the consideration and review of a potential strategic transaction with Akoya. **In light of the departure of Mr. Blaser and Ms. Olson** from the Quanterix Board earlier in 2024, **Martin Madaus** was appointed to the Quanterix Transaction Committee, which was then comprised of **Ms. Flynn, Mr. Meister and Mr. Madaus**

(QTRX S-4 Prospectus, pg. 134) ”

Source: Capital IQ, public filings, and QTRX S-4 Prospectus filed 02/13/25

Conflicts of Interest & Governance



Significant Undisclosed **Conflict of Interest**

Dr. Ivana Magovčević-Liebisch appointed to Quanterix Board on October 2, 2024 – **the same day merger discussions resumed**

- **DUAL-FIDUCIARY CONFLICTS:** Dr. Magovčević-Liebisch serves **simultaneously** as a **director at both Quanterix and Acrivon Therapeutics (ACRV)**
 - ACRV has an **INDIRECT VESTED INTEREST** in the QTRX-AKYA deal: **ACRV's flagship OncoSignature diagnostic test critically relies on Akoya's technology platform**
 - Dr. Magovčević-Liebisch personally owns stock options for ~32,500 ACRV shares (plus potentially another 20,275 granted in 2025); **total potential position value >\$1 million based on median broker ACRV price targets**
- Appointment timeline raises serious independence concerns given ACRV director role and **LACK OF RECUSAL DISCLOSURES**, as QTRX incorporated assumptions around the AKYA-ACRV partnership in financial projections (S-4, pg. 170)
 - October 2, 2024: Dr. Magovčević-Liebisch appointed to the QTRX Board; **immediately attends transaction committee discussions regarding AKYA merger**
 - October 18, 2024: ACRV and AKYA amend commercial partnership, **adding up to \$3M in new pre-commercial milestones** paid to AKYA (**~6% AKYA TTM cash burn**)¹
 - October 25, 2024: QTRX submits initial non-binding offer to acquire AKYA

Given: 1) Transaction Committee reconvened same day as Board appointment; 2) Shared directorship with ACRV; and 3) AKYA and ACRV have a VESTED-INTEREST in each other through their CDX Agreement; ACRV has an INDIRECT VESTED-INTEREST in the QTRX-AKYA deal; NON-RECUSAL introduces DUAL-FIDUCIARY CONFLICTS

Source: Capital IQ, public filings, and QTRX S-4 Prospectus filed 02/13/25

1) \$3M divided by the sum of AKYA 9/30/24 TTM cash flow from operations (-\$46.7M) and CapEx (-\$2.0M)

Why Did Dr. Madaus Remain Board Chair Despite <50% Votes Cast?

Failure to obtain majority voting should have prompted immediate resignation

- Election votes from the June 3, 2024, annual meeting show a significant number of QTRX shareholders **withheld** votes from Karen Flynn and Martin Madaus – in fact, **Dr. Madaus received only 49.9% of total votes cast**
- As a refresher, ShareholderEducation.com (Broadridge)¹ defines a “withheld” vote as:
 - *Similar to abstain, but is used in instances in which nominees run unopposed and therefore only need a single vote to earn a “plurality.” In these cases, “against” votes are meaningless, so you only have the option to vote “for” or “withhold.”*
- In 2019, nearly 90% of S&P 500 Index companies implemented resignation policies for directors who fail to receive majority voting support²
- Quanterix’s own voting results clearly signal **no confidence** in Dr. Madaus, and **undermines trust** in the current Board
- The Board should have **immediately demanded** Dr. Madaus’ resignation – instead, he continues to serve as a member of the Quanterix Transaction Committee

On October 2, 2024, the Quanterix Board met and resolved to reconvene the previously formed Quanterix Transaction Committee for the purpose of facilitating the consideration and review of a potential strategic transaction with Akoya. In light of the departure of Mr. Blaser and Ms. Olson from the Quanterix Board earlier in 2024, Martin Madaus was appointed to the Quanterix Transaction Committee, which was then comprised of **Ms. Flynn**, Mr. Meister and **Mr. Madaus**.

QTRX Shareholders Overwhelmingly Disapprove BoD Appointments

	For	Withheld	Broker Non-Votes
Karen A. Flynn	16,037,402	10,449,328	5,106,768
Martin D. Madaus, Ph.D.	13,208,474	13,278,256	5,106,768

Source: QTRX BK dated 06/07/2024 and S-4 Prospectus filed 02/13/25

1) [Proxy Voting – Frequently Asked Questions | Broadridge](#)

2) [Glass Lewis 2025 US Benchmark Policy Guideline](#)

[22]

Madaus Governance: A Willful Disregard of Shareholder Feedback

Dr. Madaus' direct involvement on the Transaction Committee AFTER Advisor and shareholder opposition further muddies the independence of the proposed AKYA deal

- The Quanterix Board **double-downed on bad governance practices** by allowing this non-independent Chair to join the Transaction Committee overseeing QTRX's **value-destructive** pursuit of AKYA
- In its 2024 Proxy Analysis, ISS explicitly classifies Dr. Madaus as a "non-independent" director, **raising significant independence and oversight concerns**
- Rather than resign after the majority WITHHOLD vote at the 2024 Annual Meeting, Dr. Madaus **remained** Chairman of the Board – **a willful disregard of shareholder feedback**



Source: ISS Proxy Analysis & Benchmark Policy Voting Recommendations – Quanterix 2024 Annual Meeting (May 10, 2024)

Strategic Rationale



Rebutting QTRX's Strategic Rationale From Their 3/3/25 PR

Scale and Profitability:

- QTRX management's ~\$1B revenue target (**~5x growth**) within five years post-merger is **highly unrealistic given 0% pro forma combined 2024 revenue**

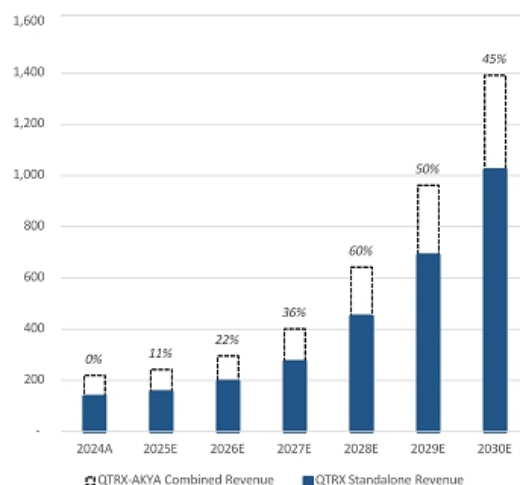
Synergy Generation:

- AKYA management explicitly anticipate revenue **dis-synergies** post-merger (S-4, pg. 153), making us skeptical of the projected \$40M annual cost synergies by 2026 – particularly in light of AKYA's recent ~35% workforce reduction

Expanded Addressable Market:

- Quanterix has promoted its "best-in-class" liquid biopsy neurology tests, but these are not yet commercially proven. **Management should first demonstrate successful execution in clinical diagnostics before pursuing dilutive high risk acquisitions**
- Quanterix is already expanding into Immunology/Oncology markets through its upcoming SIMOA ONE platform launch, making the Akoya acquisition **unnecessary**

QTRX Management's S-4 Projections¹



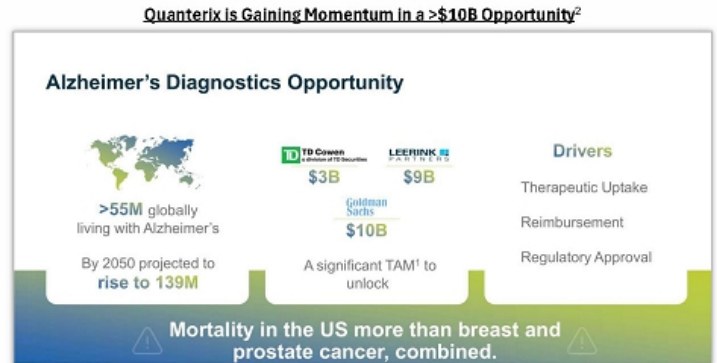
Source: ISS Proxy Analysis & Benchmark Policy Voting Recommendations – Quanterix 2024 Annual Meeting (May 10, 2024)

1) Pro forma combined 2024 revenue is the sum of QTRX pre-announced \$137.2M and AKYA Capital IQ consensus \$81.8M

Merger Distracts From Core Alzheimer's Diagnostics Opportunity

Quanterix is at a critical juncture in Alzheimer's diagnosis with its groundbreaking blood test, but the merger risks diverting cash and management attention away from this high-value opportunity

- QTRX's ultra-sensitive "Simoa" technology enables detection of early-stage, pre-symptomatic disease markers of neurodegenerative diseases, including Alzheimer's, with a simple blood draw
 - This approach has the potential to reach a broad patient population and could **dramatically** lower diagnostic costs relative to PET imagery or invasive procedures
- Alzheimer's currently affects >55M people globally and is expected to reach 139M by 2050; street analysts estimate a ~\$10B TAM opportunity¹
 - QTRX management have emphasized the need to rapidly advance its Alzheimer's diagnostic capabilities through R&D, FDA submissions, and partnerships.
- Successfully integrating Akoya's spatial imaging platform will demand significant time and attention, pulling management focus away from Quanterix's high-priority pipeline



1) [Dementia statistics | Alzheimer's Disease International \(ADI\)](#)

2) [QTRX June 2024 Investor Deck](#)

Dr. Toloue Alleges NIH Cuts Are Cause of QTRX Shares Decline

While partially true, Kent Lake believes the AKYA merger is the bigger driver of QTRX share decline.

However, if Dr. Toloue is correct and NIH issues have caused a revaluation across the sector, **that is still an excellent reason for shareholders to reject this flawed merger.**

- If QTRX's current multiple accurately reflects a standalone company valuation adjusted for NIH headwinds, **applying the same multiple to AKYA implies a negative \$20M equity value**
- Bottom line: If Dr. Toloue is correct, NIH-driven industry revaluation of QTRX shares **further strengthens the case for shareholders to vote NO on this deal**

AKYA at QTRX's Current Multiple	
AKYA Pro Forma Share Price	-\$0.41
Shares Outstanding	51.4
Market Cap	-\$21.0
Net Debt	\$47.0
Enterprise Value	\$26.0
FY24 Revenue	\$81.8
QTRX TEV / FY24 Rev	0.3x

“ *Macroeconomic pressures are expected to persist, extending sales cycles, particularly for higher CapEx life science... the recent announcements regarding NIH funding... have increased the uncertainty in the academic environment... In the near term, based on our initial conversations with customers, recently announced federal funding freezes... have added significant uncertainty* ”

– Pacific Biosciences of California, 4Q24 CC (February 13, 2025)

“ *Our team has done careful analysis and research, which we have factored into our guidance, modeling a mid-teens percentage decline in Americas academic revenue... The situation remains fluid, evolving day by day, but we anticipate that a reduction in NIH funding is likely to impact overall academia spend and priorities, particularly delaying capital equipment purchases.* ”

– Standard BioTools, 4Q24 CC (February 26, 2025)

Source: Capital IQ, public filings, and QTRX S-4 Prospectus filed 02/13/25

NIH Funding Cuts Give Shareholders Even More Reason To Reject This Deal

- A broad reduction in research spending means fewer grants to NIH-funded labs, directly shrinking the pool of instrument buyers, and likely undermining both Akoya and Quanterix's Outlook. This **intensifies** the risk for Akoya, whose core revenue is ~1/3 instrument sales (vs. QTRX at ~8%)²
- AKYA's **high cash burn and substantial leverage** leave it more exposed to a softer capital-equipment environment – **amplifying concerns over financial stability**
- Industry deterioration could jeopardize management's near-term synergy assumptions, **making the combined model look even worse**
- Overall, we believe the subsequent impact of NIH funding cuts make AKYA's balance sheet situation untenable – currently, **AKYA likely has little (or no) positive equity value**

“ NIH overhang and anxiety among researchers **remain at peak with no end in sight**. Researchers are “hoping” for overhang to resolve but **are not in control**. **Expect significant impact (layoffs/shutdowns) if NIH IDC cap of 15% holds, but for now it appears business as usual for 1Q with risk in 2Q-3Q**. In our view, **research labs and KOLs think they are walking towards a cliff but are hoping a bridge will eventually emerge as it would be too catastrophic to lose American leadership in biomedical research and innovation**. ”

– Leerink Partners Research Report (February 26, 2025)

“ A substantial portion of our revenue in the United States is derived from academic and governmental institutions, research organizations and other entities that may rely in part on academic and government funding, **including grants from the U.S. National Institutes of Health (NIH) and other U.S. government agencies**... If U.S. academic and governmental researchers experience reductions or delays in government funding, or modifications of the terms or conditions of funding, including allowable overhead rates, **they may reduce or delay their purchases of our products and services, which could have a material effect on our revenues and financial performance**. ”

– BRKR FY24 10K (Filed March 3, 2025)

Source: Capital IQ, public filings

- 1) Per QTRX CFO on 1/10/25 M&A Call: "...our split is ... between 45% and 55% between pharma biotech and academia. For Akoya, the split is slightly more skewed towards pharma a little over 60% for pharma, biotech and CROs and the rest is academic and government."
- 2) Instrument revenues as a percentage of total revenues over the trailing twelve-months 9/30/24.

AKYA's Competitive Position Is Weaker Than Management Claims

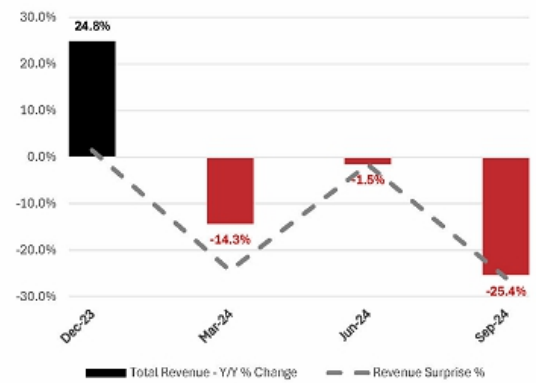
Commentary by Akoya management reveal that **market share is already slipping** – So again we ask, why do the QTRX Board want to invest more dollars in an already struggling asset?

“ And as you look at the areas where we've seen the most challenges in terms of the highest volatility, it's really been primarily in North America and really within the academic environment... I wouldn't say it's necessarily competition. **I think the macro is the overwhelming driver. That said, we do see additional competition out there from our colleagues at Lunaphore...**

So in areas like core labs, CROs, people that are providing shared services, **you are competing for dollars versus competing head-to-head on should I buy a spatial proteomic platform from Akoya or somebody else...** For us, one telltale sign is as we look at the contribution of core labs to our revenue, **that's been an area of meaningful weakness this year...** In terms of a head-to-head competitive dynamic, I would say **where we are potentially losing in the discovery side with the PhenoCycler-Fusion is where we don't have content... That is an area where the competition is probably the most fierce.**

– Brian McKelligon, AKYA President, CEO & Director (November 14, 2024) ”

AKYA – Revenue Changes and Surprise %



Source: Capital IQ, public filings

New Competitor Lunaphore (TECH) Is Taking Market Share From Akoya

AKYA's three consecutive misses and guidance cuts in 2024 coincided with Bio-Techne launching Lunaphore's competing COMET system for the first time

Competitor Quotes

“ We have a fantastic platform in COMET through the Lunaphore application, obviously, newer in the market, **which is now 3 quarters or so**. We see good traction. Initially, we couldn't keep up with demand. We fixed that. It's a very competitive instrument and promising for us.

– TECH @ Stephens Conference (November 2024) ”

“ We anticipate improvement in the Diagnostics and Spatial Biology operating margin **as Lunaphore continues to scale ... A year ago, we launched the COMET**, and we have been very happy with the uptake and the capabilities of an instrument ... we expect the pull-through of consumables on the COMET to be one of the higher pull-through instruments in our portfolio.

– TECH Q2'25 CC (January 2025) ”



Merger Deepens Quanterix's Exposure to China During Trade War

AKYA's China sales – hurt by slow demand and tightened regulations – now become QTRX's burden

- AKYA derives ~10-11% of total revenue¹ from China, while QTRX has relatively small² end-market exposure
- AKYA's relies upon its global distributor channel for Chinese sell-through, and any further international trade disputes (tariffs, export controls, etc.) could further its supply chain
- Illumina – the 800lb gorilla – is facing its own issues. On March 4, 2025, China imposed new bans on the import of genetic sequencers in response to tariffs by the Trump administration
- We caution that ILMN's issues could act as a harbinger for broader industry disruptions

The imposition of tariffs and trade restrictions as a result of international trade disputes or changes in trade policies may adversely affect our sales and profitability... Most notably, on February 1, 2025, President Donald Trump signed executive orders imposing a 25% tariff on certain imports from Mexico and Canada... and a 10% tariff on certain imports from China... These tariffs, and the related geopolitical uncertainty between the United States and China, may cause decreased demand for our products, which could have a material adverse effect on our business and results of operations.

– CTKB FY24 10K (Filed February 27, 2025)

Source: Capital IQ, public filings

- Per AKYA CEO on 1Q24 CC, China comprises between 60-70% of APAC revenue (~16% mix over the TTM 3Q24)
- Per QTRX CEO on 2Q24 CC
- [Illumina Faces New Headaches After Trump's Trade War, Cuts to Science Funding - WSI](#)

China bans imports of Illumina's gene sequencers right after Trump tariff action

By Andrew Silver

March 4, 2025 2:46 AM PST Updated 5 hours ago



Illumina HiSeq X Ten gene sequencing system is shown in this undated handout photo distributed by Reuters, September 26, 2022. (Reuters/Phil Witte via AP Photo/Justin Sullivan)

Summary	Companies
<ul style="list-style-type: none"> The export ban takes effect on March 4. Illumina says it is assessing China's announcement. Shares in ILMN's China rivals soar. 	

SHANGHAI, March 4 (Reuters) – China announced a ban on Tuesday on imports of genetic sequencers from U.S. medical equipment maker Illumina (ILMN.O), just minutes after U.S. President Donald Trump's government imposed 10% tariffs on Chinese goods.

The ban is part of a series of measures China announced in response to the extra U.S. tariffs and comes after Beijing put Illumina on its "unreliable entities" list in January.

China accounts for about 7% of Illumina's sales. Gene sequencers help determine the sequence of DNA, allowing scientists to study genetic variations associated with diseases and diagnose rare genetic conditions.

Conclusion

It's Not Too Late! Quanterix Isn't Handcuffed to This Deal!

Under the merger terms, if QTRX identifies a better opportunity or AKYA's issues prove too great, Quanterix can terminate the deal (subject to \$9M fee).

Kent Lake Calls on Quanterix's Board to consider what is best for QTRX shareholders and withdraw this deal immediately.

If this deal goes to a vote, QTRX will not pay any break fees if shareholders reject the deal.

Under the Merger Agreement, Quanterix is required to pay Akoya a termination fee of \$9 million upon any of the following events:

- a Quanterix Recommendation Change Termination;
- a Quanterix No-Shop Termination;
- a Quanterix Superior Proposal Termination, or
- if (i) the Merger Agreement is terminated by Akoya by way of a Quanterix Breach Termination or by Quanterix or Akoya by way of an Expiration Termination or a No Quanterix Vote Termination, (ii) at or prior to the Quanterix stockholders' meeting to vote on the approval of the Quanterix Share Issuance Proposal (in the case of a No Quanterix Vote Termination) or at or prior to the time of such termination (in the case of a Quanterix Breach Termination or Expiration Termination), an Acquisition Proposal with respect to Quanterix is publicly proposed, disclosed or known, and such Acquisition Proposal is not irrevocably and publicly withdrawn at or prior to such applicable time described in this clause (ii), and (iii) concurrently with or within 12 months after any such termination, Quanterix or any of its subsidiaries enters into a definitive agreement (which is ultimately consummated, whether during such 12-month period or thereafter) with respect to, or otherwise consummates, any Acquisition Proposal with respect to Quanterix (substituting 50% for 20% in the definition of "Acquisition Proposal").

22

“ **the Quanterix Board is permitted under the Merger Agreement to withdraw its recommendation if required by its fiduciary duties** ”
(QTRX S-4 Prospectus, pg. 154)

“ **Kent Lake is admittedly not an activist investor, and this is their first public campaign. However, that does not stop them from making an incredibly compelling case for abandoning the merger. Their case for abandonment made on February 13 has only been strengthened by the significant contraction in the QTRX share price, now trading below net cash implying a negative EV, and widening the valuation gap between them and AKYA** ”
- 13D Monitor (March 6, 2025)

Source: Capital IQ, public filings, and QTRX S-4 Prospectus filed 02/13/25

[33]

Questions for Quanterix

- How do you justify QTRX shareholders receiving 18% of the combined company's proforma enterprise value for the superior Quanterix business, while AKYA shareholders will receive 82% for a clearly inferior business?
- If uncertainty around NIH funding are the cause of a significant revaluation of Quanterix's stock price, then why isn't it also an opportunity to revalue Akoya's business and end this deal?
- Given the S4 discloses that multiple bidders walked away from AKYA for a variety of issues, including high cash burn and low profitability, and QTRX was bidding against itself by end of process, shouldn't that change the QTRX board's recommendation to go forward with this merger?
- How is the Transaction not a bailout for Akoya, given that the latter has disclosed a going concern risk and was facing a major cash crunch and on the cusp of needing to do a large dilutive financing for an amount nearly equal to their entire market capitalization?
- What is management's rationale for providing Akoya with \$30 million in bridge financing without any shareholder input?
- Given Akoya's 15% revenue decline in 2024, how does Quanterix know that is not due to changes Akoya's competitive position?
- Why does Quanterix continue to pursue a transaction that, since announced, has contributed to a significant decline in the QTRX share price?

Source: QTRX S-4 Prospectus filed 02/13/25

About Kent Lake PR LLC & Kent Lake Partners LP

Kent Lake Partners LP is an investment fund founded by Ben Natter in 2019 with a focus on U.S. small/mid cap public equities.



KENT LAKE PARTNERS

- Kent Lake has been investing in and following Quanterix since 2022, and currently is a top three QTRX shareholder owning over 7% of QTRX's common stock
- Historically, Kent Lake has avoided shareholder activism; this marks the Fund's **first** public stand against a decision by one of its portfolio companies
- Healthcare is the majority of Kent Lake Partners LP's investment exposure
- Ben Natter has over a decade of experience as a successful healthcare investor
- Acting in our investors' best interests, we could not let the proposed deal move forward without raising our concerns

1) Based on Kent Lake Partners ownership refer to latest form 13D/A on file with the SEC

Contact Information

Investor Contacts

Ben Natter, 415-237-0007

info@kentlakecap.com

Saratoga Proxy Consulting LLC

John Ferguson / Ann Marie Mellone

(212) 257-1311 / (888) 368-0379

info@saratogaproxy.com

Media Contact

Longacre Square Partners

KentLake@longacresquarepartners.com

Appendix

Akoya's Nov' 2024 Off-Cycle RSU Grants Introduce Accounting Risk

- **RSUs Awarded Mid-Sales Process:** AKYA management issued sizable RSU awards¹ on November 19, 2024, during active merger discussions
- AKYA's November 14 announcement of "evaluating strategic alternatives" understates material positive MNPI available at the time, and **these equity awards could be considered SPRING-LOADED GRANTS:**
 - AKYA had already received multiple bids within \$3.50 to \$5.00 per-share range
 - AKYA management had already held post-earnings investor discussions when RSUs were granted
 - Bidding process had already entered the 2nd round
- AKYA states the fair value of RSUs are calculated using the grant date stock price
 - This is **INAPPROPRIATE** for spring-loaded RSUs under SAB 120 and would result in a **MISSTATEMENT**
- **Given their own restatements, QTRX MUST NOT IGNORE additional accounting risk that could MISLEAD investors and further affect financial reporting**

Current Price of the Underlying Share (Including Considerations for Spring-Loaded Grants)?

Determining whether an adjustment to the observable market price is necessary, and if so, the magnitude of any adjustment, **requires significant judgment**. The staff acknowledges that companies generally possess non-public information when entering into share-based payment transactions. The staff believes that an observable market price on the grant date is generally a reasonable and supportable estimate of the current price of the underlying share in a share-based payment transaction, for example, when estimating the grant-date fair value of a routine annual grant to employees that is not designed to be spring-loaded.

However, companies should carefully consider whether an adjustment to the observable market price is required, for example, **when share-based payments arrangements are entered into in contemplation of or shortly before a planned release of material non-public information, and such information is expected to result in a material increase in share price**. The staff believes that **non-routine spring-loaded grants merit particular scrutiny** by those charged with compensation and financial reporting governance. Additionally, **when a company has a planned release of material non-public information within a short period of time after the measurement date of a share-based payment, the staff believes a material increase in the market price of the company's shares upon release of such information indicates marketplace participants would have considered an adjustment to the observable market price on the measurement date to determine the current price of the underlying share.**

1) See AKYA Form 4 filings dated November 21, 2024

2) [SEC.gov | Staff Accounting Bulletin No. 120](#)

Deal Offered QTRX Holders Unfair Valuation on Date of Announcement

At QTRX's pre-deal price of \$11.73, AKYA shareholders received an EV/Revenue Multiple TWICE as high as QTRX shareholders, and >50% of the combined pro forma enterprise value, showing it was an unfavorable deal for QTRX shareholders from the start

AKYA IMPLIED VALUE - UNAFFECTED (1/8/25)	QTRX	Implied AKYA
Share Price (1/8/25)	\$ 11.73	0.318x
Total Akoya FDSO		51,448
QTRX Equivalent FDSO	39,745	16,361
Market Capitalization	\$ 466,213	\$ 191,912
Replacement of AKYA Options + Frac. Shares		\$ 13,698
Total Value of Consideration		\$ 205,610
Net Cash	\$ 269,123	\$ (46,993)
Enterprise Value	\$ 197,090	\$ 252,603
% of Pro Forma TEV	44%	56%
<i>Forward Revenue (CIQ Est.)</i>		
2024E Revenue	\$ 137,200	\$ 81,906
2025E Revenue	\$ 157,607	\$ 91,584
2026E Revenue	\$ 189,924	\$ 105,541
TEV / 2024E Revenue	1.4x	3.1x
TEV / 2025E Revenue	1.3x	2.8x
TEV / 2026E Revenue	1.0x	2.4x

Source: Capital IQ, and public filings

QTRX Share Price Below \$30 is a Wealth Transfer to AKYA

For QTRX shareholders to receive an equivalent TEV/Revenue multiple, QTRX shares would need to exceed \$30, compared to the \$7.83 figure in the S-4

AKYA JUSTIFIED VALUE	QTRX	Implied AKYA
Pro Forma Share Price	\$ 30.16	0.318x
Total Akoya FDSO		51,448
QTRX Equivalent FDSO	39,745	16,361
Market Capitalization	\$ 1,198,540	\$ 493,368
Replacement of AKYA Options + Frac. Shares		\$ 13,698
Total Value of Consideration		\$ 507,066
Net Cash	\$ 269,123	\$ (46,993)
Enterprise Value	\$ 929,417	\$ 554,059
% of Pro Forma TEV	63%	37%
<i>Forward Revenue (CIQ Est.)</i>		
2024E Revenue	\$ 137,200	\$ 81,790
2025E Revenue	\$ 154,701	\$ 91,571
2026E Revenue	\$ 179,867	\$ 105,306
TEV / 2024E Revenue	6.8x	6.8x
TEV / 2025E Revenue	6.0x	6.1x
TEV / 2026E Revenue	5.2x	5.3x

Source: Capital IQ, and public filings

New Director Has **Significant Undisclosed Conflict of Interest**

Why was Dr. Magovčević-Liebisch, who sits on the Board of Acrivon Therapeutics (ACRV) – which has a material interest in Akoya's survival – appointed to the Quanterix Board on the same day transaction discussions resumed?

- On October 2, 2024, QTRX appointed Vigil Neuroscience President, CEO, and Director, to serve as a Class III director. Per the 8-K filing, **"there is no arrangement or understanding between Dr. Magovčević-Liebisch and any other person pursuant to which she was selected as a director of the Company"**
- Dr. Magovčević-Liebisch concurrently serves on the Board of Acrivon Therapeutics (ACRV) – a company whose flagship OncoSignature test is critically reliant on Akoya's platform
 - From the ACRV March 28, 2024, 10-K: **"We rely on our external companion diagnostic partner, Akoya, to perform ACR-368 OncoSignature testing in our clinical trial. If Akoya encounters delays or technical challenges, enrollment in our clinical trials may be substantially delayed"**
- The Quanterix Board met to discuss a potential transaction with Akoya **on the same day Dr. Magovčević-Liebisch was appointed** (10/2/24)... **QTRX submitted an offer three weeks later**
- Dr. Magovčević-Liebisch personally owns ACRV stock options to purchase 32,500 shares (with another 20,275 likely granted following the 2025 AGM). While ACRV's stock is only ~\$5.5 (as of February 25, 2025), the median sell-side price target is ~\$20 – **meaning these options could be worth well over \$1 million**
- Given the sudden appointment of Dr. Magovčević-Liebisch, **was the Quanterix Board's decision effectively sealed before the process even began, making any further diligence irrelevant?**

“ Also on October 18, 2024, the Quanterix Transaction Committee held a meeting to, among other things, discuss the process letter and a potential strategic transaction with Akoya. **That meeting was also attended by Ivana Magovčević-Liebisch** (a member of the Quanterix Board) ...

On October 25, 2024, the Quanterix Transaction Committee held a meeting to, among other things, further discuss ... a potential strategic transaction with Akoya ... **That meeting was also attended by William Donnelly and Ivana Magovčević-Liebisch** ... At that meeting, and after discussion, the Quanterix Transaction Committee resolved to recommend to the Quanterix Board that Quanterix submit a non-binding offer to acquire 100% of the Akoya Common Stock in an all-stock transaction.

(QTRX S-4 Prospectus, pg. 135 - 136)

”

Source: Capital IQ, public filings, and QTRX S-4 Prospectus filed 02/13/25

On June 20, 2024, Acrivon Therapeutics, Inc. (the "Company") held its 2024 annual meeting of stockholders (the "Annual Meeting"). At the Annual Meeting, a quorum of 27,422,027 shares of the Company's common stock, or 88.81% of the outstanding shares of common stock entitled to vote as of the record date of April 22, 2024, were present or represented by proxy.

Two items of business were acted upon by the stockholders at the Annual Meeting. The final results for the votes regarding each proposal are set forth below.

Proposal One: Election of Kristina Masson, Ph.D., M.B.A.; Sharon Shacham, Ph.D., M.B.A.; and Ivana Magovčević-Liebisch, Ph.D., J.D. as Class II Directors.

Kristina Masson, Ph.D., M.B.A.; Sharon Shacham, Ph.D., M.B.A.; and Ivana Magovčević-Liebisch, Ph.D., J.D. were each elected to serve as a Class II director to hold office until the Company's 2027 Annual Meeting of Stockholders and until the election and qualification of his successor. Votes were cast as follows:

	For	Withheld	Broker Non-Votes
Kristina Masson, Ph.D., M.B.A.	24,563,469	109,386	2,749,172
Sharon Shacham, Ph.D., M.B.A.	23,268,783	1,466,015	2,687,229
Ivana Magovčević-Liebisch, Ph.D., J.D.	24,498,933	237,318	2,685,776

Source: public filings

Companion Diagnostic Agreement

In June 2022, we entered into a companion diagnostic agreement with Akoya Biosciences, Inc., or Akoya, pursuant to which we agreed to co-develop, validate, and commercialize our proprietary ACR-368 OncoSignature test, the companion diagnostic that will be used to identify patients with cancer most likely to respond to ACR-368.

Pursuant to the agreement, we paid Akoya a one-time, non-refundable, non-creditable upfront payment in the amount of \$0.6 million. We are obligated to pay Akoya up to an aggregate of \$17.3 million upon the achievement of specified development milestones. In October 2024, we entered into the Fourth Amendment to the Akoya Agreement, or the Fourth Amendment, with Akoya, effective September 30, 2024. The Fourth Amendment added pre-commercialization progress milestones totaling \$3.0 million. As of November 13, 2024, development milestones totaling \$14.6 million have been achieved by Akoya under the agreement. Other than certain specified pass-through costs, each party is responsible for its own costs associated with the development of the companion diagnostic. Akoya will procure and manufacture necessary supplies to perform the ACR-368 OncoSignature test to support our clinical development and commercial requirements, in accordance with a supply agreement to be mutually agreed upon by the parties. We may terminate the agreement at our convenience, subject to the payment of a termination fee in the amount of \$1.0 million.

**Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned
(e.g., puts, calls, warrants, options, convertible securities)**

1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	3A. Deemed Execution Date, if any (Month/Day/Year)	4. Transaction Code (Instr. 8)	5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4, and 5)	6. Date Exercisable and Expiration Date (Month/Day/Year)		7. Title and Amount of Underlying Securities (Instr. 3 and 4)		8. Price of Derivative Security (Instr. 5)	9. Number of Derivative Securities Beneficially Owned Following Reported Transaction(s) (Instr. 4)	10. Ownership Form of Derivative Security: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)
						Date Exercisable	Expiration Date	Title	Amount or Number of Shares				
Stock Option (Right to Buy)	\$ 3.54	02/08/2024		A	32,500	(1)	02/07/2034	Common Stock	32,500	\$ 0	32,500	D	

Reporting Owners

Reporting Owner Name / Address	Relationships			
	Director	10% Owner	Officer	Other
Magovcevic-Liebisch Ivana C/O ACRIVON THERAPEUTICS, INC. 480 ARSENAL WAY, SUITE 100 WATERTOWN, MA 02472	X			

Peers Are Issuing New NIH Risk Disclosures

In addition, various state, federal and foreign agencies that provide grants and other funding may be subject to stringent budgetary constraints that could result in spending reductions, reduced grant making, reduced allocations or budget cutbacks, including as a result of negative or worsening conditions in the general economy, **which could jeopardize the ability of these customers, or the customers to whom they provide funding, to purchase our solutions.** For example, congressional appropriations to the National Institutes of Health (the "NIH") may experience occasional year-over-year decreases in appropriations. There is no guarantee that NIH appropriations will not decrease or halt in the future. **A decrease in the amount or halt of, or delay in the approval of, appropriations to NIH or other similar United States or foreign organizations, such as the Medical Research Council in the United Kingdom, could result in fewer grants benefiting life sciences research.** These reductions or delays could also result in a decrease in the aggregate amount of grants awarded for life sciences research or the redirection of existing funding to other projects or priorities, any of which in turn could cause our customers and potential customers to reduce or delay purchases of our solutions.

– CTKB FY24 10K (Filed February 27, 2025)

Recently, budget cuts and layoffs at various federal agencies and programs, including at the National Institute of Health (NIH), have **created uncertainty and budget cuts at various research organizations and institutes that rely on NIH funding.** Our operating results may fluctuate substantially due to any such reductions and delays. Any decrease in our customers' budgets or expenditures, or in the size, scope or frequency of their capital or operating expenditures, could materially and adversely affect our business, results of operations, financial condition and prospects.

– SEER FY24 10K (Filed March 3, 2025)

A decrease in the amount of, or delay in the approval of, appropriations to National Institutes of Health ("NIH") or other similar U.S. or international organizations, such as the Medical Research Council in the United Kingdom, could result in fewer grants benefiting life sciences research. For example, on February 7, 2025, the NIH imposed a standard indirect rate of 15% across all NIH grants for indirect costs, defined as "facilities" and "administration," in lieu of a separately negotiated rate for indirect costs in every grant. **Indirect costs represent more than 25% of total grant dollars in 2023 awarded by the NIH.** This limitation may impact research institution's ability to govern or even accept NIH grants, **thereby affecting total spending on products such as ours.** Further, as of February 2025, **much of the current period NIH funding allotments have either been slowed or frozen.** These reductions or delays could also result in a decrease in the aggregate amount of grants awarded for life sciences research or the redirection of existing funding to other projects or priorities, any of which in turn could cause our potential customers to reduce or delay purchases of our products.

– QSI FY24 10K (Filed March 3, 2025)

A substantial portion of our revenue in the United States is derived from academic and governmental institutions, research organizations and other entities that may rely in part on academic and government funding, including grants from the U.S. National Institutes of Health (NIH) and other U.S. government agencies. Government funding may fluctuate and is subject to annual appropriations and budgetary constraints, and there is no assurance that such funding will continue at current levels. **If U.S. academic and governmental researchers experience reductions or delays in government funding, or modifications of the terms or conditions of funding, including allowable overhead rates, they may reduce or delay their purchases of our products and services, which could have a material effect on our revenues and financial performance.**

– BRKR FY24 10K (Filed March 3, 2025)

Source: Capital IQ, public filings

Merger Consideration

(in thousands, except for number of shares and per share amounts):	
Akoya common stock outstanding as of September 30, 2024	49,522
Exchange Ratio	0.318
Total shares of Quanterix common stock to be issued	15,748
Quanterix stock price	\$ 7.83
Fair value of Quanterix common stock transferred to Akoya shareholders ^(a)	\$ 123,309
Akoya Restricted Stock Units outstanding	1,926
Exchange Ratio	0.318
Total shares of Quanterix common stock to be issued following acceleration of vesting up on change in control	613
Quanterix stock price	\$ 7.83
Fair value of Akoya Restricted Stock Units converted to Quanterix Restricted Stock Units ^(b)	\$ 4,796
Replacement of Akoya stock options by Quanterix ^(c)	\$ 13,649
Cash to be paid in lieu of fractional shares issued by Quanterix ^(d)	\$ 49
Total preliminary fair value of consideration transferred	\$ 141,803

Notes:

- a) Represents the preliminary estimated fair value of 15,748,228 shares of Quanterix common stock, issued to Akoya stockholders, who held 49,522,728 common shares of Akoya as of September 30, 2024. The fair value of those shares has been estimated using \$7.83 per share, which was the last reported sale price of Quanterix shares on the Nasdaq Global Market on February 10, 2025.
- b) Represents the preliminary estimated fair value of 612,538 shares of Quanterix common stock, issued to holders of Akoya's RSU as purchase consideration upon the effectiveness of the Merger. The fair value of those shares has been estimated using \$7.83 per share, which was the last reported sale price of Quanterix shares on the Nasdaq Global Market on February 10, 2025.
- c) Represents the preliminary estimated fair value of Quanterix stock options as replacement to holders of outstanding and unexercised stock option holders of Akoya. The fair value of the replacement awards has been estimated using the Black Scholes option pricing model. Of that amount, \$13.6 million was allocated to purchase consideration upon Merger based on the portion attributable to pre-combination employee services and the remaining amount of \$4 million is allocated to future employee services will be expensed as stock-based compensation on a straight-line basis over the remaining service periods of those awards.
- d) Represents the estimated fair value of the fractional shares to be paid by cash to Akoya award holders as purchase consideration upon the effectiveness of the Merger.

Source: QTRX Form S-4 Prospectus

Piper Sandler Thinks the Deal Makes Sense – But They Own >7% of AKYA

Piper Sandler owns >3.5M AKYA shares through the firm's merchant banking arm

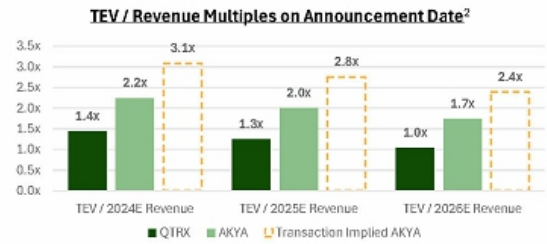
Name and Address of Beneficial Owner	Number of Shares Beneficially Owned	Percentage of Shares Beneficially Owned(1)
5% and Greater Stockholders		
Entities affiliated with Telegraph Hill Partners(2)	17,675,247	35.8%
Entities affiliated with PSC Capital Partners LLC(3)	3,537,161	7.2%
Entities affiliated with aMoon Growth Fund II, L.P.(4)	3,077,195	6.2%
Named Executive Officers and Directors		
Brian McKelligon(5)	1,083,219	2.1%
Johnny Ek(6)	106,666	*
Frederic Pla(7)	355,932	*
Myla Lai-Goldman, M.D.(8)	76,544	*
Scott Mendel(9)	82,757	*
Thomas Raffin, M.D.(10)	282,001	*
Thomas P. Schnettler(11)	47,409	*
Robert Shepler(12)	417,001	*
Matthew Winkler, Ph.D.(13)	1,031,922	2.1%
All executive officers and directors as a group (11 persons)(14)	3,726,556	7.3%

Source: Capital IQ, public filings, and QTRX S-4 Prospectus filed 02/13/25

QTRX Shareholders Deserved a Market Check

In a stock merger, Quanterix and Akoya are swapping ownership – both are effectively sellers – yet it’s unclear if the QTRX Board fully explored available alternatives

- For the merger to be accretive to Quanterix shareholders, AKYA shares (the acquiree) **MUST be undervalued relative to the acquirer (QTRX)**
- The 0.318x exchange ratio implies that the Quanterix Board believe Akoya is worth >2x more as a multiple of 2024 revenue
- QTRX formed a Transaction Committee in October 2023 to explore strategic alternatives – yet the company **provided no disclosures** regarding:
 - Specific strategic options evaluated
 - Recommendations made by the Committee
 - Timeline or conclusion of the review process
- **It remains unclear if the QTRX Board fully reviewed all available strategic alternatives.**



“ In a cash deal, the roles of the two parties are clear-cut, and the exchange of money for shares completes a simple transfer of ownership. **But in an exchange of shares, it becomes far less clear who is the buyer and who is the seller. In some cases, the shareholders of the acquired company can end up owning most of the company that bought their shares.** Companies that pay for their acquisitions with stock **share both the value and the risks** of the transaction with the shareholders of the company they acquire. The decision to use stock instead of cash **can also affect shareholder returns.** In studies covering more than 1,200 major deals, researchers have consistently found that, at the time of announcement, **shareholders of acquiring companies fare worse in stock transactions than they do in cash transactions.** What’s more, the findings show **that early performance differences between cash and stock transactions become greater—much greater—over time.** ”

– Harvard Business Review (November 1999)¹

Source: Capital IQ, and public filings

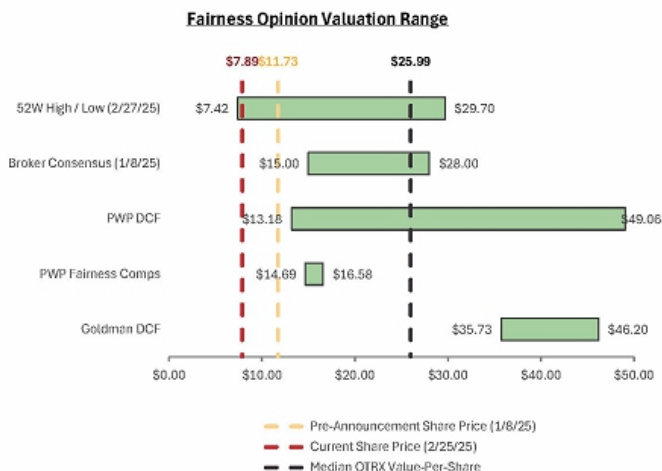
1) [Stock or Cash?: The Trade-Offs for Buyers and Sellers in Mergers and Acquisitions by Alfred Rappaport and Mark L. Sirower](#)

2) TEV/Revenue multiple calculated using the closing share price of QTRX and AKYA common stock as of 1/8/25. Net debt calculated as total debt less unrestricted cash and securities. AKYA net debt is adjusted to reflect the \$6.9M early exit fee as disclosed in the S-4 filing. [47]

S4 Disclosures Demonstrate QTRX Is Undervalued as a Standalone Company

Both transaction advisors acknowledge that QTRX is materially undervalued – A strong reason for Quanterix NOT to issue equity to buy struggling Akoya.

- Both PWP and Goldman value QTRX equity **far above its current trading price**
- The median share price across the PWP/GS fairness opinions, sell-side consensus, and QTRX's 52-week trading range is ~\$26/share
- At the 0.318x exchange ratio, QTRX is effectively valuing AKYA at nearly \$10/share – **more than 5x FY26 consensus revenue and >400% above the current share price**²
- QTRX's share price has declined by **over 30%** since the announcement, clearly signaling shareholders **oppose the transaction**
- QTRX **nearly trades for a 10% premium to its own** cash. In contrast, AKYA – flagged for **going concern risk**¹ – is being valued with an enterprise value at **>5x that of QTRX**



Source: Capital IQ, and public filings

PWP Fairness Comps include QTRX, AKYA, MASS, CTKB, ONT, PACB, QSI, SEER, and LAB

1) AKYA [3Q24 10Q](#)

2) Calculated as justified AKYA enterprise value less \$47.0M assumed net debt, divided by total AKYA shares outstanding at 9/30/24; relative to closing AKYA share price dated 2/25/25.

[48]

Akoya's Own Banker (PWP) Suggests the Ratio Is Too High

PWP's analysis of publicly-traded peers suggests the agreed upon Exchange Ratio is **nearly 2x above fair market value**

	PWP Implied Exchange Ratio Range		
	High	Low	Average
EV / 2024E Revenue	0.1618x	0.1738x	0.1678x
EV / 2025E Revenue	0.1968x	0.2140x	0.2054x
Average Implied Exchange Ratio	0.1866x		
Actual Exchange Ratio	0.3180x		
Over (Under)	1.7x		

PWP Selected Publicly-Traded Companies Analysis¹

The ranges of implied values per share derived from these calculations are summarized in the following table:

	Akoya Share Price Range	Quanterix Share Price Range
EV / 2024E Revenue	\$ 2.39 - \$2.86	\$ 14.75 - \$16.45
EV / 2025E Revenue	\$ 2.88 - \$3.53	\$ 14.62 - \$16.51

PWP compared (i) these implied ranges for Akoya to the Akoya Current Share Price of \$2.66 and the Implied Offer Price of \$3.73 and (ii) these implied ranges for Quanterix to the Quanterix Current Share Price of \$11.73.

PWP then calculated the exchange ratio ranges implied by the selected publicly-traded companies analysis. For each of the foregoing analyses, PWP calculated (i) the ratio of the highest implied value per share for Akoya derived from the selected publicly-traded companies analysis to the highest implied value per share for Quanterix derived from the selected publicly-traded companies analysis and (ii) the ratio of the lowest implied value per share for Akoya derived from the selected publicly-traded companies analysis to the lowest implied value per share for Quanterix derived from the selected publicly-traded companies analysis to calculate the following ranges of implied exchange ratios and ranges of implied pro forma ownership of the Combined Company:

	Implied Exchange Ratio Range	Implied Akoya Pro Forma Ownership	Implied Quanterix Pro Forma Ownership
EV / 2024E Revenue	0.1618x - 0.1738x	18% - 19%	81% - 82%
EV / 2025E Revenue	0.1968x - 0.2140x	21% - 22%	78% - 79%

PWP compared these implied exchange ratio ranges to (i) the Exchange Ratio of 0.3180 shares of Quanterix Common Stock to be received for each share of Akoya Common Stock as provided for in the Merger Agreement and (ii) the At Market Implied Exchange Ratio of 0.2268 shares of Quanterix Common Stock to be received for each share of Akoya Common Stock.

Source: Capital IQ, and public filings

1) QTRX Form S-4 Prospectus page 163

PWP Analysis, Continued

And undervalues Quanterix relative to both management teams' expectations

PWP Selected Publicly-Traded Companies Analysis¹

PWP then calculated the exchange ratio ranges implied by the discounted cash flow analysis. For each of the foregoing analyses, PWP calculated (i) the ratio of the highest implied value per share for Akoya derived from the discounted cash flow analysis to the highest implied value per share for Quanterix derived from the discounted cash flow analysis and (ii) the ratio of the lowest implied value per share for Akoya derived from the discounted cash flow analysis to the lowest implied value per share for Quanterix derived from the discounted cash flow analysis to calculate the following ranges of implied exchange ratios and ranges of the implied pro forma ownership of the Combined Company:

	Implied Exchange Ratio Range	Implied Akoya Pro Forma Ownership	Implied Quanterix Pro Forma Ownership
Akoya Street Estimates/ Quanterix Street Estimates	0.0814x - 0.1168x	9% - 13%	87%-91%
Akoya Mgmt. (15%-35% PTS) Case/ Quanterix Mgmt. Base Case	0.2258x - 0.3365x	23%-31%	69% - 77%
Akoya Mgmt. (15%-35% PTS) Case / Quanterix Mgmt. Upside Case	0.1487x - 0.2089x	16% - 22%	78% - 84%
Akoya Mgmt. (Excl. CDx) Case/ Quanterix Mgmt. RUO Case	0.2479x - 0.2879x	25% - 28%	72%-75%

PWP compared these implied exchange ratio ranges to (i) the Exchange Ratio of 0.3180 shares of Quanterix Common Stock to be received for each share of Akoya Common Stock as provided for in the Merger Agreement and (ii) the At Market Implied Exchange Ratio of 0.2268 shares of Quanterix Common Stock to be received for each share of Akoya Common Stock.

Source: Capital IQ, and public filings

1) QTRX Form S-4 Prospectus page 165 and 186: "For each Akoya and Quanterix, PWP performed a discounted cash flow analysis, which is a method of deriving an implied value range for a company's equity securities based on the sum of the company's unlevered free cash flows over a forecast period and the terminal value at the end of the forecast period. In connection with this analysis, PWP used the Akoya Street Estimates, Akoya Mgmt. (15%-35% PTS) Case, and the Akoya Mgmt. (Excl. CDx) Case for Akoya, and the Quanterix Street Estimates, Quanterix Mgmt. RUO Case, the Quanterix Mgmt. Base Case and the Quanterix Mgmt. Upside Case for Quanterix." [50]

AKYA Consistently Misses Wallstreet Estimates

AKYA has reported a consistent string of misses, resulting in numerous analyst downgrades

- **BTIG downgrades from Buy to Hold in August 2024**
- **Morgan Stanley downgrades from Buy to Hold in August 2024**
- Craig Hallum initiates with a Buy in June 2024
 - **Craig Hallum downgrades to Hold in November 2024**

Analyst Quotes

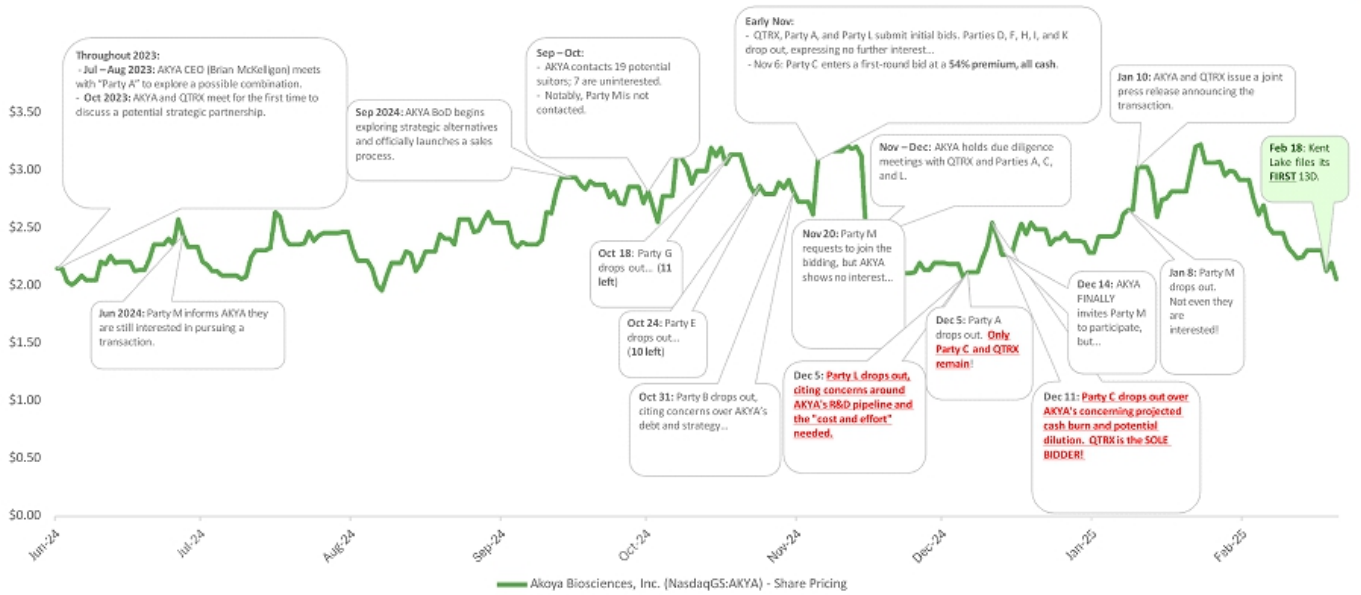
“ We believe the deal was motivated by two key issues: (1) **we think AKYA needed to raise money again** (we think it just simply didn't raise enough money when it did its last two raises); and (2) AKYA had been struggling like many other life sciences tools companies. **AKYA has now gone three-for-three on quarterly revenue misses and guide-downs this year**

– BTIG (August 2024) ”

Broker	Date	Broker Raise (Lower) %	Broker Value Current (\$M)	Broker Value Prior (\$M)
Entitlement Needed	1/28/2025	-1.1%	82.3	83.2
Entitlement Needed	1/10/2025	0.0%	83.2	83.2
Entitlement Needed	11/25/2024	-14.2%	83.2	97.0
Piper Sandler Companies	11/18/2024	-16.2%	80.7	96.3
Craig-Hallum Capital Group LLC	11/15/2024	-18.7%	80.5	99.0
Entitlement Needed	11/15/2024	-0.6%	82.5	83.0
Entitlement Needed	11/15/2024	0.0%	97.0	97.0
Entitlement Needed	11/14/2024	-17.6%	80.8	98.0
Guggenheim Securities, LLC	11/14/2024	-13.9%	82.1	95.4
Entitlement Needed	11/14/2024	-16.1%	83.0	98.9
BTIG	11/14/2024	-17.9%	83.0	101.1
Entitlement Needed	11/14/2024	0.0%	97.0	97.0
Craig-Hallum Capital Group LLC	11/14/2024	-5.2%	99.0	104.4
BTIG	11/14/2024	-2.8%	101.1	104.0
Entitlement Needed	9/16/2024	0.0%	97.0	97.0
Entitlement Needed	8/14/2024	-2.1%	97.0	99.1
Piper Sandler Companies	8/12/2024	-5.6%	96.3	102.0
Entitlement Needed	8/6/2024	-0.1%	98.9	99.0
Craig-Hallum Capital Group LLC	8/6/2024	-5.2%	99.0	104.4
Entitlement Needed	8/6/2024	-4.8%	99.1	104.1
BTIG	8/6/2024	-2.8%	101.1	104.0
Guggenheim Securities, LLC	8/5/2024	-4.3%	95.4	99.7
Capital One Securities, Inc.	8/5/2024	-8.1%	96.5	105.0
Entitlement Needed	8/5/2024	-5.9%	98.0	104.1
Entitlement Needed	8/5/2024	-6.7%	99.0	106.1
Entitlement Needed	8/5/2024	-10.3%	104.1	116.0
Craig-Hallum Capital Group LLC	8/5/2024	0.0%	104.4	-
Capital One Securities, Inc.	8/5/2024	-8.7%	105.0	115.0
Entitlement Needed	8/5/2024	0.1%	106.1	106.0
Entitlement Needed	7/22/2024	-4.8%	99.1	104.1
Craig-Hallum Capital Group LLC	6/21/2024	0.0%	104.4	-
Piper Sandler Companies	5/14/2024	-11.3%	102.00	115.00

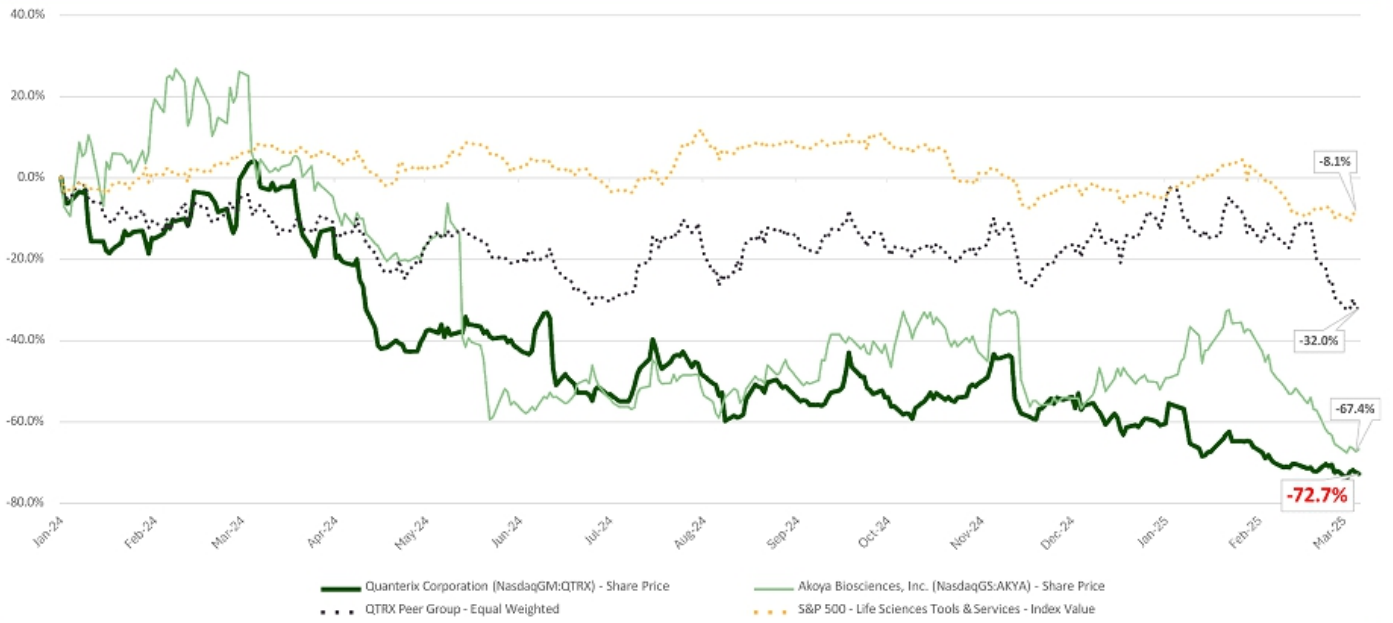
Source: Capital IQ, public filings, and analyst research reports

All AKYA Bidders Dropped Before QTRX Final Offer



Source: Capital IQ, and public filings

QTRX Relative Underperformance



"Peer Group" classified as equal weighted index of dividend-adjusted share price for MASS, ABSI, AKYA, TECH, CTKB, EVT, MXCT, NAUT, NEO, ONT, PACB, PSNL, QSI, SEER, SOPH, LAB, and TWST. Share price data as of March 7, 2025. Source: Capital IQ, and public filings

TSR Has Been Abysmal Under Quanterix's Existing Board and CEO

Total Stockholder Returns: 1-, 3-, 5-Year and CEO Tenure

	1-Year	3-Year	5-Year	CEO Tenure
Quanterix	-73.6%	-73.2%	-65.2%	-75.7%
Russell 2000 Index (^RUT)	2.8%	7.8%	61.8%	1.9%
S&P 500 Health Care (Sector) (^HCX)	3.5%	20.5%	67.3%	13.4%

Total Stockholder Returns: Director Tenure

	Masoud Toloue (3 Years)	William P. Donnelly (2 Years)	Jeffrey T. Elliot (1 Year)	Karen A. Flynn (3 Years)	Sarah E. Hlavinka (6 Years)	Martin D. Madaus (10+ Years)	Ivana Magovčević-Liebisch (<1 Year)	Paul M. Meister (10+ Years)	David R. Walt (10+ Years)
Quanterix	-75.7%	-70.3%	-47.3%	-55.1%	-79.1%	-57.6%	-40.8%	-57.6%	-57.6%
Russell 2000 Index (^RUT)	1.9%	6.8%	-5.1%	9.6%	45.1%	44.5%	-4.8%	44.5%	44.5%
S&P 500 Health Care (^HCX)	13.4%	14.8%	-0.4%	20.1%	77.3%	107.3%	-2.1%	107.3%	107.3%

Source: Capital IQ. Total stockholder return data as of 3/7/25.