#### UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

#### FORM 8-K

#### **CURRENT REPORT**

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

Date of Report (Date of earliest event reported): May 2, 2024

### **QUANTERIX CORPORATION**

(Exact name of registrant as specified in its charter)

**Delaware** (State or other jurisdiction of incorporation)

001-38319 (Commission File Number) **20-8957988** (IRS Employer Identification No.)

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

**01821** (Zip Code)

(617) 301-9400

(Registrant's telephone number, including area code)

Check the appropriate box below if the Foregistrant under any of the following prov		ly satisfy the filing obligation of the
□Written communications pursuant to Ru	le 425 under the Securities Act (17 CFR 2	230.425)
☐ Soliciting material pursuant to Rule 14a	1-12 under the Exchange Act (17 CFR 240	.14a-12)
☐ Pre-commencement communications pu	ursuant to Rule 14d-2(b) under the Exchan	ge Act (17 CFR 240.14d-2(b))
☐ Pre-commencement communications pu	ursuant to Rule 13e-4(c) under the Exchan	ge Act (17 CFR 240.13e-4(c))
Securities registered pursuant to Section 1	2(b) of the Act:	
Title of each class	Trading symbol(s)	Name of each exchange on which registered
Common Stock, \$0.001 par value per share	QTRX	The Nasdaq Global Market
Indicate by check mark whether the regist of 1933 (17 CFR §230.405) or Rule 12b-2		
Emerging Growth Company $\square$		
If an emerging growth company, indicate period for complying with any new or re Exchange Act. □		

#### Item 2.02 Results of Operations and Financial Condition.

On May 7, 2024, Quanterix Corporation ("Quanterix") issued a press release announcing its financial results for the first quarter ended March 31, 2024 (the "Earnings Release"). A copy of the Earnings Release is furnished as Exhibit 99.1 and is incorporated herein by reference.

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

## Item 5.02 Departure of Directors or Certain Officers; Election of Directors; Appointment of Certain Officers; Compensatory Arrangements of Certain Officers.

(b) On May 2, 2024, Brian J. Blaser notified Quanterix of his resignation from its Board of Directors, effective as of May 3, 2024. His resignation from all public company boards on which he served was a condition of his appointment as President and Chief Executive Officer, and as a director, of QuidelOrtho Corporation as of May 6, 2024. His resignation was not because of any disagreement with Quanterix on any matter relating to its operations, policies or practices.

#### Item 9.01 Financial Statements and Exhibits.

(d) Exhibits

Exhibit	
No.	Description
99.1	Earnings Release dated May 7, 2024.
104	Cover Page Interactive Data File (embedded within the inline XBRL document)

#### SIGNATURES

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

#### QUANTERIX CORPORATION

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

Date: May 7, 2024

#### Quanterix Releases Financial Results for the First Ouarter of 2024

**BILLERICA**, Mass. – May 7, 2024 - Quanterix Corporation (NASDAQ: QTRX), a company fueling scientific discovery through ultrasensitive biomarker detection, today announced financial results for the first quarter ended March 31, 2024.

"Building on the foundation our team developed last year, Quanterix is entering a new innovation and product delivery phase," said Masoud Toloue, Chief Executive Officer of Quanterix. "We've announced the first of several new assays that could re-define the way Tau is measured, from the brain versus peripheral sources. We will continue to lead in bringing novel biomarker detection solutions to market for the research community; developing tests to study and enable the early detection of neuro-pathologies such as Multiple Sclerosis, Parkinson's, and Alzheimer's disease."

#### First Quarter Financial Highlights

- Revenue of \$32.1 million, an increase of 12.7% compared to \$28.5 million in the prior year period.
- GAAP gross margin of 61.2%, an increase of approximately 170 basis points compared to the prior year period and non-GAAP gross margin of 54.5%, an increase of approximately 140 basis points compared to the prior year period.
- Net loss of \$10.1 million, an increase of \$4.0 million compared to the corresponding prior year period.
- Net cash usage in the quarter was \$19.4 million, and the Company ended the period with \$304.5 million of cash, cash equivalents, marketable securities, and restricted cash.

#### **Operational and Business Highlights**

- The Company continues to execute against its transformation initiatives, successfully commercializing five new assays in the quarter.
- The Company announced its Simoa® phospho-Tau 217 (P-tau217) blood test was granted Breakthrough Device designation by the U.S. Food and Drug Administration (FDA) as an aid in diagnostic evaluation of Alzheimer's Disease. The FDA's Breakthrough Device designation is granted to products that have the potential to offer more effective diagnosis of life-threatening diseases with an unmet medical need.
- Eli Lilly Clinical Diagnostics Laboratory launched <u>CertuitAD</u>, an *in vitro* immunoassay that measures plasma tau protein fragments phosphorylated at threonine 217 (P-tau217), using the Quanterix SP-X Imaging and Analysis System<sup>TM</sup>. CertuitAD is intended to be used in patients aged 60 years and older who present with cognitive impairment and who are being evaluated for Alzheimer's disease (AD) and other causes of cognitive decline. CertuitAD is a laboratory developed test (LDT) and is not intended to be used as a screening or stand-alone diagnostic test and is not intended for therapeutic monitoring.

#### 2024 Full Year Business Outlook

The Company's 2024 financial guidance is unchanged. Management continues to expect full-year 2024 revenue in the range of \$139 to \$144 million. This guide is for the Company's research-only business, and does not include revenues from Diagnostics testing, which to date have not been material. GAAP gross margin percentage is expected to be in the range of 57-61%, and non-GAAP gross margin percentage is expected to be in the range of \$1-55%. Finally, the Company continues to anticipate 2024 cash usage in the range of \$25 to \$30 million.

For additional information on the non-GAAP financial measures included in this press release, please see "Use of Non-GAAP Financial Measures" and "Reconciliation of GAAP to Non-GAAP Financial Measures" below.

#### **Conference Call**

In conjunction with this announcement, the Company will host a conference call on May 08, 2024 at 8:30 a.m. E.T. <u>Click here to preregister for the conference call and obtain your dial-in number and passcode.</u>

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

# Quanterix Corporation Consolidated Statements of Operations (Unaudited and in thousands, except per share data)

	 Three Months Ended March 31,			
	2024		2023	
Revenues:				
Product revenue	\$ 19,670	\$	19,287	
Service and other revenue	11,967		8,579	
Collaboration and license revenue	155		368	
Grant revenue	 274		222	
Total revenues	32,066		28,456	
Costs of goods sold and services:				
Cost of product revenue	7,145		7,033	
Cost of service and other revenue	 5,295		4,497	
Total costs of goods sold and services	 12,440		11,530	
Gross profit	19,626		16,926	
Operating expenses:				
Research and development	6,675		4,720	
Selling, general, and administrative	25,993		20,850	
Other lease costs	924		776	
Total operating expenses	33,592		26,346	
Loss from operations	 (13,966)		(9,420)	
Interest income, net	3,948		3,449	
Other income, net	206		8	
Loss before income taxes	 (9,812)		(5,963)	
Income tax expense	(260)		(140)	
Net loss	\$ (10,072)	\$	(6,103)	
Net loss per common share, basic and diluted	\$ (0.26)	\$	(0.16)	
•	 <u> </u>		` '	
Weighted-average common shares outstanding, basic and diluted	 38,126		37,327	

#### Quanterix Corporation Consolidated Balance Sheets (Unaudited and in thousands)

	March 31, 2024		December 31, 2023		
ASSETS					
Current assets:					
Cash and cash equivalents	\$ 45,281	\$	174,422		
Marketable securities	256,640		146,902		
Accounts receivable	29,276		25,414		
Inventory	26,015		22,365		
Prepaid expenses and other current assets	9,551		9,291		
Total current assets	 366,763		378,394		
Restricted cash	2,605		2,604		
Property and equipment, net	17,492		17,926		
Intangible assets, net	5,339		6,034		
Operating lease right-of-use assets	17,748		18,251		
Other non-current assets	1,802		1,802		
Total assets	\$ 411,749	\$	425,011		
LIABILITIES AND STOCKHOLDERS' EQUITY	 				
Current liabilities:					
Accounts payable	\$ 3,914	\$	5,048		
Accrued compensation and benefits	6,706		13,659		
Accrued expenses and other current liabilities	7,021		6,041		
Deferred revenue	10,234		9,468		
Operating lease liabilities	4,366		4,241		
Total current liabilities	 32,241		38,457		
Deferred revenue, net of current portion	933		1,227		
Operating lease liabilities, net of current portion	36,084		37,223		
Other non-current liabilities	1,053		1,177		
Total liabilities	 70,311		78,084		
Total stockholders' equity	 341,438		346,927		
Total liabilities and stockholders' equity	\$ 411,749	\$	425,011		

#### Quanterix Corporation Consolidated Statements of Cash Flows (Unaudited and in thousands)

		Three Months Ended March 31,		
		2024		2023
Cash flows from operating activities:				
Net loss	\$	(10,072)	\$	(6,103)
Adjustments to reconcile net loss to net cash used in operating activities:				
Depreciation and amortization expense		1,523		1,439
Credit losses on accounts receivable		176		178
Accretion of marketable securities		(1,657)		_
Operating lease right-of-use asset amortization		478		334
Stock-based compensation expense		5,265		3,902
Other operating activity		55		270
Changes in assets and liabilities:				
Accounts receivable		(4,233)		(3,741)
Inventory		(3,670)		(89)
Prepaid expenses and other current assets		(254)		(422)
Other non-current assets		(21)		(33)
Accounts payable		(1,122)		(1,271)
Accrued compensation and benefits, accrued expenses, and other current liabilities		(6,126)		(5,983)
Deferred revenue		472		2,041
Operating lease liabilities		(988)		179
Other non-current liabilities		10		(203)
Net cash used in operating activities		(20,164)		(9,502)
Cash flows from investing activities:				
Purchases of marketable securities		(137,889)		_
Proceeds from maturities of marketable securities		29,200		_
Purchases of property and equipment		(506)		(136)
Net cash used in investing activities		(109,195)		(136)
Cash flows from financing activities:				
Proceeds from common stock issued under stock plans		2,037		564
Payments for employee taxes withheld on stock-based compensation awards		(1,438)		(13)
Net cash provided by financing activities		599		551
Net decrease in cash, cash equivalents, and restricted cash		(128,760)		(9,087)
Effect of exchange rate changes on cash, cash equivalents, and restricted cash		(380)		24
Cash, cash equivalents, and restricted cash at beginning of period		177,026		341,337
Cash, cash equivalents, and restricted cash at end of period	\$	47,886	\$	332,274
Cubit, cubit equitations, and restricted cubit at one of period	Ψ	17,000	<del>)</del>	332,211

#### **Use of Non-GAAP Financial Measures**

To supplement our financial statements presented on a U.S. GAAP basis, we present non-GAAP gross profit, non-GAAP gross margin, non-GAAP total operating expenses, and non-GAAP loss from operations, which are calculated by including shipping and handling costs for product sales within cost of goods sold instead of within selling, general, and administrative expenses. Management uses these non-GAAP measures to evaluate our operating performance in a manner that allows for meaningful period-to-period comparison and analysis of trends in our business and our competitors. Management believes that presentation of these non-GAAP measures provides useful information to investors in assessing our operating performance within our industry and in order to allow comparability to the presentation of other companies in our industry where shipping and handling costs are included in cost of goods sold for products. The non-GAAP financial information presented here should be considered in conjunction with, and not as a substitute for, the financial information presented in accordance with U.S. GAAP. Set forth below is a reconciliation of non-GAAP gross profit, non-GAAP gross margin, non-GAAP total operating expenses, and non-GAAP loss from operations to their most directly comparable GAAP financial measures.

#### Reconciliation of GAAP to Non-GAAP Financial Measures

## Quanterix Corporation Reconciliation of GAAP Financial Measures to Non-GAAP Financial Measures (Unaudited and in thousands, except percentages)

		Three Months Ended March 31,				
		2024		2023		
GAAP gross profit	\$	19,626	\$	16,926		
Shipping and handling costs (1)		(2,142)		(1,829)		
Non-GAAP gross profit	\$	17,484	\$	15,097		
GAAP revenue	\$	32,066	\$	28,456		
GAAP gross margin (gross profit as % of revenue)		61.2%		59.5%		
Non-GAAP gross margin (non-GAAP gross profit as % of revenue)		54.5%		53.1%		
GAAP total operating expenses	S	33,592	\$	26,346		
Shipping and handling costs (1)	•	(2,142)		(1,829)		
Non-GAAP total operating expenses	\$	31,450	\$	24,517		
GAAP loss from operations	\$	(13,966)	\$	(9,420)		
Non-GAAP loss from operations	\$	(13,966)	\$	(9,420)		

(1) Shipping and handling costs, which include freight and other activities costs associated with product shipments, are captured within operating expenses in our consolidated statements of operations. During the three months ended March 31, 2024 and 2023, we incurred \$2.1 million and \$1.8 million, respectively, of shipping and handling costs recorded within operating expenses.

#### **About Quanterix**

From discovery to diagnostics, Quanterix's ultrasensitive biomarker detection is driving breakthroughs only made possible through its unparalleled sensitivity and flexibility. The Company's Simoa technology has delivered the gold standard for earlier biomarker detection in blood, serum or plasma, with the ability to quantify proteins that are far lower than the Level of Quantification of conventional analog methods. Its industry-leading precision instruments, digital immunoassay technology and CLIA-certified Accelerator laboratory have supported research that advances disease understanding and management in neurology, oncology, immunology, cardiology and infectious disease. Quanterix has been a trusted partner of the scientific community for nearly two decades, powering research published in more than 2,900 peer-reviewed journals. Find additional information about the Billerica, Massachusetts-based company at <a href="https://www.quanterix.com">https://www.quanterix.com</a> or follow us on <a href="mailto:Twitter">Twitter</a> and <a href="mailto:LinkedIn">LinkedIn</a>.

#### Forward-Looking Statements

Quanterix's current financial results, as discussed in this press release, are preliminary and unaudited, and subject to adjustment. This press release contains forward-looking statements within the meaning of the Private Securities Litigation Reform Act of 1995. Words such as "may," "will," "expect," "plan," "anticipate," "estimate," "intend" and similar expressions (as well as other words or expressions referencing future events, conditions or circumstances) are intended to identify forward-looking statements. These forward-looking statements include, but are not limited to, statements about Quanterix's financial performance, including statements under the header "2024 Full Year Business Outlook" set forth above, and are subject to a number of risks, uncertainties and assumptions. Forward-looking statements in this press release are based on Quanterix's expectations and assumptions as of the date of this press release. Each of these forward-looking statements involves risks and uncertainties. Factors that may cause Quanterix's actual results to differ from those expressed or implied in the forward-looking statements in this press release include, but are not limited to, those described in our periodic reports filed with the U.S. Securities and Exchange Commission, including the "Risk Factors" sections contained therein. Except as required by law, Quanterix assumes no obligation to update any forward-looking statements contained herein to reflect any change in expectations, even as new information becomes available.

#### **Contacts**

#### Media:

PAN Communications Maya Nimnicht 510-334-6273 quanterix@pancomm.com

Investor Relations: Francis Pruell, Quanterix (508)-789-1725 ir@quanterix.com