
**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): June 3, 2022

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)

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

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

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

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

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

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

Emerging Growth Company

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

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

On June 3, 2022, the board of directors (the “Board”) of Quanterix Corporation (the “Company”) appointed Karen A. Flynn to serve on the Board, effective as of June 6, 2022. Ms. Flynn will serve as a Class I director with a term that will expire at the Company’s 2024 annual meeting of stockholders.

There is no arrangement or understanding between Ms. Flynn and any other person pursuant to which she was selected as a director of the Company, and there is no family relationship between Ms. Flynn and any of the Company’s other directors or executive officers. Ms. Flynn does not have a material interest in any transaction that is required to be disclosed under Item 404(a) of Regulation S-K.

In accordance with the Company’s non-employee director compensation policy (the “Policy”), Ms. Flynn will receive an equity award valued at \$400,000 comprised of (i) 60% non-qualified stock options to purchase the Company’s common stock at an exercise price equal to the closing price of the common stock on June 6, 2022 and (ii) 40% RSUs. The options vest over three years from the date of grant, with one-third vesting on the first anniversary and the remainder vesting over the following two years in 24 successive equal monthly installments. The RSUs vest over a three-year period, with one-third vesting on each of the first, second, and third anniversaries of the grant date. Ms. Flynn will also receive annual compensation and annual equity awards as further provided in the Policy. A description of the Policy is available in the Company’s proxy statement filed with the Securities and Exchange Commission on April 29, 2022 (File No. 001-38319) under the heading “Executive Officer and Director Compensation – Director Compensation – Non-Employee Director Compensation Policy.”

A copy of the press release dated June 6, 2022 announcing the appointment of Ms. Flynn is filed as Exhibit 99.1 hereto and is incorporated by reference herein.

Item 9.01 Financial Statements and Exhibits.

(d) Exhibits

Exhibit No.	Description
99.1	Press Release dated June 6, 2022.
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/ John Fry

John Fry
General Counsel and Secretary

Date: June 6, 2022

Quanterix

Quanterix Announces Appointment of Karen A. Flynn to Board of Directors

BILLERICA, Mass. – June 6, 2022 – Quanterix Corporation (NASDAQ: QTRX), a company digitizing biomarker analysis with the goal of advancing the science of precision health, today announced that Karen A. Flynn, Senior Vice President and Chief Commercial Officer of Catalent Pharma Solutions (NYSE: CTLT), has been appointed to Quanterix' Board of Directors effective June 6, 2022.

Ms. Flynn is an accomplished healthcare leader with over 35 years of commercial, operations and general management experience in the pharmaceutical services industry. She currently serves as Chief Commercial Officer of Catalent Pharma Solutions, a global provider of development and manufacturing solutions for drugs, biologics, cell and gene therapies and consumer health products. She previously served as the President of Catalent's Biologics business. Prior to Catalent, she served as Senior Vice President and Chief Commercial Officer for West Pharmaceutical Services and as President of West's Pharmaceutical Packaging Systems business. In addition to her experience leading high growth businesses, Ms. Flynn brings an extensive background in commercial strategy, strategic planning and innovation and quality management. She previously served on the Board of Directors of Recro Pharma, Inc., a specialty pharma and contract development and manufacturing organization focused on developing non-opioid products for acute pain and on solving complex formulation and manufacturing challenges primarily in small molecule therapeutic development.

"We are pleased to welcome Karen to the Quanterix Board, bringing proven expertise that will be invaluable as a Director," said Kevin Hrusovsky, Executive Chairman of the Board of Quanterix. "As Quanterix continues executing on its next phase of growth, we are confident that Karen's successful track record of developing and overseeing commercial and market strategies for global businesses will complement the skill sets already on our Board. We look forward to Karen's guidance and perspective as the Company continues to expand the impact it has in powering advances in healthcare research and diagnostics."

"I am excited to join the Quanterix Board of Directors and work alongside a group of talented and distinguished board members during such an exciting time at the Company," said Karen Flynn. "Quanterix is driving innovation that has the potential to deliver life-changing solutions in support of earlier disease detection, better prognoses and enhanced treatment methods. With the Company's strong foundation and focused vision for the future, Quanterix is well-positioned to further scale and strengthen its business and drive long-term value for shareholders. I look forward to supporting Quanterix' continued success."

Ms. Flynn received her Master of Science in Engineering from University of Pennsylvania and Master of Science in Business Administration from Boston University. She received her Bachelor of Science in Pre-Professional Studies from the University of Notre Dame. Ms. Flynn has completed the Women's Senior Leadership Program and the Corporate Governance, Effectiveness and Accountability in the Board Room Program at Kellogg School of Management at Northwestern University and has been an active leader in the EcSell Institute Advisory Board, the Forum of Executive Women, the Healthcare Businesswomen's Association and the Advisory Board for Downingtown STEM Academy. She has also received numerous awards, including the Silver Stevie Award for Women in Business and the Ellen Ann Roberts Business Award.

Quanterix

About Quanterix

Quanterix is a company that's digitizing biomarker analysis with the goal of advancing the science of precision health. The company's digital health solution, Simoa, has the potential to change the way in which healthcare is provided today by giving researchers the ability to closely examine the continuum from health to disease. Quanterix' technology is designed to enable much earlier disease detection, better prognoses and enhanced treatment methods to improve the quality of life and longevity of the population for generations to come. The technology is currently being used for research applications in several therapeutic areas, including oncology, neurology, cardiology, inflammation and infectious disease. The company was established in 2007 and is located in Billerica, Massachusetts. For additional information, please visit <https://www.quanterix.com>.

Forward-Looking Statements

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. Forward-looking statements in this news release are based on Quanterix' 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' actual results to differ from those expressed or implied in the forward-looking statements in this press release are discussed in Quanterix' filings 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 Contact:

PAN Communications
Lauren Force, 617-502-4366
pan.quanterix@pancomm.com

Investor Relations Contact:

Stephen Hrusovsky
(774) 278-0496
shrusovsky@quanterix.com
