
**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): April 27, 2021

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 April 27, 2021, the board of directors (the “Board”) of Quanterix Corporation (the “Company”) appointed Laura J. Olson to serve on the Board, effective immediately. Ms. Olson will serve as a Class I director with a term that will expire at the Company’s 2021 annual meeting of stockholders.

There is no arrangement or understanding between Ms. Olson and any other person pursuant to which she or he was selected as a director of the Company, and there is no family relationship between Ms. Olson and any of the Company’s other directors or executive officers. Ms. Olson 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. Olson was granted an option for 15,800 shares of the Company’s common stock and 4,540 restricted stock units (“RSUs”) subject to the vesting and other terms as provided therein. Ms. Olson will also receive cash compensation and annual equity awards as further provided in the Policy.

A copy of the press release dated April 28, 2021 announcing the appointment of Ms. Olson is filed as Exhibit 99.1 hereto and is incorporated by reference herein.

Item 9.01 Financial Statements and Exhibits

(d) Exhibits

Exhibit

<u>No.</u>	<u>Description</u>
99.1	Press Release dated April 28, 2021.
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: April 28, 2021

Quanterix

Quanterix Appoints Laurie Olson to Board of Directors

Pharmaceutical veteran adds deep expertise in commercial and corporate strategy

Billerica, Mass. – April 28, 2021 — Quanterix Corporation (NASDAQ: QTRX), a company digitizing biomarker analysis with the goal of advancing the science of precision health, today announced the appointment of Laurie Olson to its Board of Directors, effective April 27, 2021.

Olson is a seasoned pharmaceutical executive who brings more than three decades of experience in commercial and corporate strategy from Pfizer, Inc. to the Quanterix Board. She most recently served as executive vice president, strategy and commercial operations at Pfizer until 2018, where she led a team of approximately 2,000 people across 100 markets, with a scope including corporate strategy, market access and reimbursement, digital strategy, portfolio management, business analytics and global commercial operations. She also served as a member of Pfizer’s executive leadership team.

Olson also serves on the Board of Directors of Karuna Therapeutics, a clinical stage biopharmaceutical company committed to developing novel therapies for neuropsychiatric disorders.

“We are pleased to welcome Laurie to our board as we continue to scale the business and capitalize on the tremendous market opportunity ahead,” said Kevin Hrusovsky, Chairman, Chief Executive Officer, and President of Quanterix. “Laurie’s philosophy that the needs of patients are best realized via deep customer insights and data-driven decisions integrates seamlessly with Quanterix’ broader precision health mission. We believe her business acumen and far-reaching expertise in commercial and corporate strategy will be a valuable addition to our very skilled and knowledgeable board.”

During her career at Pfizer, Olson also served as a member of the Pfizer Board of Directors Science and Technology Committee, where she informed therapeutic area strategies and the utilization of new technologies and analytics to drive R&D productivity. As a member of the Company’s Business Development, Portfolio, Disclosure, Diversity and Compliance Committees, Olson developed M&A strategies, ensured rigorous capital allocation and maintained Pfizer’s transparency and compliance commitments with stakeholders.

“I am honored to be appointed to Quanterix’ Board of Directors at such a pivotal moment for biomarker research, when technologies like Simoa are increasingly being called upon to detect, measure and inform disease treatment,” said Olson. “I look forward to working alongside Kevin and the board, utilizing the expertise I’ve gained across a number of roles within the pharmaceutical industry, to advance Quanterix’ mission of transforming the way healthcare is provided.”

Olson is also a member of the Board of Trustees at the Mystic Seaport Museum in Mystic, CT. She received her BA in economics from State University of New York at Stony Brook and earned her MBA in marketing from Hofstra University.

To learn more about Quanterix, visit www.quanterix.com/about.

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.quoterix.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.

Media Contact:

PAN Communications

Staci Didner, (407) 734-7325

quoterix@pancomm.com
