
**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 24, 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))

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.

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

Item 5.07 Submission of Matters to a Vote of Security Holders.

At the 2021 Annual Meeting of Stockholders of Quanterix Corporation (the “**Company**”) held on June 24, 2021 (the “**Annual Meeting**”), the stockholders of the Company voted on and approved the following matters, which are described in detail in the Company’s Definitive Proxy Statement filed with the Securities and Exchange Commission on April 29, 2021 (File No. 001-38319): (1) to elect Martin D. Madaus, Ph.D. and Laurie J. Olson as Class I directors to each serve for a three-year term expiring at the Company’s annual meeting of stockholders in 2024, and until their successors have been elected and qualified, or until their earlier death, resignation, retirement or removal (“**Proposal 1**”); and (2) to ratify the appointment of Ernst & Young LLP as the Company’s independent registered public accounting firm for the fiscal year ending December 31, 2021 (“**Proposal 2**”).

The tabulation of votes with respect to the proposals at the Annual Meeting was as follows:

Proposal 1 — Election of Directors:

	<u>For</u>	<u>Withheld</u>	<u>Broker Non-Votes</u>
Martin D. Madaus, Ph.D.	14,802,359	12,563,292	4,842,888
Laurie J. Olson	27,308,589	57,062	4,842,888

Proposal 2 — Ratification of Independent Registered Public Accounting Firm:

<u>For</u>	<u>Against</u>	<u>Abstain</u>
32,117,134	64,079	27,326

Item 9.01 Financial Statements and Exhibits

(d) Exhibits

<u>Exhibit No.</u>	<u>Description</u>
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 25, 2021
