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 7, 2023

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

is intended to simultaneously satisfy the filing	obligation of the registrant under any of the			
 □ 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:				
Trading symbol(s)	Name of each exchange on which registered			
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).				
α if the registrant has elected not to use the extension to Section 13(a) of the Exchange Act. \Box	ended transition period for complying with any new			
	tr the Securities Act (17 CFR 230.425) the Exchange Act (17 CFR 240.14a-12) the 14d-2(b) under the Exchange Act (17 CFR tule 13e-4(c) under the Exchange Act (17 CFR tt: Trading symbol(s) QTRX Trading growth company as defined in Rule 405 CFR §240.12b-2).			

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

At the 2023 Annual Meeting of Stockholders of Quanterix Corporation (the "Company") held on June 7, 2023 (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 14, 2023 (File No. 001-38319) (the "Proxy Statement"): (1) to elect Brian J. Blaser and Paul M. Meister as Class III directors to each serve for a three-year term expiring at the Company's annual meeting of stockholders in 2026, and until their successors have been elected and qualified, or until their earlier death, resignation, retirement or removal ("Proposal 1"); (2) an advisory vote to approve the compensation of the Company's named executive officers, as disclosed in the Proxy Statement ("Proposal 2"); and (3) to ratify the appointment of Ernst & Young LLP as the Company's independent registered public accounting firm for the fiscal year ending December 31, 2023 ("Proposal 3").

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

Proposal 1 — Election of Directors:

	For	Withheld	Broker Non-Votes
Brian J. Blaser	26,710,969	231,019	3,771,384
Paul M. Meister	17,804,718	9,137,270	3,771,384

Proposal 2 — Advisory Vote to Approve the Compensation of the Company's Named Executive Officers:

For	Against	Abstain	Broker Non-Votes
26,690,270	239.915	11.803	3,771,384

<u>Proposal 3 — Ratification of Independent Registered Public Accounting Firm:</u>

For	Against	Abstain
30,669,815	37,672	5,885

Item 9.01 Financial Statements and Exhibits.

Description

(d) Exhibits

104

Exhibit No.

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/ Michael A. Doyle

Michael A. Doyle Chief Financial Officer

Date: June 8, 2023