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): September 4, 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 following provisions:	; is intended to simultaneously satis	fy the filing obligation of the registrant under any of the
☐ Written communications pursuant to Rule 425 under the	ne Securities Act (17 CFR 230.425)	
☐ Soliciting material pursuant to Rule 14a-12 under the I	Exchange Act (17 CFR 240.14a-12)	
☐ Pre-commencement communications pursuant to Rule	14d-2(b) under the Exchange Act (17	7 CFR 240.14d-2(b))
☐ Pre-commencement communications pursuant to Rule	13e-4(c) under the Exchange Act (17	7 CFR 240.13e-4(c))
Indicate by check mark whether the registrant is an emergi Rule 12b-2 of the Securities Exchange Act of 1934 (17 CF)		ale 405 of the Securities Act of 1933 (17 CFR §230.405) or
Emerging Growth Company ⊠		
If an emerging growth company, indicate by check mark is or revised financial accounting standards provided pursuan	•	the extended transition period for complying with any new ct. \square
Securities registered pursuant to Section 12(b) of the Act:		
Title of each class Common Stock, \$0.001 par value per share	Trading symbol(s) QTRX	Name of each exchange on which registered The Nasdaq Global Market

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

(a) On September 4, 2021, Quanterix Corporation (the "Company") received notice from Marijn Dekkers that he was resigning as a member of the Company's Board of Directors and all committees thereof effective December 31, 2021. His resignation was not because of any disagreement with the Company on matters relating to its operations, policies or practices.

Item 9.01 Financial Statements and Exhibits.

(d) Exhibits

Exhibit

No. Description

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: September 10, 2021