# UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

#### **SCHEDULE 14A**

Proxy Statement Pursuant to Section 14(a) of the Securities Exchange Act of 1934 (Amendment No. )

Filed by the Registrant  $\,x\,$ 

Filed by a Party other than the Registrant o

Check the appropriate box:

- o Preliminary Proxy Statement
- o Confidential, for Use of the Commission Only (as permitted by Rule 14a-6(e)(2))
- o Definitive Proxy Statement
- x Definitive Additional Materials
- Soliciting Material under §240.14a-12

## **Quanterix Corporation**

(Name of Registrant as Specified In Its Charter)

(Name of Person(s) Filing Proxy Statement, if other than the Registrant)

Payment of Filing Fee (Check the appropriate box):

- x No fee required.
- o Fee computed on table below per Exchange Act Rules 14a-6(i)(1) and 0-11.
  - (1) Title of each class of securities to which transaction applies:
  - (2) Aggregate number of securities to which transaction applies:
  - (3) Per unit price or other underlying value of transaction computed pursuant to Exchange Act Rule 0-11 (set forth the amount on which the filing fee is calculated and state how it was determined):
  - (4) Proposed maximum aggregate value of transaction:
  - (5) Total fee paid:
- o Fee paid previously with preliminary materials.
- Check box if any part of the fee is offset as provided by Exchange Act Rule 0-11(a)(2) and identify the filing for which the offsetting fee was paid previously. Identify the previous filing by registration statement number, or the Form or Schedule and the date of its filing.
  - (1) Amount Previously Paid:
  - (2) Form, Schedule or Registration Statement No.:
  - (3) Filing Party:
  - (4) Date Filed:



000004

ENDORSEMENT LINE SACKPACK Սլիրմյանորդորդյիրիմ**իցիՍԱլի**նդինվյինիլցինդներաննան

MR A SAMPLE DESIGNATION (IF ANY)

ADD 1

ADD 2

ADD 3 ADD 4

ADD 5



in the shaded bar below.

#### Online



Go to www.investorvote.com/QTRX or scan the QR code - login details are located



Votes submitted electronically must be received by 1:00 a.m., Eastern Time, on July 25, 2019.

# 2019 Stockholder Meeting Notice

1234 5678 9012 345

Important Notice Regarding the Availability of Proxy Materials for the Quanterix Corporation Stockholder Meeting to be held on July 25, 2019.

Under Securities and Exchange Commission rules, you are receiving this notice that the proxy materials for the annual stockholders' meeting are available on the Internet. Follow the instructions below to view the materials and vote online or request a copy. The items to be voted on and location of the annual meeting are on the reverse side. Your vote is important!

This communication presents only an overview of the more complete proxy materials that are available to you on the Internet. We encourage you to access and review all of the important information contained in the proxy materials before voting. The proxy statement and 2018 annual report to stockholders are available at:

# www.investorvote.com/QTRX



# Easy Online Access — View your proxy materials and vote.

- Step 1: Go to www.investorvote.com/QTRX.
- Step 2: Click on the icon on the right to view meeting materials.
- Step 3: Return to the investorvote.com window and follow the instructions on the screen to log in.
- Step 4: Make your selections as instructed on each screen for your delivery preferences.
- Step 5: Vote your shares.

When you go online, you can also help the environment by consenting to receive electronic delivery of future materials.



Obtaining a Copy of the Proxy Materials - If you want to receive a copy of the proxy materials, you must request one. There is no charge to you for requesting a copy. Please make your request as instructed on the reverse side on or before July 15, 2019 to facilitate timely delivery.







## 2019 Stockholder Meeting Notice

Quanterix Corporation's Annual Meeting of Stockholders will be held on July 25, 2019 at Quanterix Corporation, 900 Middlesex Turnpike, Building 1, Billerica, MA 01821, at 10:00 a.m. Eastern Time.

Proposals to be voted on at the meeting are listed below along with the Board of Directors' recommendations.

The Board of Directors recommends a vote FOR all the nominees listed and FOR Proposal 2:

- 1. Election of Class II Directors for a three-year term expiring in 2022:
  - 01 Marijn Dekkers, Ph.D.
  - 02 Sarah E. Hlavinka
  - 03 David R. Walt, Ph.D.
- Ratification of the appointment of Ernst & Young LLP as the independent registered public accounting firm for the fiscal year ending December 31, 2019.
- Such other business as may properly come before the meeting or an adjournment or postponement thereof.

PLEASE NOTE - YOU CANNOT VOTE BY RETURNING THIS NOTICE. To vote your shares you must go online or request a paper copy of the proxy materials to receive a proxy card. If you wish to attend and vote at the meeting, please bring this notice with you.



### Here's how to order a copy of the proxy materials and select delivery preferences:

Current and future delivery requests can be submitted using the options below.

If you request an email copy, you will receive an email with a link to the current meeting materials.

PLEASE NOTE: You must use the number in the shaded bar on the reverse side when requesting a copy of the proxy materials.

- Internet Go to www.investorvote.com/QTRX.
- Phone Call us free of charge at 1-866-641-4276.
- Email Send an email to investorvote@computershare.com with "Proxy Materials Quanterix Corporation" in the subject line. Include
  your full name and address, plus the number located in the shaded bar on the reverse side, and state that you want a paper copy of
  the meeting materials.

To facilitate timely delivery, requests for a paper copy of proxy materials must be received by July 15, 2019.