
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

Filed by a Party other than the Registrant

Check the appropriate box:

- Preliminary Proxy Statement
 Confidential, for Use of the Commission Only (as permitted by Rule 14a-6(e)(2))
 Definitive Proxy Statement
 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):

- No fee required.
 Fee paid previously with preliminary materials.
 Fee computed on table in exhibit required by Item 25(b) per Exchange Act Rules 14a-6(i)(1) and 0-11.
-
-



C 1234567890



000004

ENDORSEMENT_LINE _____ SACKPACK _____

MR A SAMPLE
DESIGNATION (IF ANY)
ADD 1
ADD 2
ADD 3
ADD 4
ADD 5
ADD 6

Online



Go to www.investorvote.com/QTRX
or scan the QR code – login details are located
in the shaded bar below.

2022 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 June 23, 2022.**

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 at the virtual 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 2022 proxy statement and our 2021 annual report to stockholders are available for viewing, printing and downloading 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 June 13, 2022 to facilitate timely delivery.



2 N O T

C O Y



03NIVC

2022 Stockholder Meeting Notice

Quanterix Corporation's Annual Meeting of Stockholders will be held on June 23, 2022, at 10:00 a.m. Eastern Time, virtually via the internet at meetnow.global/MD2UGTW. To access the virtual meeting, you must have the 15 digit number that is printed in the shaded bar located on the reverse side of this form.

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, **FOR** Proposals 2 and 4 and **FOR 1 YEAR** on Proposal 3:

1. Election of Class II Directors for a three-year term expiring in 2025:
01 - Sarah E. Hlavinka
02 - Masoud Toloue, Ph.D.
03 - David R. Walt, Ph.D.
2. To approve, by a non-binding advisory vote, the compensation of our named executive officers, as disclosed in the proxy statement.
3. To approve, by a non-binding advisory vote, the frequency of holding an advisory vote on compensation of our named executive officers.
4. Ratification of the appointment of Ernst & Young LLP as the independent registered public accounting firm for the fiscal year ending December 31, 2022.

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.



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 June 13, 2022.