

Use these links to rapidly review the document

[TABLE OF CONTENTS](#)

[Table of Contents](#)

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 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: _____
 - 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: _____
-

Quanterix™

June 5, 2019

To Our Stockholders:

You are cordially invited to attend the 2019 annual meeting of stockholders of Quanterix Corporation to be held at 10:00 a.m. ET on Thursday, July 25, 2019 at our offices, located at 900 Middlesex Turnpike, Building 1, Billerica, Massachusetts 01821.

Details regarding the meeting, the business to be conducted at the meeting, and information about Quanterix Corporation that you should consider when you vote your shares are described in the accompanying proxy statement.

At the annual meeting, three persons will be elected to our board of directors. In addition, we will ask stockholders to ratify the appointment of Ernst & Young LLP as our independent registered public accounting firm for our fiscal year ending December 31, 2019. The board of directors recommends the approval of each of these proposals. Such other business will be transacted as may properly come before the annual meeting.

Under Securities and Exchange Commission rules that allow companies to furnish proxy materials to stockholders over the Internet, we have elected to deliver our proxy materials to the majority of our stockholders over the Internet. This delivery process allows us to provide stockholders with the information they need, while at the same time conserving natural resources and lowering the cost of delivery. On or about June 6, 2019, we intend to begin sending to our stockholders a Notice of Internet Availability of Proxy Materials (the "Notice") containing instructions on how to access our proxy statement for our 2019 Annual Meeting of Stockholders and our 2018 annual report to stockholders. The Notice also provides instructions on how to vote online or by telephone and how to receive a paper copy of the proxy materials by mail.

We hope you will be able to attend the annual meeting. Whether you plan to attend the annual meeting or not, it is important that you cast your vote either in person or by proxy. You may vote over the Internet as well as by telephone or by mail. When you have finished reading the proxy statement, you are urged to vote in accordance with the instructions set forth in the proxy statement. We encourage you to vote by proxy so that your shares will be represented and voted at the meeting, whether or not you can attend.

Thank you for your continued support of Quanterix Corporation. We look forward to seeing you at the annual meeting.

Sincerely,

/s/ E. KEVIN HRUSOVSKY

E. Kevin Hrusovsky
Chairman, President and Chief Executive Officer

QUANTERIX CORPORATION

900 Middlesex Turnpike
Billerica, Massachusetts 01821

June 5, 2019

NOTICE OF 2019 ANNUAL MEETING OF STOCKHOLDERS

TIME: 10:00 a.m. ET

DATE: Thursday, July 25, 2019

PLACE: Quanterix Corporation, 900 Middlesex Turnpike, Building 1, Billerica, Massachusetts 01821

PURPOSES:

1. To elect three directors to serve three-year terms expiring in 2022;
2. To ratify the appointment of Ernst & Young LLP as our independent registered public accounting firm for the fiscal year ending December 31, 2019; and
3. To transact such other business that is properly presented at the annual meeting and any adjournments or postponements thereof.

WHO MAY VOTE:

You may vote if you were the record owner of Quanterix Corporation common stock at the close of business on June 3, 2019. A list of stockholders of record will be available at the annual meeting and, during the 10 days prior to the annual meeting, at our principal executive offices, located at 900 Middlesex Turnpike, Building 1, Billerica, Massachusetts 01821.

All stockholders are cordially invited to attend the annual meeting. **Whether you plan to attend the annual meeting or not, we urge you to vote by following the instructions in the Notice of Internet Availability of Proxy Materials that you previously received and submit your proxy by the Internet, telephone or mail in order to ensure the presence of a quorum.** You may change or revoke your proxy at any time before it is voted at the meeting.

BY ORDER OF THE BOARD OF DIRECTORS

/s/ BRIAN P. KEANE

Brian P. Keane
General Counsel and Secretary

TABLE OF CONTENTS

	<u>Page</u>
Important Information About the Annual Meeting and Voting	1
Security Ownership of Certain Beneficial Owners and Management	7
Management and Corporate Governance	10
Executive Officer and Director Compensation	21
Equity Compensation Plan Information	26
Report of Audit Committee	27
Section 16(a) Beneficial Ownership Reporting Compliance	28
Certain Relationships and Related Person Transactions	29
Election of Directors	34
Ratification of the Appointment of the Independent Registered Public Accounting Firm	35
Corporate Code of Conduct and Ethics	37
Other Matters	37
Stockholder Proposals and Nominations for Director	37

QUANTERIX CORPORATION

900 Middlesex Turnpike
Billerica, Massachusetts 01821

PROXY STATEMENT FOR QUANTERIX CORPORATION 2019 ANNUAL MEETING OF STOCKHOLDERS TO BE HELD ON JULY 25, 2019

This proxy statement, along with the accompanying Notice of 2019 Annual Meeting of Stockholders, contains information about the 2019 annual meeting of stockholders of Quanterix Corporation, including any adjournments or postponements of the annual meeting. We are holding the annual meeting at 10:00 a.m. ET on Thursday, July 25, 2019 at our offices, located at 900 Middlesex Turnpike, Building 1, Billerica, Massachusetts 01821.

In this proxy statement, we refer to Quanterix Corporation as "Quanterix," "the Company," "we" and "us."

This proxy statement relates to the solicitation of proxies by our board of directors for use at the annual meeting.

On or about June 6, 2019, we intend to begin sending the Important Notice Regarding the Availability of Proxy Materials to stockholders entitled to vote at the annual meeting.

IMPORTANT NOTICE REGARDING THE AVAILABILITY OF PROXY MATERIALS FOR THE STOCKHOLDER MEETING TO BE HELD ON JULY 25, 2019

This proxy statement and our 2018 annual report to stockholders are available for viewing, printing and downloading at www.investorvote.com/qtrx. To view these materials, please have your 15-digit control number(s) available that appears on your Notice or proxy card. On this website, you can also elect to receive future distributions of our proxy statements and annual reports to stockholders by electronic delivery.

Additionally, you can find a copy of our Annual Report on Form 10-K, which includes our financial statements, for the fiscal year ended December 31, 2018 on the website of the Securities and Exchange Commission, or the SEC, at www.sec.gov, or in the "SEC Filings" section of the "Financial Information" section of the "Investors" section of our website at www.quanterix.com. You may also obtain a printed copy of our Annual Report on Form 10-K, including our financial statements, free of charge, from us by sending a written request to: Quanterix Corporation, Attn: Investor Relations, 900 Middlesex Turnpike, Building 1, Billerica, Massachusetts 01821. Exhibits will be provided upon written request and payment of an appropriate processing fee.

IMPORTANT INFORMATION ABOUT THE ANNUAL MEETING AND VOTING

Why is the Company Soliciting My Proxy?

Our board of directors is soliciting your proxy to vote at the 2019 annual meeting of stockholders to be held at our offices, located at 900 Middlesex Turnpike, Building 1, Billerica, Massachusetts 01821, on Thursday, July 25, 2019, at 10:00 a.m. ET, and any adjournments or postponements of the meeting, which we refer to as the annual meeting. This proxy statement, along with the accompanying Notice of Annual Meeting of Stockholders, summarizes the purposes of the meeting and the information you need to know to vote at the annual meeting.

We have made available to you on the Internet or have sent you this proxy statement, the Notice of Annual Meeting of Stockholders, the proxy card and a copy of our Annual Report on Form 10-K for the fiscal year ended December 31, 2018 because you owned shares of our common stock on the record date. We intend to commence distribution of the Important Notice Regarding the Availability of Proxy Materials, which we refer to throughout this proxy statement as the Notice, and, if applicable, proxy materials to stockholders on or about June 6, 2019.

Why Did I Receive a Notice in the Mail Regarding the Internet Availability of Proxy Materials Instead of a Full Set of Proxy Materials?

As permitted by the rules of the U.S. Securities and Exchange Commission, or the SEC, we may furnish our proxy materials to our stockholders by providing access to such documents on the Internet, rather than mailing printed copies of these materials to each stockholder. Most stockholders will not receive printed copies of the proxy materials unless they request them. We believe that this process should expedite stockholders' receipt of proxy materials, lower the costs of the annual meeting and help to conserve natural resources. If you received a Notice by mail or electronically, you will not receive a printed or email copy of the proxy materials, unless you request one by following the instructions included in the Notice. Instead, the Notice instructs you as to how you may access and review all of the proxy materials and submit your proxy on the Internet. If you requested a paper copy of the proxy materials, you may authorize the voting of your shares by following the instructions on the proxy card, in addition to the other methods of voting described in this proxy statement.

Who Can Vote?

Only stockholders who owned our common stock at the close of business on June 3, 2019 are entitled to vote at the annual meeting. On this record date, there were 24,896,311 shares of our common stock outstanding and entitled to vote. Our common stock is our only class of voting stock.

You do not need to attend the annual meeting to vote your shares. Shares represented by valid proxies, received in time for the annual meeting and not revoked prior to the annual meeting, will be voted at the annual meeting. For instructions on how to change or revoke your proxy, see "May I Change or Revoke My Proxy?" below.

How Many Votes Do I Have?

Each share of our common stock that you own entitles you to one vote.

How Do I Vote?

Whether you plan to attend the annual meeting or not, we urge you to vote by proxy. All shares represented by valid proxies that we receive through this solicitation, and that are not revoked, will be voted in accordance with your instructions on the proxy card or as instructed via Internet or telephone. You may specify whether your shares should be voted for or withheld for each nominee for director and whether your shares should be voted for, against or abstain with respect to the other proposal. If you properly submit a proxy without giving specific voting instructions, your shares will be voted in accordance with the board of director's recommendations as noted below. Voting by proxy will not affect your right to attend the annual meeting. If your shares are registered directly in your name through our stock transfer agent, Computershare Trust Company N.A., or you have stock certificates registered in your name, you may vote:

- **By Internet or by telephone.** Follow the instructions included in the Notice or, if you received printed materials, in the proxy card to vote over the Internet or by telephone.
- **By mail.** If you received a proxy card by mail, you can vote by mail by completing, signing, dating and returning the proxy card as instructed on the card. If you sign the proxy card but do not specify how you want your shares voted, they will be voted in accordance with the board of director's recommendations as noted below.
- **In person at the meeting.** If you attend the meeting, you may deliver a completed proxy card in person or you may vote by completing a ballot, which will be available at the meeting.

Telephone and Internet voting facilities for stockholders of record will be available 24 hours a day and will close at 1:00 a.m. ET on July 25, 2019.

If your shares are held in "street name" (held in the name of a bank, broker or other holder of record), you will receive instructions from the holder of record. You must follow the instructions of the holder of record in order for your shares to be voted. Telephone and Internet voting also will be offered to stockholders owning shares through certain banks and brokers. If your shares are not registered in your own name and you plan to vote your shares in person at the annual meeting, you should contact your broker or agent to obtain a legal proxy or broker's proxy card and bring it to the annual meeting in order to vote.

How Does the Board of Directors Recommend That I Vote on the Proposals?

The board of directors recommends that you vote as follows:

- "FOR" the election of the three nominees for director; and
- "FOR" the ratification of the appointment of Ernst & Young LLP as our independent registered public accounting firm for our fiscal year ending December 31, 2019.

If any other matter is presented at the annual meeting, your proxy provides that your shares will be voted by the proxy holder listed in the proxy in accordance with his best judgment. At the time this proxy statement was first made available, we knew of no matters that needed to be acted on at the annual meeting, other than those discussed in this proxy statement.

May I Change or Revoke My Proxy?

If you give us your proxy, you may change or revoke it at any time before the annual meeting. You may change or revoke your proxy in any one of the following ways:

- if you received a proxy card, by signing a new proxy card with a date later than your previously delivered proxy and submitting it as instructed above;
- by re-voting by Internet or by telephone as instructed above;
- by notifying our Corporate Secretary in writing before the annual meeting that you have revoked your proxy; or
- by attending the annual meeting in person and voting in person. Attending the annual meeting in person will not in and of itself revoke a previously submitted proxy. You must specifically request at the annual meeting that it be revoked.

Your most current vote, whether by telephone, Internet or proxy card is the one that will be counted.

What if I Receive More Than One Notice or Proxy Card?

You may receive more than one Notice or proxy card if you hold shares of our common stock in more than one account, which may be in registered form or held in street name. Please vote in the manner described above under "How Do I Vote?" for each account to ensure that all of your shares are voted.

Will My Shares be Voted if I Do Not Vote?

If your shares are registered in your name or if you have stock certificates, they will not be counted if you do not vote as described above under "How Do I Vote?" If your shares are held in street name and you do not provide voting instructions to the bank, broker or other nominee that holds

your shares as described above, the bank, broker or other nominee that holds your shares has the authority to vote your unvoted shares only on the ratification of the appointment of our independent registered public accounting firm without receiving instructions from you. Therefore, we encourage you to provide voting instructions to your bank, broker or other nominee. This ensures your shares will be voted at the annual meeting and in the manner you desire. A "broker non-vote" will occur if your broker cannot vote your shares on a particular matter because it has not received instructions from you and does not have discretionary voting authority on that matter or because your broker chooses not to vote on a matter for which it does have discretionary voting authority.

What Vote is Required to Approve Each Proposal and How are Votes Counted?

Proposal 1: Elect Directors	The three nominees for director who receive the most votes (also known as a "plurality" of the votes cast) will be elected. You may vote either FOR all of the nominees, WITHHOLD your vote from all of the nominees or WITHHOLD your vote from any one or more of the nominees. Votes that are withheld will not be included in the vote tally for the election of the directors. Brokerage firms do not have authority to vote customers' unvoted shares held by the firms in street name for the election of the directors. As a result, any shares not voted by a customer will be treated as a broker non-vote. Such broker non-votes will have no effect on the results of this vote.
--	---

Proposal 2: Ratify Appointment of Independent Registered Public Accounting Firm	The affirmative vote of a majority of the votes cast affirmatively or negatively for this proposal is required to ratify the appointment of our independent registered public accounting firm. Abstentions will have no effect on the results of this vote. Brokerage firms have authority to vote customers' unvoted shares held by the firms in street name on this proposal. If a broker does not exercise this authority, such broker non-votes will have no effect on the results of this vote. We are not required to obtain the approval of our stockholders to appoint our independent registered public accounting firm. However, if our stockholders do not ratify the appointment of Ernst & Young LLP as our independent registered public accounting firm for 2019, the audit committee of our board of directors will reconsider its appointment.
--	---

Is Voting Confidential?

We will keep all the proxies, ballots and voting tabulations private. We only let our Inspector of Election, Computershare Trust Company, N.A., examine these documents. Management will not know how you voted on a specific proposal unless it is necessary to meet legal requirements. We will, however, forward to management any written comments you make on the proxy card or that you otherwise provide.

Where Can I Find the Voting Results of the Annual Meeting?

The preliminary voting results will be announced at the annual meeting, and we will publish preliminary results, or final results if available, in a Current Report on Form 8-K within four business days of the annual meeting. If final results are unavailable at the time we file the Form 8-K, then we will file an amended Current Report on Form 8-K to disclose the final voting results within four business days after the final voting results are known.

What Are the Costs of Soliciting these Proxies?

We will pay all of the costs of soliciting these proxies. Our directors and employees may solicit proxies in person or by telephone, fax or email. We will pay these employees and directors no additional compensation for these services. We will ask banks, brokers and other institutions, nominees and fiduciaries to forward these proxy materials to their principals and to obtain authority to execute proxies. We will then reimburse them for their expenses.

What Constitutes a Quorum for the Annual Meeting?

The presence, in person or by proxy, of the holders of a majority of the voting power of all outstanding shares of our common stock entitled to vote at the annual meeting is necessary to constitute a quorum at the annual meeting. Votes of stockholders of record who are present at the annual meeting in person or by proxy, abstentions, and broker non-votes are counted for purposes of determining whether a quorum exists.

Attending the Annual Meeting

The annual meeting will be held at 10:00 a.m. ET on Thursday, July 25, 2019 at our offices, located at 900 Middlesex Turnpike, Building 1, Billerica, Massachusetts 01821. When you arrive, signs will direct you to the appropriate meeting rooms. You need not attend the annual meeting in order to vote.

Householding of Annual Disclosure Documents

SEC rules concerning the delivery of annual disclosure documents allow us or your broker to send a single Notice or, if applicable, a single set of our proxy materials to any household at which two or more of our stockholders reside, if we or your broker believe that the stockholders are members of the same family. This practice, referred to as "householding," benefits both you and us. It reduces the volume of duplicate information received at your household and helps to reduce our expenses. The rule applies to our Notices, annual reports, proxy statements and information statements. Once you receive notice from your broker or from us that communications to your address will be "household," the practice will continue until you are otherwise notified or until you revoke your consent to the practice. Stockholders who participate in householding will continue to have access to and utilize separate proxy voting instructions.

If your household received a single Notice or, if applicable, a single set of proxy materials this year, but you would prefer to receive your own copy, please contact our transfer agent, Computershare Trust Company, N.A., by calling their toll free number, 1-800-736-3001.

If you do not wish to participate in "householding" and would like to receive your own Notice or, if applicable, set of our proxy materials in future years, follow the instructions described below. Conversely, if you share an address with another stockholder and together both of you would like to receive only a single Notice or, if applicable, set of proxy materials, follow these instructions:

- If your shares are registered in your own name, please contact our transfer agent, Computershare Trust Company, N.A., and inform them of your request by calling them

at 1-800-736-3001 or writing them at Computershare Trust Company, N.A., P.O. Box 505000, Louisville, KY 40233-5000.

- If a broker or other nominee holds your shares, please contact the broker or other nominee directly and inform them of your request. Be sure to include your name, the name of your brokerage firm and your account number.

Electronic Delivery of Company Stockholder Communications

Most stockholders can elect to view or receive copies of future proxy materials over the Internet instead of receiving paper copies in the mail.

You can choose this option and save the Company the cost of producing and mailing these documents by:

- following the instructions provided on your Notice or proxy card;
- following the instructions provided when you vote over the Internet; or
- going to www.computershare.com and logging into the Investor Center to follow the instructions provided.

SECURITY OWNERSHIP OF CERTAIN BENEFICIAL OWNERS AND MANAGEMENT

The following table sets forth certain information with respect to the beneficial ownership of our common stock as of May 15, 2019 for (a) the executive officers named in the Summary Compensation Table on page 21 of this proxy statement, (b) each of our directors and director nominees, (c) all of our current directors and executive officers as a group and (d) each stockholder known by us to own beneficially more than 5% of our common stock. Beneficial ownership is determined in accordance with the rules of the SEC and includes voting or investment power with respect to the securities. We deem shares of common stock that may be acquired by an individual or group within 60 days of May 15, 2019 pursuant to the exercise of options, warrants or the vesting of restricted stock units to be outstanding for the purpose of computing the percentage ownership of such individual or group, but are not deemed to be outstanding for the purpose of computing the percentage ownership of any other person shown in the table. Except as indicated in footnotes to this table, we believe that the stockholders named in this table have sole voting and investment power with respect to all shares of common stock shown to be beneficially owned by them based on information provided to us by these stockholders. Percentage of ownership is based on 22,608,199 shares of common stock outstanding on May 15, 2019.

<u>Name and Address of Beneficial Owner</u>	<u>Shares Beneficially Owned(1)</u>	
	<u>Number</u>	<u>Percent</u>
<i>Principal Stockholders</i>		
ARCH Venture Partners and affiliated funds(1) 8755 West Higgins Avenue, Suite 1025 Chicago, IL 60631	3,949,684	17.5%
Bain Capital Ventures Entities(2) 200 Clarendon Street Boston, MA 02116	2,136,085	9.4%
bioMérieux SA(3) 376 Chemin de l'Orme, 69280 Marcy l'Étoile, France	2,055,632	9.1%
Flagship Ventures Fund 2004, L.P.(4) 55 Cambridge Parkway, Suite 800E Cambridge, MA 02142	2,030,495	9.0%
<i>Directors, Director Nominees and Named Executive Officers</i>		
E. Kevin Hrusovsky(5)	1,146,482	5.0%
Douglas G. Cole, M.D.(6)	10,170	*
John M. Connolly(7)	12,870	*
Keith L. Crandell(8)	3,960,356	17.5%
Marijn Dekkers, Ph.D.(9)	137,916	*
Martin D. Madaus, Ph.D.(10)	322,028	1.4%
Sarah E. Hlavinka	—	—
Paul M. Meister(11)	34,092	*
David Walt, Ph.D.(12)	1,225,670	5.4%
Dawn R. Mattoon, Ph.D.(13)	25,826	*
Jackson Streeter, M.D.	—	—
All directors and current executive officers as a group (13 persons)(14)	7,178,438	30.8%

* Represents beneficial ownership of less than 1% of the outstanding shares of our common stock.

(1) Based on a Schedule 13G/A filed by ARCH Venture Fund VI, L.P. ("ARCH Venture Fund VI"); ARCH Venture Fund VIII Overage, L.P. ("AVF VIII Overage"); ARCH Venture Partners VI, L.P.

("AVP VI LP"); ARCH Venture Partners VI, LLC ("AVP VI LLC"); ARCH Venture Partners VIII, LLC ("AVP VIII LLC"); and Keith Crandell ("Crandell"), Clinton Bybee ("Bybee") and Robert Nelsen ("Nelsen") (collectively, the "Managing Directors" and individually, each a "Managing Director") with the SEC on February 14, 2019, which reported ownership as of December 31, 2018. ARCH Venture Fund VI is the record owner of 2,412,000 shares of common stock, which includes 10,333 shares of common stock issuable upon the exercise of warrants (the "ARCH VI Shares") and AVF VIII Overage is the record holder of 1,537,684 shares of common stock (the "ARCH Overage Shares"); combined with ARCH VI Shares, the "Arch Shares") as of December 31, 2018. AVP VI LP, as the sole general partner of ARCH Venture Fund VI, may be deemed to beneficially own the AVF VI Shares. AVP VI LLC, as the sole general partner of AVP VI LP, may be deemed to beneficially own the AVF VI Shares. AVP VIII LLC, as the sole general partner of AVF VIII Overage, may be deemed beneficially own the ARCH Overage Shares. As managing directors of AVP VI LLC and AVP VIII LLC, each Managing Director may also be deemed to share the power to direct the disposition and vote of the ARCH Shares. The Managing Directors disclaim beneficial ownership of the ARCH Shares, except to the extent of any pecuniary interest therein.

- (2) Based on a Schedule 13G filed by (1) Bain Capital Venture Fund 2005, L.P. ("Fund 2005"), (2) BCIP Associates III, LLC ("BCIP III"), and (3) BCIP Associates III-B, LLC ("BCIP III-B" and together with Fund 2005 and BCIP III, the "Bain Capital Entities") with the SEC on February 14, 2018, which reported ownership as of December 31, 2017. Bain Capital Venture Investors, LLC, a Delaware limited liability company ("BCVI"), is the general partner of Bain Capital Venture Partners 2005, L.P. ("BCVP"), which is the general partner of Fund 2005. Boylston Coinvestors, LLC ("Boylston"), is the managing partner of (i) BCIP Associates III ("BCIPA III"), which is the manager of BCIP III, and (ii) BCIP Associates III-B ("BCIPA III-B"), which is the manager of BCIP III-B. The governance, investment strategy and decision-making process with respect to the investments held by all of the Bain Capital Entities is directed by the Executive Committee of BCVI, which consists of Michael A. Krupka and Ajay Agarwal. As a result, BCVI and Messrs. Krupka and Agarwal may be deemed to share voting and dispositive power with respect to all of the shares of common stock held by all of the Reporting Persons. As of December 31, 2017, Fund 2005 held 1,865,010 shares of common stock, which includes 8,931 shares of common stock issuable upon the exercise of warrants, BCIP III held 265,727 shares of common stock, which includes 1,272 shares of common stock issuable upon the exercise of warrants, and BCIP III-B held 5,348 shares of common stock, which includes 25 shares of common stock issuable upon the exercise of warrants. As a result of the foregoing, the Bain Capital Entities may collectively be deemed to beneficially own in the aggregate 2,136,085 shares of common stock.
- (3) Based on a Schedule 13G filed by bioMérieux SA ("bioMérieux") with the SEC on February 14, 2018, which reported ownership as of December 31, 2017. The board of directors of bioMérieux consists of Jean-Luc Belingard, Alexandre Mérieux, Alain Mérieux, Marie-Hélène Habert, Michele Palladino, Philippe Gillet, Agnès Lemarchand, Philippe Archinard, and Harold Boël, who have shared investment and voting control with respect to the shares held by bioMérieux and may exercise such control only with the support of the majority of the members of the board of directors of bioMérieux. No individual member of the board of directors of bioMérieux is deemed to hold any beneficial ownership or reportable pecuniary interest in the shares held by bioMérieux.
- (4) Based on a Schedule 13G filed by Flagship Ventures Fund 2004, L.P. ("Fund 2004"), Flagship Ventures General Partner LLC ("Fund 2004 GP"), Noubar B. Afeyan, Ph.D. ("Dr. Afeyan"), and Edwin M. Kania, Jr. ("Mr. Kania") with the SEC on February 14, 2018, which reported ownership as of December 31, 2017. As the general partner of Fund 2004, Fund 2004 GP may be deemed to beneficially own shares of common stock directly held by Fund 2004. Messrs. Afeyan and Kania, as Managers of Fund 2004 GP, may be deemed to beneficially own shares of common stock

beneficially owned by Fund 2004 GP. Douglas G. Cole, M.D., a member of our board of directors, is a managing partner of Flagship Pioneering and a member of the Flagship Ventures General Partner LLC ("GP"), the sole general partner of Fund 2004. Dr. Cole does not have either voting or investment control over Fund 2004's shares and he disclaims beneficial ownership of such shares, except to the extent of his pecuniary interest therein.

- (5) Consists of 951,180 shares of common stock (which includes 39,815 shares of restricted stock subject to performance-based vesting), 181,480 shares of common stock issuable upon the exercise of stock options exercisable within 60 days of May 15, 2019 and 13,822 shares of common stock issuable upon the vesting of RSUs within 60 days of May 15, 2019 held by Mr. Hrusovsky.
- (6) Consists of 2,270 shares of common stock and 7,900 shares of common stock issuable upon the exercise of stock options exercisable within 60 days of May 15, 2019 held by Dr. Cole. Dr. Cole is a managing partner of Flagship Pioneering but has no voting or investment power with respect to the securities described in footnote 4.
- (7) Consists of 4,970 shares of common stock and 7,900 shares of common stock issuable upon the exercise of stock options exercisable within 60 days of May 15, 2019 held by Mr. Connolly. Mr. Connolly is a senior advisor for Bain Capital Ventures but has no voting or investment power with respect to the securities described in footnote 2.
- (8) See footnote 1. Also includes 2,772 shares of common stock and 7,900 shares of common stock issuable upon the exercise of stock options exercisable within 60 days of May 15, 2019 held by Mr. Crandell. Mr. Crandell disclaims beneficial ownership of the securities described in footnote 1 except to the extent of his pecuniary interest therein, if any.
- (9) Consists of 119,648 shares of common stock and 18,268 shares of common stock issuable upon the exercise of stock options exercisable within 60 days of May 15, 2019 held by Dr. Dekkers.
- (10) Consists of 2,991 shares of common stock held by Dr. Madaus, 94,071 shares of common stock issuable upon the exercise of stock options exercisable within 60 days of May 15, 2019 held by Dr. Madaus and 224,966 shares of common stock held by a trust for which Dr. Madaus is a trust advisor who shares voting and investment power over the shares held by the trust.
- (11) Consists of 2,857 shares of common stock and 31,235 shares of common stock issuable upon the exercise of stock options exercisable within 60 days of May 15, 2019 held by Mr. Meister. Does not include 120,035 shares of common stock held by Woburn Abbey March 2009 Trust, U.D.T. March 31, 2009 as Mr. Meister does not have voting or investment power over the shares held by the trust.
- (12) Consists of 1,217,770 shares of common stock and 7,900 shares of common stock issuable upon the exercise of stock options exercisable within 60 days of May 15, 2019 held by Dr. Walt. Does not include 200,684 shares of common stock held by The David R. Walt 2008 Irrevocable Family Trust (the "Walt Trust"), of which his spouse is trustee and has sole voting and investment power over the shares held by the Walt Trust.
- (13) Consists of 3,290 shares of common stock, 21,800 shares of common stock issuable upon the exercise of stock options exercisable within 60 days of May 15, 2019, and 736 shares of common stock issuable upon the vesting of RSUs within 60 days of May 15, 2019 held by Dr. Mattoon.
- (14) See footnotes 5 through 13. Also includes an aggregate of 36,673 shares of common stock, 265,830 shares of common stock issuable upon the exercise of stock options exercisable within 60 days of May 15, 2019 and 525 shares of common stock issuable upon the vesting of RSUs within 60 days of May 15, 2019 held by David C. Duffy, Ph.D., our Senior Vice President, Research & Development and Chief Technology Officer, Mark T. Roskey, Ph.D., our Senior Vice President, Commercial and Accelerator, and Amol Chaubal, our Chief Financial Officer.

MANAGEMENT AND CORPORATE GOVERNANCE

The Board of Directors

Our restated certificate of incorporation and our restated bylaws provide that our business is to be managed by or under the direction of our board of directors. Our board of directors is divided into three classes for purposes of election. One class is elected at each annual meeting of stockholders to serve for a three-year term. Our board of directors currently consists of eight members, classified into three classes as follows: (1) Douglas G. Cole, M.D., Marijn Dekkers, Ph.D. and David R. Walt, Ph.D. constitute Class II with a term ending at the 2019 annual meeting; (2) Keith L. Crandell, E. Kevin Hrusovsky and Paul M. Meister constitute Class III with a term ending at the 2020 annual meeting; and (3) John M. Connolly and Martin D. Madaus, Ph.D. constitute Class I with a term ending at the 2021 annual meeting. Sarah E. Hlavinka has been nominated for election at the 2019 annual meeting, along with Dr. Dekkers and Dr. Walt, to serve as the Class II directors.

Our board of directors accepted the recommendation of the nominating and governance committee and voted to nominate Marijn Dekkers, Ph.D., Sarah E. Hlavinka and David R. Walt, Ph.D. for election at the annual meeting for a term of three years to serve until the 2022 annual meeting of stockholders, and until their respective successors have been elected and qualified.

Set forth below are the names of the persons nominated as directors and directors whose terms do not expire this year, their ages, their offices in the Company, if any, their principal occupations or employment for at least the past five years, the length of their tenure as directors and the names of other public companies in which such persons hold or have held directorships during the past five years, as of May 15, 2019. Additionally, information about the specific experience, qualifications, attributes or skills that led to our board of directors' conclusion at the time of filing of this proxy statement that each person listed below should serve as a director is set forth below:

<u>Name</u>	<u>Age</u>	<u>Position with the Company</u>
E. Kevin Hrusovsky	58	Chairman, President and Chief Executive Officer
John M. Connolly(1)(3)	67	Director
Keith L. Crandell(2)	59	Director
Marijn Dekkers, Ph.D.(3)	61	Director
Sarah E. Hlavinka	54	Director Nominee
Martin D. Madaus, Ph.D.(1)(3)	59	Director
Paul M. Meister(1)	66	Director
David R. Walt, Ph.D.(2)	66	Director

- (1) Member of our audit committee
- (2) Member of our compensation committee
- (3) Member of our nominating and governance committee

Our board of directors has reviewed the materiality of any relationship that each of our directors and director nominees has with Quanterix Corporation, either directly or indirectly. Based upon this review, our board of directors has determined that the following directors and director nominees are "independent directors" as defined by The Nasdaq Stock Market: Douglas G. Cole, M.D., John M. Connolly, Keith L. Crandell, Marijn Dekkers, Ph.D., Sarah E. Hlavinka, Martin D. Madaus, Ph.D., Paul M. Meister and David R. Walt, Ph.D.

E. Kevin Hrusovsky has been a member of our board of directors and our Chairman since June 2014 and our President and Chief Executive Officer since January 2015. Prior to joining us, Mr. Hrusovsky served as Senior Vice President of PerkinElmer, Inc., a publicly traded company that produces analytical instruments, genetic testing and diagnostic tools, medical imaging components,

software, instruments and consumables for multiple end markets, from February 2012 to May 2013 and served as President of the Life Sciences and Technology business unit of PerkinElmer, Inc. from November 2011 to May 2013. From May 2013 through September 2013, he served as a consultant to PerkinElmer. Previously, Mr. Hrusovsky served as Chief Executive Officer and President of Caliper Life Sciences, Inc., a life sciences company that developed and sold lab automation equipment, from July 2003 to November 2011 when it was acquired by PerkinElmer, Inc. Prior to that, he served as Chief Executive Officer and President of Zymark, a company that provided laboratory automation, robotics, and liquid handling solutions, and Director of International Business, Agricultural Chemical Division, and President of the Pharmaceutical Division for FMC Corporation, a publicly traded chemical manufacturing company. He also held several management positions at E.I. DuPont de Nemours, a publicly traded chemical manufacturing company. Mr. Hrusovsky serves on the boards of directors of several private companies, including BioreclamationIVT, LLC, Cellaria, Cell Signaling Technology, Inc., 908 Devices Inc., and Solect Energy Development, LLC, and is the founder of the Powering Precision Health Summit, an executive think tank of researchers, scientists, physicians and innovators focused on the latest research on new biomarkers that have the potential to advance personalized medicine. Mr. Hrusovsky also serves on the Educational Board of the Massachusetts Biotech Council, the Advisory Committee for the Center for Biomedical Engineering at Brown University, the Association for Laboratory Automation, the JALA Editorial Board and the Strategy Committee of Children's Hospital Boston. He formerly served on the boards of SynapDx Corp., SeraCare, Caliper Life Sciences, Xenogen—XGEN and Alliant Medical Technology. He was selected as the 2013 Entrepreneur of the Year from Ohio State University and holds an Honorary Doctorate degree from Framingham State University for contributions in life sciences and personal medicine. Mr. Hrusovsky has a B.S. in mechanical engineering from Ohio State University and an M.B.A. from Ohio University. Our board of directors has concluded that Mr. Hrusovsky possesses specific attributes that qualify him to serve as a member of our board of directors, including the perspective and experience he brings as our President and Chief Executive Officer, which provides our board with historic knowledge, operational expertise and continuity.

John M. Connolly has served as a member of our board of directors since December 2012. Since December 2015, Mr. Connolly has served as a senior advisor for Bain Capital Ventures, a venture capital firm he joined in 2009. Mr. Connolly was previously a managing director at Bain Capital Ventures from September 2009 to January 2015. Mr. Connolly also served as Interim Chief Executive Officer of Education Holdings 1, Inc. from May 2012 to September 2012, Interim President and Chief Executive Officer of The Princeton Review from March 2011 to April 2012, Chairman of the board of directors of EDGAR Online, Inc. from March 2011 to August 2012 and Interim President and Chief Executive Officer of EDGAR Online, Inc. from September 2010 to March 2011. Prior to that, Mr. Connolly served as President and Chief Executive Officer of M-C Communications, Inc., a leading provider of continuing medical education in the United States, from March 2007 to July 2009, and as Chairman of the board of directors of M-C Communications, Inc. from August 2009 to August 2010. Prior to that, from March 2004 to March 2007, Mr. Connolly served as President and Chief Executive Officer of Institutional Stockholder Services Inc., which provides proxy voting, corporate governance, compliance, and risk management solutions. Education Holdings 1, Inc. filed a voluntary chapter 11 bankruptcy petition in January 2013. Mr. Connolly has a B.A. from St. Norbert College and an Executive Education Degree from the Executive Education Program at INSEAD which is located in Fontainebleau, France. Our board of directors has concluded that Mr. Connolly possesses specific attributes that qualify him to serve as a member of our board of directors, including his substantial experience as an investor in emerging companies as well as his significant managerial and corporate governance experience.

Keith L. Crandell has served as a member of our board of directors since June 2007. Since July 1994, Mr. Crandell has served as a managing director of ARCH Venture Partners, a venture capital firm focused on early-stage technology companies. He serves on the board of directors of Adesto

Technologies Corporation, a publicly traded provider of application-specific and ultra-low power non-volatile memory products, and of Twist Bioscience Corporation, a publicly traded synthetic biology company. He is also a director of several private companies and he also serves as a director of the Illinois Venture Capital Association. Mr. Crandell has a B.S. degree in chemistry and mathematics from St. Lawrence University, an M.S. degree in chemistry from the University of Texas at Arlington and an M.B.A. from the University of Chicago. Our board of directors has concluded that Mr. Crandell possesses specific attributes that qualify him to serve as a member of our board of directors, including his financial expertise and his substantial experience as an investor in emerging companies.

Marijn Dekkers, Ph.D. has served as a member of our board of directors since March 2017. He is the founder and Chairman of Novalis LifeSciences LLC, an investment and advisory firm for the life science industry. Since April 2016, he has served as Chairman of the board of directors of Unilever. From 2010 to 2016, Dr. Dekkers was Chief Executive Officer of Bayer AG in Leverkusen, Germany. From 2000 to 2002, he was the Chief Operating Officer at Thermo Electron Corporation (later renamed Thermo Fisher Scientific Inc.). In 2002, he became Thermo Fisher's President and Chief Executive Officer, where he served until 2009. Dr. Dekkers began his career in 1985 as a research scientist at General Electric Company. He holds a degree in chemistry from Radboud University in Nijmegen, the Netherlands, and a Ph.D. in chemical engineering from the University of Eindhoven. Our board of directors has concluded that Dr. Dekkers possesses specific attributes that qualify him to serve as a member of our board of directors, including his substantial experience leading commercial stage healthcare companies and his significant corporate governance experience.

Sarah E. Hlavinka has been nominated by our board of directors for election at the annual meeting. Since August 2018, Ms. Hlavinka has served as Senior Vice President, General Counsel and Secretary of Itron, Inc., a publicly-traded technology and services company focused on critical infrastructure solutions. Prior to joining Itron, from January 2017 to July 2018, Ms. Hlavinka served as Executive Vice President, General Counsel and Secretary of Xerox Corporation, a publicly-traded document management systems and solutions company. Prior to joining Xerox Corporation, from September 2007 to January 2017, Ms. Hlavinka served as Executive Vice President, General Counsel and Secretary of ABM Industries Incorporated, a publicly-traded integrated facility services provider. Ms. Hlavinka holds a J.D. from the University of Texas School of Law and a B.A. in history from Texas A&M University. Our board of directors has concluded that Ms. Hlavinka possesses specific attributes that qualify her to serve as a member of our board of directors, including her significant experience as a senior executive of global companies operating in a variety of industries and her substantial expertise in corporate governance and risk management.

Martin D. Madaus, Ph.D. has served as a member of our board of directors since November 2010. Dr. Madaus previously served as our Executive Chairman from November 2010 to June 2014, as our Chief Executive Officer from October 2011 to July 2012 and as our President from June 2011 to July 2012. Since February 2019, he has served as Operating Executive to The Carlyle Group, a multinational private equity, alternative asset management and financial services corporation. From June 2014 to February 2019, Dr. Madaus served as Chairman and Chief Executive Officer at Ortho-Clinical Diagnostics, Inc., a diagnostics company that makes products and diagnostic equipment for blood testing. Previously, Dr. Madaus was the Chairman, President and Chief Executive Officer of Millipore Corporation (MIL), a life sciences company serving the bioscience research and biopharmaceutical manufacturing industry, from January 2005 to July 2010, when Millipore was acquired by Merck KGaA. From July 2009 to May 2015, Dr. Madaus served as a member of the board of directors of Mettler Toledo International, a manufacturer of scales and analytical instruments. Dr. Madaus received a Doctor of Veterinary Medicine from the University of Munich in Germany and a Ph.D. in Veterinary Medicine from the Veterinary School of Hanover in Germany. Dr. Madaus has extensive public and private company board experience and our board of directors has concluded that Dr. Madaus possesses

specific attributes that qualify him to serve as a member of our board of directors, including his substantial knowledge of and managerial experience in the diagnostics industry.

Paul M. Meister has served as a member of our board of directors since September 2013. Mr. Meister is Co-Founder, and since 2008, Chief Executive Officer, of Liberty Lane Partners, LLC, a private investment company with diverse investments in healthcare, technology and distribution-related industries, and is Vice Chairman and Co-Founder of Perspecta Trust, a New Hampshire-based trust company. Mr. Meister served as President of MacAndrews & Forbes from 2014 to 2018. Mr. Meister also served as Executive Vice Chairman of Revlon, Inc., a leading beauty products company, on an interim basis from January 2018 to November 2018, when the Chief Executive Officer of Revlon, Inc. resigned. Mr. Meister previously served as Chairman and Chief Executive Officer of inVentiv Health (now known as Syneos Health, Inc.), a provider of commercial, consulting and clinical research services to the pharmaceutical and biotech industries, from 2010 to 2015. Mr. Meister was Chairman of Thermo Fisher Scientific, Inc., a scientific instruments equipment and supplies company, from November 2006 to April 2007. He was previously Vice Chairman of Fisher Scientific International, Inc., a predecessor of Thermo Fisher Scientific, Inc., from 2001 to 2006, and Chief Financial Officer of Fisher Scientific International, Inc. from 1991 to 2001. Prior to joining Fisher Scientific International, Inc., Mr. Meister held executive positions with the Henley Group, Wheelabrator Technologies and Abex, Inc. Mr. Meister has served as a director of Scientific Games Corporation, which provides customized, end-to-end solutions to the gaming industry, since 2012. He also previously served as a director of LKQ Corporation, a distributor of vehicle products, from 1999 to 2018, and vTv Therapeutics, Inc., a clinical-stage biopharmaceutical company, from 2015 to 2018 and Revlon, Inc. from 2016 to 2019. Mr. Meister is Co-Chair of the University of Michigan's Life Sciences Institute External Advisory Board and Chair of the Provost's Advisory Committee. Mr. Meister has an M.B.A. from Northwestern University and a B.A. from the University of Michigan. Our board of directors has concluded that Mr. Meister possesses specific attributes that qualify him to serve as a member of our board of directors, including his financial and investment expertise and his extensive knowledge of the life sciences industry.

David R. Walt, Ph.D. has served as a member of our board of directors since April 2007. Dr. Walt was our founding scientist and serves as chairman of our scientific advisory board. Dr. Walt currently serves as Hansjörg Wyss Professor of Biologically Inspired Engineering and Professor of Pathology at Harvard Medical School in the Department of Pathology at the Brigham and Women's Hospital and is a core faculty member of the Wyss Institute for Biologically Inspired Engineering. He is also a Howard Hughes Medical Institute Professor. He previously served as University Professor, Professor of Chemistry, Professor of Biomedical Engineering, Professor of Genetics, Professor of Neuroscience, Professor of Cell and Molecular Biology, and Professor of Oral Medicine at Tufts University, from 1981-2017. Dr. Walt was also the founding scientist of Illumina, Inc. and served as a member of the board of its directors from 1998-2016. He served on the board of directors of Cerulean Pharma Inc. (which was acquired by Daré Bioscience, Inc.) from August 2016 to July 2017. Dr. Walt was also a founder and currently serves as a member of the board of directors of Ultivue, Inc., Arbor Biotechnologies, Inc. and Sherlock Biosciences, Inc. and serves as a member of the board of directors of Exicure, Inc. He has received numerous national and international awards and honors for his fundamental and applied work in the field of optical sensors, microwell arrays and single molecule detection. He is a member of the U.S. National Academy of Engineering, U.S. National Academy of Medicine, American Academy of Arts and Sciences, a fellow of the American Institute for Medical and Biological Engineering, a fellow of the National Academy of Inventors, and a fellow of the American Association for the Advancement of Science. Dr. Walt has a B.S. in Chemistry from the University of Michigan and a Ph.D. in Chemical Biology from Stony Brook University. Our board of directors has concluded that Dr. Walt possesses specific attributes that qualify him to serve as a member of our board of directors, including his experience in developing life sciences companies and his expertise in chemistry, diagnostics technologies and biomedical engineering.

There are no family relationships between or among any of our directors or director nominees. The principal occupation and employment during the past five years of each of our directors and director nominees was carried on, in each case except as specifically identified above, with a corporation or organization that is not a parent, subsidiary or other affiliate of us. There is no arrangement or understanding between any of our directors or nominees and any other person or persons pursuant to which he or she is to be selected as a director or nominee.

There are no legal proceedings to which any of our directors is a party adverse to us or any of our subsidiaries or in which any such person has a material interest adverse to us or any of our subsidiaries.

Committees of the Board of Directors and Meetings

Meeting Attendance. During the fiscal year ended December 31, 2018, our board of directors met four times and acted by written consent on four occasions. During the fiscal year ended December 31, 2018, the various committees of the board of directors met a total of six times and acted by written consent on five occasions. No director attended fewer than 75% of the total number of meetings of the board and of committees of the board on which such director served during fiscal 2018. Our board of directors has adopted a policy under which each member of the board is encouraged but not required to attend each annual meeting of our stockholders. The directors did not attend our 2018 annual meeting of stockholders.

Audit Committee. Our audit committee met four times and acted by written consent on one occasion during the fiscal year ended December 31, 2018. This committee currently has three members, Paul M. Meister (Chairman), John M. Connolly and Martin D. Madaus, Ph.D. Our audit committee's role and responsibilities are set forth in the audit committee's written charter and include:

- selecting a firm to serve as the independent registered public accounting firm to audit our financial statements;
- ensuring the independence of the independent registered public accounting firm;
- discussing the scope and results of the audit with the independent registered public accounting firm, and reviewing, with management and that firm, our interim and year-end operating results;
- establishing procedures for employees to anonymously submit concerns about questionable accounting or audit matters;
- considering the effectiveness of our internal controls and, if any, our internal audit function;
- reviewing material related-party transactions or those that require disclosure; and
- approving or, as permitted, pre-approving all audit and non-audit services to be performed by the independent registered public accounting firm.

Except as noted below, all members of the audit committee satisfy the current independence standards promulgated by the Securities and Exchange Commission, or the SEC, and by The Nasdaq Stock Market, as such standards apply specifically to members of audit committees. Our board of directors has determined that Paul M. Meister is an "audit committee financial expert," as the SEC has defined that term in Item 407 of Regulation S-K.

A copy of the audit committee's written charter is publicly available on our website at www.quanterix.com.

Compensation Committee. Our compensation committee met two times and acted by written consent on three occasions during the fiscal year ended December 31, 2018. This committee currently has three members, Keith L. Crandell (Chairman), Douglas G. Cole, M.D. and David R. Walt, Ph.D.

Our compensation committee's role and responsibilities are set forth in the compensation committee's written charter and includes:

- reviewing and approving, or recommending that our board of directors approve, the compensation of our executive officers;
- reviewing and recommending to our board of directors the compensation of our directors;
- reviewing and recommending to our board of directors the terms of any compensatory agreements with our executive officers;
- administering our 2017 Employee, Director and Consultant Equity Incentive Plan, 2007 Stock Option and Grant Plan, as amended, and 2017 Employee Stock Purchase Plan;
- reviewing and approving, or making recommendations to our board of directors with respect to, incentive compensation and equity plans; and
- reviewing all overall compensation policies and practices.

The compensation committee is also responsible for the determination of the compensation of our chief executive officer, and will conduct its decision making process with respect to that issue without the chief executive officer present. Our board of directors has determined that all members of the compensation committee qualify as "independent" under the definition promulgated by The Nasdaq Stock Market.

The compensation committee has the authority to retain or obtain the advice of such compensation consultants, legal counsel, experts and other advisors as it may deem appropriate in its sole discretion. The compensation committee is directly responsible for the appointment, compensation and oversight of its consultants, legal counsel, experts and advisors and has sole authority to approve their fees and retention terms, and the Company will provide funding for such fees and related expenses. Before selecting any such consultant, legal counsel, expert or advisor, the compensation committee must consider the following independence factors:

- The provision of other services to the Company by the entity that employs the consultant, legal counsel, expert or advisor.
- The amount of fees received from the Company by the entity that employs the consultant, legal counsel, expert or advisor as a percentage of the total revenue of such entity.
- The policies and procedures of the entity that employs the consultant, legal counsel, expert or advisor that are designed to prevent conflicts of interest.
- Any business or personal relationship of the consultant, legal counsel, expert or advisor with a member of the compensation committee or an executive officer of the company.
- Any of our stock owned by the consultant, legal counsel, expert or advisor.
- Any business or personal relationship the consultant, legal counsel, expert or advisor has with an executive officer of the company.

The compensation committee engaged Pay Governance as an independent advisor to the compensation committee providing executive compensation consulting services for the fiscal year ended December 31, 2018 and for the fiscal year ending December 31, 2019. Pay Governance was engaged by and reports to the compensation committee. Pay Governance is not expected to provide any services to the Company other than executive compensation consulting services during the fiscal year ending December 31, 2019. In compliance with the SEC and the corporate governance rules of The Nasdaq Stock Market, Pay Governance has provided the compensation committee with a letter addressing each of the six independence factors noted above. Their responses affirm the independence of Pay

Governance and the partners, consultants, and employees who service the compensation committee on executive compensation matters and governance issues.

A copy of the compensation committee's written charter is publicly available on our website at www.quanterix.com.

Nominating and Governance Committee. Our nominating and governance committee, or "nominating committee", acted by written consent on one occasion during the fiscal year ended December 31, 2018 and has three members, Marijn Dekkers, Ph.D. (Chairman), John M. Connolly and Martin D. Madaus, Ph.D. Our board of directors has determined that all members of the nominating committee qualify as independent under the definition promulgated by The Nasdaq Stock Market. The nominating committee's responsibilities are set forth in the nominating committee's written charter and include:

- identifying and recommending candidates for membership on our board of directors;
- recommending directors to serve on board committees;
- reviewing and recommending our corporate governance guidelines and policies;
- reviewing proposed waivers of the code of conduct for directors and executive officers;
- evaluating, and overseeing the process of evaluating, the performance of our board of directors and individual directors; and
- assisting our board of directors on corporate governance matters.

Generally, our nominating committee considers candidates recommended by stockholders as well as from other sources such as other directors or officers, third party search firms or other appropriate sources. Once identified, the nominating committee will evaluate a candidate's qualifications in accordance with our Nominating and Governance Committee Policy Regarding Qualifications of Directors appended to our nominating committee's written charter. Threshold criteria include: personal integrity and sound judgment, business and professional skills and experience, independence, knowledge of our industry, possible conflicts of interest, diversity, the extent to which the candidate would fill a present need on the board of directors, and concern for the long-term interests of our stockholders. Our nominating committee has not adopted a formal diversity policy in connection with the consideration of director nominations or the selection of nominees. However, the nominating committee will consider issues of diversity among its members in identifying and considering nominees for director, and strive where appropriate to achieve a diverse balance of backgrounds, perspectives, experience, age, gender, ethnicity and country of citizenship on the board and its committees.

If a stockholder wishes to propose a candidate for consideration as a nominee for election to the Board, it must follow the procedures described in our restated bylaws and in "Stockholder Proposals and Nominations For Director" at the end of this proxy statement. In general, persons recommended by stockholders will be considered in accordance with our Policy on Shareholder Recommendation of Candidates for Election as Directors appended to our nominating committee's written charter. Any such recommendation should be made in writing to the nominating committee, care of our Corporate Secretary at our principal office and should be accompanied by the following information concerning each recommending stockholder and the beneficial owner, if any, on whose behalf the nomination is made:

- all information relating to such person that would be required to be disclosed in a proxy statement;
- certain biographical and share ownership information about the stockholder and any other proponent, including a description of any derivative transactions in our securities;

- a description of certain arrangements and understandings between the proposing stockholder and any beneficial owner and any other person in connection with such stockholder nomination; and
- a statement whether or not either such stockholder or beneficial owner intends to deliver a proxy statement and form of proxy to holders of voting shares sufficient to carry the proposal.

The recommendation must also be accompanied by the following information concerning the proposed nominee:

- certain biographical information concerning the proposed nominee;
- all information concerning the proposed nominee required to be disclosed in solicitations of proxies for election of directors;
- certain information about any other security holder of the Company who supports the proposed nominee;
- a description of all relationships between the proposed nominee and the recommending stockholder or any beneficial owner, including any agreements or understandings regarding the nomination; and
- additional disclosures relating to stockholder nominees for directors, including completed questionnaires and disclosures required by our restated bylaws.

A copy of the nominating committee's written charter, including its appendices, is publicly available on our website at www.quanterix.com.

Compensation Committee Interlocks and Insider Participation

Our compensation committee has three members, Keith L. Crandell (Chairman), Douglas G. Cole, M.D. and David R. Walt, Ph.D. None of the members of our compensation committee has at any time during the prior three years been one of our officers or employees. None of our executive officers currently serves, or in the past fiscal year has served, as a member of the board of directors or compensation committee of any entity that has one or more executive officers serving on our board of directors or compensation committee. For a description of transactions between us and members of our compensation committee and affiliates of such members, please see "Certain Relationships and Related Person Transactions."

Board Leadership Structure and Role in Risk Oversight

Our board of directors does not have a policy regarding the separation of the roles of Chief Executive Officer and Chairman of the board of directors, as our board of directors believes it is in the best interest of the Company to make that determination based on the position and direction of the Company and the membership of the board of directors. Our board of directors has determined that having an employee director serve as Chairman is in the best interest of our stockholders at this time because of the efficiencies achieved in having the role of Chief Executive Officer and Chairman combined, and because the detailed knowledge of our day-to-day operations and business that the Chief Executive Officer possesses greatly enhances the decision-making processes of our board of directors as a whole. We have a strong governance structure in place, including independent directors, to ensure the powers and duties of the dual role are handled responsibly. Martin D. Madaus, Ph.D. serves as our lead independent director.

The Chairman of the board of directors, the lead independent director and the other members of the board of directors work in concert to provide oversight of our management and affairs. Our board of directors encourages communication among its members and between management and the board of

directors to facilitate productive working relationships. Working with the other members of the board of directors, our Chairman also strives to ensure that there is an appropriate balance and focus among key board responsibilities such as strategic development, review of operations and risk oversight.

Risk is inherent with every business, and how well a business manages risk can ultimately determine its success. Our board of directors is actively involved in oversight of risks that could affect the Company. This oversight is conducted primarily by our full board of directors, which has responsibility for general oversight of risks. For example, in determining whether and under what circumstances we will engage in financing transactions or enter into licensing, collaboration or similar arrangements, the board of directors is involved in our management of risks related to our financial condition or of the risks inherent in product development and commercialization.

As part of its oversight, our board of directors receives reports by each committee chair regarding the committee's considerations and actions. In particular, the audit committee is responsible for discussing the adequacy of our risk management activities with management and our independent registered public accounting firm. The audit committee's primary emphasis is financial risk. In addition, the compensation committee is responsible for considering whether our compensation programs and practices are reasonably likely to have a material adverse effect on us.

At each of its meetings, the board of directors receives business updates from various members of management. These updates may identify matters that have emerged within that member of management's scope of responsibility that involve operational, financial, legal or regulatory risks and, in these cases, the board of directors provides guidance to management. Our board of directors believes that full and open communication between management and the board of directors is essential for effective risk management and oversight.

Stockholder Communications to the Board

Generally, stockholders who have questions or concerns should contact our Investor Relations department at ir@quanterix.com. However, any stockholders who wish to address questions regarding our business directly with our board of directors, or any individual director, should direct his or her questions in writing to the Chairman of the Board at 900 Middlesex Turnpike, Building 1, Billerica, Massachusetts 01821 or using the "IR Contacts" page of our website at <http://ir.quanterix.com/contact-ir>. Communications will be distributed to the board of directors, or to any individual director or directors as appropriate, depending on the facts and circumstances outlined in the communications. Items that are unrelated to the duties and responsibilities of the board may be excluded, such as:

- junk mail and mass mailings;
- resumes and other forms of job inquiries;
- surveys; and
- solicitations or advertisements.

In addition, any material that is unduly hostile, threatening, or illegal in nature may be excluded, in which case it will be made available to any outside director upon request.

Executive Officers

The following table sets forth certain information regarding our executive officers who are not also directors, as of May 15, 2019. We have entered into agreements with our executive officers, and all of them are at-will employees.

<u>Name</u>	<u>Age</u>	<u>Position(s) with the Company</u>
Amol Chaubal	43	Chief Financial Officer
David C. Duffy, Ph.D.	48	Senior Vice President, Research & Development and Chief Technology Officer
Dawn R. Mattoon, Ph.D.	46	Senior Vice President, Research Products
Mark T. Roskey, Ph.D.	59	Senior Vice President, Commercial and Accelerator
Jackson Streeter, M.D.	57	Senior Vice President, Corporate Development and Strategy

Amol Chaubal has been our Chief Financial Officer since April 2019. Prior to joining us, Mr. Chaubal served as the Chief Financial Officer of Global Operations at Smith & Nephew plc, a global medical technology company, from October 2017 to April 2019. Prior to his time at Smith & Nephew plc, he served as Corporate Vice President and Head of Finance for the Clinical Research Services business segment at PAREXEL International Corporation from June 2015 to October 2017. Between May 2013 and May 2015, Mr. Chaubal served as CFO North America and later Regional Chief Operating Officer at Brookfield Renewable Energy Group, which is part of Brookfield Asset Management Inc. Prior to joining Brookfield, Mr. Chaubal was CFO North America at Novartis AG, where he spent 11 years in positions of increasing responsibility across Switzerland, Canada, and the United States. He has an M.B.A. (with distinction) from INSEAD (France). He began his career at Procter & Gamble Company after completing Masters degree in Chemical Engineering from the Indian Institute of Technology, Mumbai.

David C. Duffy, Ph.D. is Senior Vice President, Research & Development, and has been our Chief Technology Officer since March 2013. Prior to this role, Dr. Duffy served as our Vice President, Research from November 2011 to February 2013, and Senior Director, Platform Research from July 2007 to October 2011. Prior to joining us, Dr. Duffy served as Director, Pharmacomer Technology at Surface Logix, Inc., a biomedical development company. Prior to that, Dr. Duffy was a Senior Scientist at Gamera Bioscience Corporation, a developer of automated biotechnology test equipment, from December 1998 to January 2000. Previously, Dr. Duffy was a postdoctoral research fellow in the Department of Chemistry and Chemical Biology at Harvard University. He was the first Sir Alan Wilson Research Fellow of Emmanuel College, University of Cambridge. Dr. Duffy has a Ph.D. in physical chemistry from the Department of Chemistry at the University of Cambridge, and B.A. and M.A. degrees in Natural Sciences from the University of Cambridge.

Dawn R. Mattoon, Ph.D., has been our Senior Vice President, Research Products since March 2019 and served as Senior Vice President of Strategic Marketing and Assay Technologies from February 2018 to March 2019. Dr. Mattoon completed both her Ph.D. and postdoctoral work at Yale University, studying tyrosine kinase signaling pathways. Following her postdoctoral fellowship in 2004, she joined Invitrogen, where she was recognized as one of the company's best new scientists in 2005. Dr. Mattoon went on to serve in a variety of leadership roles in R&D, Program & Portfolio Management, Strategy, and General Management as Invitrogen evolved to become Life Technologies with the acquisition of Applied Biosystems in 2008, and was subsequently acquired by Thermo Fisher in 2013. Until her departure from Thermo Fisher in 2014, Dr. Mattoon also served on global leadership teams focused on mentoring and developing female leaders. Dr. Mattoon served as the Vice President for Product Development at Cell Signaling Technologies, a leading antibody provider, from July 2014 to February 2018, when she transitioned to her current role with Quanterix.

Mark T. Roskey, Ph.D., has been our Senior Vice President, Commercial and Accelerator since March 2019 and served as Senior Vice President and General Manager of Applications and Reagents from September 2014 to March 2019. Prior to joining us, Dr. Roskey served as Vice President and General Manager of the America's Sales and Service organization in PerkinElmer's Life Science and Technologies Division from December 2011 to September 2014. PerkinElmer is a global life science company. Prior to PerkinElmer, Dr. Roskey served as Senior Vice President of Biology Research and Development at Caliper Life Sciences, Inc. from March 2009 to November 2011, when it was acquired by PerkinElmer. At Caliper Life Sciences, Inc., Dr. Roskey served as Senior Vice President of Applied Biology Research from 2008 to 2011 and Vice President of Worldwide Marketing from July 2003 to 2008. Prior to that, Dr. Roskey served as Vice President of Worldwide Marketing for Zymark Corporation, a laboratory automation company, from December 2001 to August 2003 when it was acquired by Caliper Life Sciences, Inc. Prior to that, Dr. Roskey served as Director of Marketing at Applied Biosystems, Inc. Dr. Roskey completed a postdoctoral fellowship in molecular immunobiology at the Harvard Medical School and has a Ph.D. in microbiology from the University of Notre Dame and a B.S. in biology from Framingham State University.

Jackson Streeter, M.D. has been our Senior Vice President, Corporate Development and Strategy, since November 2018. From March 2016 to November 2018, Dr. Streeter was the CEO and Executive Director of the Florida Technology Seed Capital Fund, a seed stage venture fund investing in companies commercializing technologies developed from State of Florida universities, colleges or research institutes. Since inception, the fund has created over 50 companies with private investment matching Florida state funding at a 9:1 ratio. From January 2010 to March 2016, Dr. Streeter served as the CEO and Chief Medical Officer of Banyan Biomarkers. During his time the company received over \$110 million in contracts and grants with the NIH and Department of Defense and raised \$20 million in private capital for the development of a first ever blood test for traumatic brain injury, which received FDA clearance in February 2018. Prior to serving as CEO of Banyan Biomarkers, Dr. Streeter founded and served as CEO for PhotoThera, Inc. where he raised over \$120 million in private capital and was responsible for the design and management of five U.S. and international clinical trials for the treatment of stroke. He was also the co-founder of American Veterinary Laser, which was acquired in a private transaction. Dr. Streeter is the inventor of over twenty patents and author on multiple scientific publications. Dr. Streeter has served on the following boards and executive committees: Florida Research Consortium, Governor's Life Science Task Force for the State of Florida, FBI counter terrorism south east working group, Walter H. Coulter Center Committee at the University of Miami, Department of Defense Grey Team in Afghanistan, Sid Martin Biotechnology Incubator advisory board and the Gainesville Chamber of Commerce Board of Directors. Dr. Streeter served as an Officer in the U.S. Navy, completing a general surgery internship at Naval Regional Medical Center Portsmouth, Virginia and then primary Flight School at Naval Air Station Pensacola, Florida. During his military service, he was the first Naval Flight Surgeon selected as a TOPGUN staff instructor, served on several overseas deployments and was awarded the Navy Achievement Medal and the Navy Commendation Medal. He earned his B.S. degree in Biology from the University of Nevada Reno and his M.D. from the University of Nevada School of Medicine.

EXECUTIVE OFFICER AND DIRECTOR COMPENSATION

Summary Compensation Table

The following table shows the total compensation paid or accrued during the fiscal years ended December 31, 2018 and December 31, 2017, to our Chief Executive Officer and President and our two next most highly compensated executive officers who earned more than \$100,000 during the fiscal year ended December 31, 2018 and were serving as executive officers as of such date.

Name and Principal Position	Year	Salary (\$)	Bonus \$(1)	Stock Awards \$(2)	Option Awards \$(2)	All Other Compensation \$(3)	Total (\$)
E. Kevin Hrusovsky	2018	497,857	609,377	4,440,000	—	346	5,547,580
<i>Chairman, President and Chief Executive Officer</i>	2017	426,978	276,682		1,430,000	—	2,133,660
Dawn R. Mattoon, Ph.D.(4)	2018	239,077	133,986	368,042	558,215	3,102	1,302,422
<i>Senior Vice President, Research Products</i>							
Jackson Streeter, M.D.(5)	2018	41,538	20,769	464,442	409,357	1,246	937,352
<i>Senior Vice President, Corporate Development and Strategy</i>							

- (1) Amounts for 2018 represent cash bonuses earned for the 12-month period from January 1, 2018 to December 31, 2018, and exclude payments made in 2018 for 2017 bonuses. Amounts for 2017 represent cash bonuses earned for the 12-month period from January 1, 2017 to December 31, 2017, and exclude payments made in 2017 for 2016 bonuses.
- (2) These amounts represent the aggregate grant date fair value for restricted stock unit awards (RSUs) and option awards granted during our fiscal year ended December 31, 2018, computed in accordance with FASB ASC Topic 718. A discussion of the assumptions used in determining grant date fair value for the option awards may be found in Note 2 to our consolidated financial statements for the year ended December 31, 2018.
- (3) Amounts represent the dollar value of matching contributions to the qualified 401(k) plan.
- (4) Dr. Mattoon joined Quanterix on February 23, 2018.
- (5) Dr. Streeter joined Quanterix on November 12, 2018.

Narrative Disclosure to Summary Compensation Table

E. Kevin Hrusovsky

We entered into an agreement with Mr. Hrusovsky with respect to his service as President and Chief Executive Officer on January 1, 2015. Under the terms of the agreement, Mr. Hrusovsky received a one-time signing bonus of \$250,000 and was entitled to an initial annual base salary of \$400,000. His current base salary is \$536,000. Under the agreement, Mr. Hrusovsky is eligible to receive an annual bonus of up to 85% of his base salary based on the compensation committee's assessment of his and the Company's performance against goals established by the committee. For 2018, our compensation committee awarded Mr. Hrusovsky a discretionary bonus of \$609,377. In connection with his employment, in February 2015 Mr. Hrusovsky was granted 781,060 shares of restricted stock. The shares underlying this grant vest, subject to continued service, as follows: (1) with respect to 168,246 of the shares, 25% of the shares vested on July 1, 2015, with the remainder vesting over the next three years in equal monthly installments on the first day of each succeeding calendar month thereafter, (2) with respect to 549,113 of the shares, 25% of the shares vested on September 1, 2015, with the

remainder vesting over the next three years in equal monthly installments on the first day of each succeeding calendar month thereafter and (3) 63,701 shares vest based upon the achievement of certain milestones.

Pursuant to the agreement, Mr. Hrusovsky is entitled to continuation of his then-current base salary and health insurance benefits for six months in the event we terminate his employment without Cause, as defined in the agreement, or Mr. Hrusovsky terminates his employment with us for Good Reason, as defined in the agreement, subject to Mr. Hrusovsky's execution of a release satisfactory to us following such termination. If such termination occurs within 12 months of the sale of the Company, he is entitled to continuation of his then-current base salary and health insurance benefits for 12 months. If we terminate his employment without Cause, his employment terminates due to his death or disability or he terminates his employment with us for Good Reason within three months prior to the end of a year, he is entitled to a pro rata portion of his bonus. In addition, Mr. Hrusovsky has entered into an employee non-competition, non-solicitation, confidentiality and assignment agreement pursuant to which he has agreed (1) not to engage in any competitive business for six months following his termination of employment with us (12 months if the termination is within 12 months of the sale of the Company), (2) not to solicit our employees, customers or suppliers for six months following his termination of employment with us (12 months if the termination is within 12 months of the sale of the Company) and (3) to assign any inventions conceived or developed during the course of his employment with us.

Dawn R. Mattoon, Ph.D.

We entered into an agreement with Dr. Mattoon with respect to her service as Senior Vice President, Strategic Marketing & Assay Operations on February 1, 2018. In March 2019, her title was changed to Senior Vice President, Research Products. Under the terms of the agreement, Dr. Mattoon was entitled to an initial annual base salary of \$280,000. Her current base salary is \$320,000. Dr. Mattoon is also eligible to receive an annual bonus of up to 40% of her then-current base salary based on her performance and our business conditions in the sole discretion of our board of directors. Pursuant to the agreement, Dr. Mattoon was granted an option to purchase up to 40,000 shares of our common stock at an exercise price of \$22.12 per share, which was the fair market value on the date she began her employment with us (February 23, 2018). She was also granted 10,000 RSUs. The option and RSUs vested as to 25% of the shares on February 23, 2019, with the remainder vesting over the subsequent three years in equal monthly installments. In addition, Dr. Mattoon has entered into an employee non-solicitation, confidentiality and assignment agreement pursuant to which she has agreed (1) not to solicit our employees, customers or suppliers for one year following voluntary termination of her employment and (2) to assign any inventions conceived or developed during the course of her employment with us.

Jackson Streeter, M.D.

We entered into an agreement with Dr. Streeter with respect to his service as Senior Vice President, Corporate Development and Strategy on October 22, 2018. Under the terms of the agreement, Dr. Streeter was entitled to an initial annual base salary of \$300,000. Dr. Streeter is also eligible to receive an annual bonus of up to 50% of his then-current base salary based on his performance and our business conditions in the sole discretion of our board of directors. Pursuant to the agreement, Dr. Streeter was granted an option to purchase up to 66,143 shares of our common stock at an exercise price of \$16.33 per share, which was the fair market value on the date he began his employment with us (November 12, 2018). He was also granted 28,441 RSUs. The option and RSUs will vest as to 25% of the shares on November 12, 2019, with the remainder vesting over the subsequent three years in equal monthly installments. In addition, Dr. Streeter has entered into an employee non-competition, non-solicitation, confidentiality and assignment agreement pursuant to

which he has agreed (1) not to engage in any business with certain competitors for one year following his voluntary termination of employment with us, (2) not to solicit our employees, customers or suppliers for one year following voluntary termination of his employment and (3) to assign any inventions conceived or developed during the course of his employment with us.

Outstanding Equity Awards at 2018 Fiscal Year-End

The following table shows grants of stock options and grants of unvested stock awards outstanding on the last day of the fiscal year ended December 31, 2018, including both awards subject to performance conditions and non-performance-based awards, to each of the executive officers named in the Summary Compensation Table.

Name	Option Awards(1)				Stock Awards(1)			
	Number of Securities Underlying Unexercised Options (#) Exercisable	Number of Securities Underlying Unexercised Options (#) Unexercisable	Option Exercise Price (\$)	Option Expiration Date	Number of Shares or Units of Stock That Have Not Vested (#)	Market Value of Shares or Units of Stock That Have Not Vested (\$)(2)	Equity Incentive Plan Awards: Number of Unearned Shares, Units or Other Rights That Have Not Vested (#)	Equity Incentive Plan Awards: Market or Payout Value of Unearned Shares, Units or Other Rights That Have Not Vested (\$)(2)
E. Kevin Hrusovsky <i>Chairman, President and Chief Executive Officer</i>	142,594(4)	168,544(4)	\$ 8.16	3/31/2027	—	—	39,815(3)	729,013
	—	—	—	—	225,000(5)	4,119,750	—	—
Dawn R. Mattoon, Ph.D. <i>Senior Vice President, Research Products</i>	—	40,000(6)	\$ 22.12	2/23/2028	—	—	—	—
	—	26,181(7)	\$ 19.12	5/10/2028	10,000(6)	183,100	—	—
	—	—	—	—	7,680(7)	140,621	—	—
Jackson Streeter, M.D. <i>Senior Vice President, Corporate Development and Strategy</i>	—	66,143(8)	\$ 16.33	11/12/2028	—	—	—	—
	—	—	—	—	28,441(8)	520,755	—	—

- (1) Each of the outstanding equity awards in the table above was granted pursuant to our 2007 Stock Option and Grant Plan, as amended, or our 2017 Employee, Director and Consultant Equity Incentive Plan.
- (2) The market value of the stock awards is determined by multiplying the number of shares by \$18.31, the closing price of our common stock on the Nasdaq Global Market on December 31, 2018, the last day of our fiscal year.
- (3) On February 13, 2015, Mr. Hrusovsky was granted 781,060 shares of restricted stock in connection with his appointment as our President and Chief Executive Officer. The shares underlying this grant vest, subject to continued service, as follows: (i) with respect to 168,246 of the shares, 25% of the shares vested on July 1, 2015, with the remainder vesting over the next three years in equal monthly installments on the first day of each succeeding calendar month thereafter, (ii) with respect to 549,113 of the shares, 25% of the shares vested on September 1, 2015, with the remainder vesting over the next three years in equal monthly installments on the first day of each succeeding calendar month thereafter and (iii) 63,701 shares vest based upon the achievement of certain milestones. Upon the sale of the Company, the vesting is accelerated such that no more than 25% of the shares shall remain unvested. If Mr. Hrusovsky is terminated without Cause (as defined in his employment agreement) or he terminates his service for Good Reason (as defined in his employment agreement) within 12 months following a sale of the Company, all of the unvested shares will become vested.
- (4) Represents an option to purchase shares of our common stock granted on March 31, 2017. The shares underlying this option vest, subject to continued service, as follows: 25% of the shares vested on March 31, 2018, with the remainder vesting over the next three years in equal monthly installments on the last day of each succeeding calendar month thereafter.
- (5) On August 7, 2018, Mr. Hrusovsky was granted 300,000 RSUs. The RSUs vest upon continued service and vested as to 50,000 shares on August 31, 2018, and vest with respect to 6,250 shares at the end of each month thereafter for 40 months.

- (6) On February 23, 2018, Dr. Mattoon was granted (i) an option to purchase 40,000 shares of our common stock and (ii) 10,000 RSUs. The shares underlying the option and RSUs vest, subject to continued service, as follows: 25% of the shares vested on February 23, 2019, with the remainder vesting over the next three years in equal monthly installments on the last day of each calendar month thereafter.
- (7) On May 10, 2018, Dr. Mattoon was granted (i) an option to purchase 26,181 shares of our common stock and (ii) 7,680 RSUs. The shares underlying the option and RSUs vest, subject to continued service, as follows: 25% of the shares vested on May 10, 2019, with the remainder vesting over the next three years in equal monthly installments on the last day of each calendar month thereafter.
- (8) On November 12, 2018, Dr. Streeter was granted (i) an option to purchase 66,143 shares of our common stock and (ii) 28,441 RSUs. The shares underlying the option and RSUs vest, subject to continued service, as follows: 25% of the shares vest on November 12, 2019, with the remainder vesting over the next three years in equal monthly installments on the last day of each calendar month thereafter.

Director Compensation

The following table shows the total compensation paid or accrued during the fiscal year ended December 31, 2018 to each of our non-employee directors. Directors who are employed by us are not compensated for their service on our board of directors.

Name	Fees Earned or Paid in Cash (\$)	Stock Awards (\$)(1)	Option Awards (\$)(1)	All Other Compensation (\$)	Total (\$)
Douglas G. Cole, M.D.	41,000	33,596	40,062	—	114,658
John M. Connolly	40,000	33,596	40,062	—	113,658
Keith L. Crandell	54,500	33,596	40,062	—	128,158
Marijn Dekkers, Ph.D.	45,000	33,596	40,062	—	118,658
Martin D. Madaus, Ph.D.	57,500	33,596	40,062	—	131,158
Paul M. Meister	55,000	33,596	40,062	—	128,658
David R. Walt, Ph.D.	41,000	33,596	40,062	—	114,658

- (1) These amounts represent the aggregate grant date fair value for RSUs and option awards granted to each director in the fiscal year ended December 31, 2018, computed in accordance with FASB ASC Topic 718. A discussion of the assumptions used in determining grant date fair value for the option awards may be found in Note 2 to our consolidated financial statements for the year ended December 31, 2018. The following table shows the aggregate number of stock options held by each of our non-employee directors as of December 31, 2018.

Name	Aggregate Number of Shares Subject to Stock Options
Douglas G. Cole, M.D.	7,900
John M. Connolly	7,900
Keith L. Crandell	7,900
Marijn Dekkers, Ph.D.	39,013
Martin D. Madaus, Ph.D.	94,071
Paul M. Meister	31,235
David R. Walt, Ph.D.	7,900

Our current non-employee director compensation policy provides for the annual payment of: \$35,000 to each of our nonemployee directors, or \$55,000 in the case of the chair; \$7,500 to each member of our audit committee, or \$20,000 in the case of the chair; \$6,000 to each member of our

compensation committee, or \$12,000 in the case of the chair; and \$5,000 to each member of our nominating and governance committee, or \$10,000 in the case of the chair; in each case quarterly in arrears. Each non-employee director may elect to receive this annual payment in the form of common stock in lieu of cash.

Our current non-employee director compensation policy also provides that each non-employee director be granted, on the first trading day of each fiscal year, (1) a non-qualified stock option exercisable for 7,900 shares of common stock and (2) 2,270 restricted stock units (each restricted stock unit relating to one share of common stock). The annual stock options and restricted stock units vest in full on December 31 of the year in which such awards were granted, provided that the non-employee director is a director of the Company on the applicable vesting date.

In addition, each new non-employee director receives under the policy (1) a non-qualified stock option grant exercisable for 15,800 shares of common stock and (2) 4,540 restricted stock units (each restricted stock unit relating to one share of common stock) upon his or her initial election or appointment. The initial stock options vest over three years from the date of grant, with one third vesting on the first anniversary of the applicable grant date and the remainder vesting over the following two years in 24 successive equal monthly installments at the end of each month until the third anniversary of such grant date, provided that the non-employee director is a director of the Company on the applicable vesting date. The restricted stock units granted to each new non-employee director vest over a three-year period, with one third vesting on each of the first, second, and third anniversaries of the applicable grant date, provided that the non-employee director is a director of the Company on the applicable vesting date.

We reimburse in full each non-employee director for all reasonable out-of-pocket expenses incurred in attending each meeting of our board of directors or its committees. Directors may be reimbursed for travel, food, lodging and other expenses directly related to their service as directors. Directors are also entitled to the protection provided by their indemnification agreements, and the indemnification provisions in our restated certificate of incorporation and restated bylaws.

EQUITY COMPENSATION PLAN INFORMATION

The following table provides certain aggregate information with respect to all of our equity compensation plans in effect as of December 31, 2018.

<u>Plan category</u>	<u>(a)</u> Number of securities to be issued upon exercise of outstanding options, warrants and rights	<u>(b)</u> Weighted-average exercise price of outstanding options, warrants and rights	<u>(c)</u> Number of securities remaining available for future issuance under equity compensation plans (excluding securities reflected in column (a))
Equity compensation plans approved by security holders(1)	2,476,911	\$ 9.65	4,393
Equity compensation plans not approved by security holders	—	—	—
Total	2,476,911	\$ 9.65	4,393

- (1) These plans consist of our 2017 Employee, Director and Consultant Equity Incentive Plan (the "2017 Plan"), our 2007 Stock Option and Grant Plan, as amended, and our 2017 Employee Stock Purchase Plan (the "ESPP"). Does not include an additional 894,761 shares of common stock reserved for future issuance under the 2017 Plan effective January 1, 2019 by operation of the 2017 Plan's "evergreen" provision and an additional 223,690 shares of common stock reserved for future issuance under the ESPP effective January 1, 2019 by operation of the ESPP's "evergreen" provision.

REPORT OF AUDIT COMMITTEE

The audit committee of the board of directors, which consists entirely of directors who meet the independence and experience requirements of The Nasdaq Stock Market, has furnished the following report:

The audit committee assists the board of directors in overseeing and monitoring the integrity of our financial reporting process, compliance with legal and regulatory requirements and the quality of internal and external audit processes. This committee's role and responsibilities are set forth in our charter adopted by the board of directors, which is available on our website at www.quanterix.com. This committee reviews and reassesses our charter annually and recommends any changes to the board of directors for approval. The audit committee is responsible for overseeing our overall financial reporting process, and for the appointment, compensation, retention, and oversight of the work of Ernst & Young LLP. In fulfilling its responsibilities for the financial statements for the fiscal year ended December 31, 2018, the audit committee took the following actions:

- Reviewed and discussed the audited financial statements for the fiscal year ended December 31, 2018 with management and Ernst & Young LLP, our independent registered public accounting firm;
- Discussed with Ernst & Young LLP the matters required to be discussed in accordance with Auditing Standard No. 1301—*Communications with Audit Committees*; and
- Received written disclosures and the letter from Ernst & Young LLP regarding its independence as required by applicable requirements of the Public Company Accounting Oversight Board regarding Ernst & Young LLP's communications with the audit committee and the audit committee further discussed with Ernst & Young LLP their independence. The audit committee also considered the status of pending litigation, taxation matters and other areas of oversight relating to the financial reporting and audit process that the committee determined appropriate.

Based on the audit committee's review of the audited financial statements and discussions with management and Ernst & Young LLP, the audit committee recommended to the board of directors that the audited financial statements be included in our Annual Report on Form 10-K for the fiscal year ended December 31, 2018 for filing with the SEC.

Members of the Audit Committee:

Paul M. Meister (*Chairman*)
John M. Connolly
Martin D. Madaus, Ph.D.

SECTION 16(a) BENEFICIAL OWNERSHIP REPORTING COMPLIANCE

Section 16(a) of the Securities Exchange Act of 1934, as amended, or the Exchange Act, requires the Company's officers and directors, and persons who beneficially own 10% or greater of a registered class of the Company's equity securities, to file reports of ownership on Form 3 and reports of changes in ownership on Form 4 or Form 5 with the SEC. Such officers, directors and 10% or greater stockholders are also required by SEC rules to furnish the Company with copies of all Section 16(a) forms they file. To our knowledge, based solely on a review of our records and representations made by our directors and officers regarding their filing obligations, all reports which were required to be filed pursuant to Section 16(a) of the Exchange Act in 2018 were filed on a timely basis.

CERTAIN RELATIONSHIPS AND RELATED PERSON TRANSACTIONS

Since January 1, 2017, we have engaged in the following transactions with our directors, director nominees, executive officers and holders of more than 5% of our common stock, which we refer to as our principal stockholders, and affiliates or immediate family members of our directors, executive officers and principal stockholders. We believe that all of these transactions were on terms as favorable as could have been obtained from unrelated third parties.

Some of our directors are or were recently affiliated with our principal stockholders as indicated below:

<u>Director</u>	<u>Affiliation with principal stockholder</u>
Keith L. Crandell	Managing Director of ARCH Venture Partners VI, LLC, the ultimate general partner of ARCH Venture Fund VI, L.P., and Managing Director of ARCH Venture Partners VIII, LLC, the general partner of ARCH Venture Fund VIII Overage, L.P.
Douglas G. Cole, M.D.	Managing Partner of Flagship Pioneering and a member of Flagship Ventures General Partner LLC, the sole general partner of Flagship Ventures Fund 2004, L.P.
John M. Connolly	Senior Advisor to Bain Capital Ventures, the ultimate general partner of Bain Capital Venture Fund 2005, L.P., BCIP Associates III, LLC and BCIP Associates III-B, LLC.

Participation in Our Initial Public Offering

In December 2017, we issued an aggregate of 4,916,480 shares of our common stock at a purchase price of \$15.00 per share, which included 641,280 shares of common stock issued upon exercise by the underwriters of their option to purchase additional shares. Gross proceeds from the offering were approximately \$73 million, prior to deducting underwriting discounts and commissions and offering expenses payable by the Company.

The table below sets forth the aggregate number of shares of our common stock issued to our directors, executive officers or holders of more than 5% of our capital stock, or an affiliate or immediate family member thereof, at the time of the transaction:

	<u>Shares</u>	<u>Aggregate Purchase Price</u>
David R. Walt, Ph.D.	200,000	\$ 3,000,000
Marijn Dekkers, Ph.D.	40,000	\$ 600,000
E. Kevin Hrusovsky	34,000	\$ 510,000
Arch Venture Fund VIII Overage	266,000	\$ 3,990,000
Cormorant Global Healthcare Fund, LP	100,000	\$ 1,500,000
Trinitas Innovation-Q Investment Co., Ltd.	55,000	\$ 825,000

Series D-1 Preferred Stock Financing

In June 2017, we issued an aggregate of 2,113,902 shares of our Series D-1 preferred stock at \$4.021 per share for an aggregate purchase price of approximately \$8.5 million, which we refer to as the Series D-1 preferred stock financing. Each share of our Series D-1 preferred stock converted automatically into 0.3111 shares of our common stock immediately prior to the completion of our initial public offering on December 11, 2017.

The following table summarizes the Series D-1 preferred stock purchased by our directors, executive officers and principal stockholders.

<u>Name of Purchaser</u>	<u>Number of Shares of Series D-1 Preferred Stock Purchased</u>	<u>Aggregate Purchase Price</u>
<u>Marijn Dekkers, Ph.D.(1)</u>	248,694	\$ 999,999

(1) Dr. Dekkers is a director of the Company.

License Agreement with Tufts University

We entered into a License Agreement with Tufts University pursuant to which we obtained an exclusive, worldwide license to research, develop, commercialize, use, make, or have made, import or have imported, distribute or have distributed, offer or have offered, and sell or have sold products and services covered by patent rights to the Simoa technology owned by Tufts University, as well as a non-exclusive license to related know-how. The rights licensed to us are for all fields of use and are sublicensable for a fee. One of our directors and our founding scientist, David R. Walt, Ph.D. previously served as University Professor, Professor of Biomedical Engineering, Professor of Genetics, Professor of Neuroscience, Professor of Cell and Molecular Biology, and Professor of Oral Medicine at Tufts University. Under that agreement, we are required to pay Tufts University royalties on all net sales of products and services as well as a portion of any sublicensing revenues. Tufts University pays a portion of the royalties received from us to Dr. Walt, the amount of which is controlled solely by Tufts University.

Registration Rights

Pursuant to our registration rights agreement to which certain of our principal stockholders, directors and their affiliates are parties, the holders of certain shares of common stock and warrants to purchase shares of our common stock are entitled to registration rights with respect to the shares of common stock held by them. These shares include the shares held (and shares issuable upon the exercise of warrants held) by ARCH Venture Partners and affiliated funds, Bain Capital Venture entities, Flagship Ventures Fund 2005, L.P., and our directors Marijn Dekkers, Ph.D. and David R. Walt, Ph.D., and an affiliate of our director Paul M. Meister. See "Security Ownership of Certain Beneficial Owners and Management."

Under our registration rights agreement, holders of registrable shares can demand that we file a registration statement or request that their shares be included on a registration statement that we are otherwise filing, in either case, registering the resale of their shares of common stock. The registration rights will terminate at the earlier of the closing of a Change of Control Event, as such term is defined in our restated certificate of incorporation, (b) such time as Rule 144 under the Securities Act of 1933, as amended, or the Securities Act, or another similar exemption under the Securities Act is available for the sale of all of such registrable shares without limitation during a three-month period without registration (and without the requirement for us to be in compliance with the current public information required under Rule 144(c)(1)), or (3) the fifth anniversary of the completion of our initial public offering.

Demand Registration Rights. The holders of a majority of the registrable shares may require us to file a registration statement under the Securities Act at our expense, subject to certain exceptions, with respect to the resale of their registrable shares having an aggregate offering price (net of underwriting discounts and commissions, if any) of at least \$5 million, and we are required to use our reasonable best efforts to effect the registration. The holders of a majority of the registrable shares may require us to effect up to two such demand registrations for the holders of registrable shares as a group. In the

event we are required to effect such a demand registration, we may not effect any other registration of securities for sale for our own account (other than a registration effected solely to implement an employee benefit plan or in certain business combination transactions) within 120 days following the effective date of the demand registration.

We are also required to use our reasonable best efforts to qualify and remain qualified to register securities pursuant to a registration statement on Form S-3 under the Securities Act. Any holder or holders of registrable shares anticipated to have an aggregate sale price (net of underwriting discounts and commissions, if any) in excess of \$1 million will have the right to require us to file, at our expense, an unlimited number of registration statements on Form S-3 for the registrable shares held by such requesting holder or holders, and we are required to use our reasonable best efforts to effect such registrations.

Piggyback Registration Rights. If we propose to register any of our securities under the Securities Act for sale to the public (except with respect to registration statements on Form S-4, Form S-8 or another form not available for registering the registrable shares for sale to the public), the holders of registrable shares are entitled to notice of such registration and to request that we include registrable shares for resale on such registration statement, subject to the right of any underwriter to limit the number of shares included in such registration. The requisite holders of these "piggyback" registration rights have waived their rights in connection with the filing of our registration statement on Form S-3 (File No. 333-230399).

We will pay all registration expenses, other than underwriting fees, commissions or discounts, related to any demand or piggyback registration, including up to \$50,000 of fees and disbursements of counsel for the holders of registrable shares. The Registration Rights Agreement contains customary cross-indemnification provisions, pursuant to which we are obligated to indemnify the selling stockholders in the event of misstatements or omissions in the registration statement attributable to us and they are obligated to indemnify us for misstatements or omissions attributable to them.

Agreements with Directors

On March 20, 2017, we entered into a letter agreement with our director Marijn Dekkers, Ph.D. pursuant to which Dr. Dekkers agreed to serve as a member of our board of directors. Pursuant to the letter agreement, we granted Dr. Dekkers an option to purchase 100,000 shares of our common stock vesting over four years. In addition, we agreed to provide Dr. Dekkers with the opportunity to participate in any private placement financing transaction we entered into prior to completion of our initial public offering on the same terms as other investors in such transaction. We also agreed to pay Dr. Dekkers compensation of \$25,000 per year for service on our board of directors and to reimburse Dr. Dekkers for his reasonable out-of-pocket expenses in attending board and committee meetings. We also agreed to provide indemnification of Dr. Dekkers pursuant to our certificate of incorporation, an indemnification agreement and our director and officer insurance. This agreement terminated upon adoption of our non-employee director compensation policy in 2018.

On August 7, 2013, we entered into a letter agreement with our director Paul M. Meister pursuant to which Mr. Meister agreed to serve as a member of our board of directors. Pursuant to the letter agreement, we granted Mr. Meister an option to purchase 75,000 shares of our common stock vesting over four years. We also agreed to pay Mr. Meister compensation of \$25,000 per year for service on our board of directors and to reimburse Mr. Meister for his reasonable out-of-pocket expenses in attending board and committee meetings. We also agreed to provide indemnification of Mr. Meister pursuant to our certificate of incorporation, an indemnification agreement and our director and officer insurance. This agreement terminated upon adoption of our non-employee director compensation policy in 2018.

On January 1, 2014, we entered into a letter agreement with our director David Walt, Ph.D. pursuant to which we agreed to pay Dr. Walt compensation of \$25,000 per year for service on our board of directors, commencing as of January 1, 2014 and for so long as he continues to serve as a non-employee member of our board of directors, and to reimburse Dr. Walt for his reasonable out-of-pocket expenses in attending board and committee meetings. This agreement terminated upon adoption of our non-employee director compensation policy in 2018.

Indemnification Agreements with Officers and Directors and Directors' and Officers' Liability Insurance

We have entered into indemnification agreements with each of our executive officers and directors. The indemnification agreements, our restated certificate of incorporation and our restated bylaws require us to indemnify our directors to the fullest extent not prohibited by Delaware law. Subject to certain limitations, our restated bylaws also require us to advance expenses incurred by our directors and officers.

Our 2017 Employee, Director and Consultant Equity Incentive Plan and our 2007 Stock Option and Grant Plan, as amended, also provide that our directors (and in the case of the 2017 plan, our employees) will not be liable for any act, omission, interpretation, construction or determination made in good faith in connection with the respective plan, and the members of our board of directors and any committee administering the plan will be entitled to indemnification and reimbursement by us in respect of any claim, loss, damage or expense (including reasonable attorneys' fees) arising or resulting therefrom to the fullest extent permitted by law and/or under any directors' and officers' liability insurance coverage which may be in effect from time to time.

We also maintain a general liability insurance policy which covers certain liabilities of directors and officers of our company arising out of claims based on acts or omissions in their capacities as directors or officers.

Policies and Procedures for Related Party Transactions

Pursuant to its charter, our audit committee is responsible for reviewing and approving, prior to our entry into any such transaction, all transactions reportable by us under Item 404 of Regulation S-K in which we are a participant and in which any parties related to us, including our executive officers, our directors, beneficial owners of more than 5% of our securities, immediate family members of the foregoing persons and any other persons whom our board of directors determines may be considered related persons under Item 404 of Regulation S-K, has or is expected to have a direct or indirect material interest.

In reviewing and approving such transactions, the audit committee will obtain, or will direct our management to obtain on its behalf, all information that the committee believes to be relevant and important to a review of the transaction prior to its approval. Following receipt of the necessary information, a discussion will be held of the relevant factors if deemed to be necessary by the committee prior to approval. If a discussion is not deemed to be necessary, approval may be given by written consent of the committee. This approval authority may also be delegated to the chair of the audit committee in some circumstances. No related person transaction will be entered into prior to the completion of these procedures.

The audit committee or its chair, as the case may be, will approve only those related person transactions that are determined to be in, or not inconsistent with, the best interests of us and our stockholders, taking into account all available facts and circumstances as the committee or the chair determines in good faith to be necessary in accordance with principles of Delaware law generally applicable to directors of a Delaware corporation. These facts and circumstances will typically include, but not be limited to, the benefits of the transaction to us; the impact on a director's independence in

the event the related person is a director, an immediate family member of a director or an entity in which a director is a partner, stockholder or executive officer; the availability of other sources for comparable products or services; the terms of the transaction; and the terms of comparable transactions that would be available to unrelated third parties or to employees generally. No member of the audit committee will participate in any review, consideration or approval of any related person transaction with respect to which the member or any of his or her immediate family members has an interest.

ELECTION OF DIRECTORS

(Notice Item 1)

Our board of directors nominated Marijn Dekkers, Ph.D., Sarah E. Hlavinka and David R. Walt, Ph.D. for election at the annual meeting. Our board of directors currently consists of eight members, classified into three classes as follows: (1) Douglas G. Cole, M.D., Marijn Dekkers, Ph.D., and David R. Walt, Ph.D. constitute Class II with a term ending at the upcoming annual meeting of stockholders; (2) E. Kevin Hrusovsky, Keith L. Crandell and Paul M. Meister constitute Class III with a term ending at the 2020 annual meeting of stockholders; and (3) John M. Connolly and Martin D. Madaus, Ph.D. constitute Class I with a term which expires at the 2021 annual meeting of stockholders. At each annual meeting of stockholders, directors are elected for a full term of three years to succeed those directors whose terms are expiring.

Our board of directors has voted to nominate Marijn Dekkers, Ph.D., Sarah E. Hlavinka and David R. Walt, Ph.D. for election at the annual meeting as Class II directors for a term of three years to serve until the 2022 annual meeting of stockholders, and until their respective successors are elected and qualified. The Class III directors (E. Kevin Hrusovsky, Keith L. Crandell and Paul M. Meister) and the Class I directors (John M. Connolly and Martin D. Madaus, Ph.D.) will serve until the annual meetings of stockholders to be held in 2020 and 2021, respectively, and until their respective successors have been elected and qualified.

Unless authority to vote for any of these nominees is withheld, the shares represented by the enclosed proxy will be voted **FOR** the election of Marijn Dekkers, Ph.D., Sarah E. Hlavinka and David R. Walt, Ph.D. as directors. In the event that any of the nominees becomes unable or unwilling to serve, the shares represented by the enclosed proxy will be voted for the election of such other person as the board of directors may recommend in that nominee's place. We have no reason to believe that any of the nominees will be unable or unwilling to serve as a director.

A plurality of the shares voted **FOR** each nominee is required to elect each nominee as a director.

THE BOARD OF DIRECTORS RECOMMENDS THE ELECTION OF MARIJN DEKKERS, PH.D., SARAH E. HLAVINKA AND DAVID R. WALT, PH.D. AS DIRECTORS, AND PROXIES SOLICITED BY THE BOARD WILL BE VOTED IN FAVOR THEREOF UNLESS A STOCKHOLDER HAS INDICATED OTHERWISE ON THE PROXY.

**RATIFICATION OF THE APPOINTMENT OF THE INDEPENDENT
REGISTERED PUBLIC ACCOUNTING FIRM**

(Notice Item 2)

The audit committee has appointed Ernst & Young LLP as our independent registered public accounting firm, to audit our financial statements for the fiscal year ending December 31, 2019. The board of directors proposes that the stockholders ratify this appointment. Ernst & Young LLP audited our financial statements for the fiscal years ended December 31, 2018 and 2017. We expect that representatives of Ernst & Young LLP will be present at the annual meeting, will be able to make a statement if they so desire, and will be available to respond to appropriate questions.

In deciding to appoint Ernst & Young LLP, the audit committee reviewed auditor independence issues and existing commercial relationships with Ernst & Young LLP and concluded that Ernst & Young LLP has no commercial relationship with the Company that would impair its independence for the fiscal year ending December 31, 2019.

The following table presents fees for professional audit services rendered by Ernst & Young LLP, our independent registered public accounting firm, for the audit of our annual financial statements for the years ended December 31, 2018 and 2017, and fees billed for other services rendered by Ernst & Young LLP during those periods.

	<u>2018</u>	<u>2017</u>
Audit fees:(1)	\$ 892,738	\$ 1,498,921
Tax fees:(2)	\$ 59,523	\$ 48,295
Total	\$ 952,261	\$ 1,547,216

- (1) Audit fees consisted of audit work performed in the preparation of financial statements, as well as work generally only the independent registered public accounting firm can reasonably be expected to provide, such as statutory audits.
- (2) Tax fees consist principally of assistance with matters related to tax compliance and reporting and tax consulting services primarily related to international tax planning strategies.

Policy on Audit Committee Pre-Approval of Audit and Permissible Non-audit Services of Independent Public Accountant

Consistent with SEC policies regarding auditor independence, the audit committee has responsibility for appointing, setting compensation and overseeing the work of our independent registered public accounting firm. In recognition of this responsibility, the audit committee has established a policy to pre-approve all audit and permissible non-audit services provided by our independent registered public accounting firm; provided however, that de minimis non-audit services may instead be approved in accordance with applicable SEC rules.

Prior to engagement of an independent registered public accounting firm for the next year's audit, management will submit an aggregate of services expected to be rendered during that year for each of four categories of services to the audit committee for approval.

1. **Audit** services include audit work performed in the preparation of financial statements, as well as work that generally only an independent registered public accounting firm can reasonably be expected to provide, including comfort letters, statutory audits, and attest services and consultation regarding financial accounting and/or reporting standards.

2. **Audit-Related** services are for assurance and related services that are traditionally performed by an independent registered public accounting firm, including due diligence related to mergers and acquisitions, employee benefit plan audits, and special procedures required to meet certain regulatory requirements.
3. **Tax** services include all services performed by an independent registered public accounting firm's tax personnel except those services specifically related to the audit of the financial statements, and includes fees in the areas of tax compliance, tax planning, and tax advice.
4. **Other Fees** are those associated with services not captured in the other categories. The Company generally does not request such services from our independent registered public accounting firm.

Prior to engagement, the audit committee pre-approves these services by category of service. The fees are budgeted and the audit committee requires our independent registered public accounting firm and management to report actual fees versus the budget periodically throughout the year by category of service. During the year, circumstances may arise when it may become necessary to engage our independent registered public accounting firm for additional services not contemplated in the original pre-approval. In those instances, the audit committee requires specific pre-approval before engaging our independent registered public accounting firm.

The audit committee may delegate pre-approval authority to one or more of its members. The member to whom such authority is delegated must report, for informational purposes only, any pre-approval decisions to the audit committee at its next scheduled meeting.

In the event the stockholders do not ratify the appointment of Ernst & Young LLP as our independent registered public accounting firm, the audit committee will reconsider its appointment.

The affirmative vote of a majority of the votes cast affirmatively or negatively for this proposal is required to ratify the appointment of the independent registered public accounting firm.

THE BOARD OF DIRECTORS RECOMMENDS A VOTE TO RATIFY THE APPOINTMENT OF ERNST & YOUNG LLP AS OUR INDEPENDENT REGISTERED PUBLIC ACCOUNTING FIRM, AND PROXIES SOLICITED BY THE BOARD WILL BE VOTED IN FAVOR OF SUCH RATIFICATION UNLESS A STOCKHOLDER INDICATES OTHERWISE ON THE PROXY.

CORPORATE CODE OF CONDUCT AND ETHICS

We have adopted a corporate code of conduct and ethics that applies to all of our employees, including our chief executive officer and chief financial and accounting officers. The text of the code of conduct and ethics is posted on our website at www.quanterix.com and will be made available to stockholders without charge, upon request, in writing to the Corporate Secretary at 900 Middlesex Turnpike, Building 1, Billerica, Massachusetts 01821. Disclosure regarding any amendments to, or waivers from, provisions of the code that apply to our directors, principal executive officer or principal financial officer will be included in a Current Report on Form 8-K within four business days following the date of the amendment or waiver, unless website posting or the issuance of a press release of such amendments or waivers is then permitted by the rules of The Nasdaq Stock Market.

OTHER MATTERS

The board of directors knows of no other business which will be presented to the annual meeting. If any other business is properly brought before the annual meeting, proxies will be voted in accordance with the judgment of the persons named therein.

STOCKHOLDER PROPOSALS AND NOMINATIONS FOR DIRECTOR

To be considered for inclusion in the proxy statement relating to our 2020 annual meeting of stockholders, we must receive stockholder proposals (other than for director nominations) no later than February 7, 2020. To be considered for presentation at the 2020 annual meeting, although not included in the proxy statement, proposals (including director nominations that are not requested to be included in our proxy statement) must be received no earlier than March 27, 2020 and no later than April 26, 2020. Proposals that are not received in a timely manner will not be voted on at the 2020 annual meeting. If a proposal is received on time, the proxies that management solicits for the meeting may still exercise discretionary voting authority on the proposal under circumstances consistent with the proxy rules of the SEC. All stockholder proposals should be marked for the attention of Secretary, Quanterix Corporation, 900 Middlesex Turnpike, Building 1, Billerica, Massachusetts 01821.

Billerica, Massachusetts
June 5, 2019

2019 Annual Meeting Admission Ticket

2019 Annual Meeting of Quanterix Corporation Stockholders

July 25, 2019 10:00 AM ET

Quanterix Corporation
900 Middlesex Turnpike, Building 1
Billerica, MA 01821

Upon arrival, please present this admission ticket and photo identification at the registration desk.

Important notice regarding the Internet availability of proxy materials for the Annual Meeting of Stockholders.
The material is available at: www.investorvote.com/QTRX



▼ IF VOTING BY MAIL, SIGN, DETACH AND RETURN THE BOTTOM PORTION IN THE ENCLOSED ENVELOPE. ▼

Quanterix Corporation



Notice of 2019 Annual Meeting of Stockholders

Proxy Solicited by Board of Directors for Annual Meeting – July 25, 2019

E. Kevin Hrusovsky and Amol Chaubal, or any of them, each with the power of substitution, are hereby authorized to represent and vote the shares of the undersigned, with all the powers which the undersigned would possess if personally present, at the Annual Meeting of Stockholders of Quanterix Corporation to be held on July 25, 2019 or at any postponement or adjournment thereof.

Shares represented by this proxy will be voted as directed by the stockholder. If no such directions are indicated, the Proxies will have authority to vote FOR the election of the nominees as directors and FOR Proposal 2.

In their discretion, the Proxies are authorized to vote upon such other business as may properly come before the meeting.

(Items to be voted appear on reverse side)

C Non-Voting Items

Change of Address – Please print new address below.

Comments – Please print your comments below.

