

**UNITED STATES
SECURITIES AND EXCHANGE COMMISSION**
Washington, D.C. 20549

FORM 8-K

CURRENT REPORT
Pursuant to Section 13 or 15(d)
of the Securities Exchange Act of 1934

Date of Report (Date of earliest event reported): August 3, 2023

QUANTERIX CORPORATION

(Exact name of registrant as specified in its charter)

Delaware
(State or other jurisdiction
of incorporation)

001-38319
(Commission File Number)

20-8957988
(IRS Employer
Identification No.)

900 Middlesex Turnpike
Billerica, MA
(Address of principal executive offices)

01821
(Zip Code)

Registrant's telephone number, including area code: (617) 301-9400

Check the appropriate box below if the Form 8-K filing is intended to simultaneously satisfy the filing obligation of the registrant under any of the following provisions:

- Written communications pursuant to Rule 425 under the Securities Act (17 CFR 230.425)
- Soliciting material pursuant to Rule 14a-12 under the Exchange Act (17 CFR 240.14a-12)
- Pre-commencement communications pursuant to Rule 14d-2(b) under the Exchange Act (17 CFR 240.14d-2(b))
- Pre-commencement communications pursuant to Rule 13e-4(c) under the Exchange Act (17 CFR 240.13e-4(c))

Securities registered pursuant to Section 12(b) of the Act:

Title of each class	Trading symbol(s)	Name of each exchange on which registered
Common Stock, \$0.001 par value per share	QTRX	The Nasdaq Global Market

Indicate by check mark whether the registrant is an emerging growth company as defined in Rule 405 of the Securities Act of 1933 (17 CFR §230.405) or Rule 12b-2 of the Securities Exchange Act of 1934 (17 CFR §240.12b-2).

Emerging Growth Company

If an emerging growth company, indicate by check mark if the registrant has elected not to use the extended transition period for complying with any new or revised financial accounting standards provided pursuant to Section 13(a) of the Exchange Act.

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

(b) Effective as of August 21, 2023, Michael Doyle will cease serving as the Chief Financial Officer and Treasurer of Quanterix Corporation (the “Company”), as previously contemplated by the terms of his amended and restated employment agreement dated March 27, 2023 and effective with the appointment of Vandana Sriram (as described below). Mr. Doyle will remain employed by the Company as Executive Director, Finance to support finance department operations and the Company until his retirement on March 31, 2024, as disclosed in a Current Report on Form 8-K filed by the Company on March 27, 2023.

(c) On August 9, 2023, the Company issued a press release announcing the appointment of Vandana Sriram as Chief Financial Officer of the Company, effective August 21, 2023. A copy of the press release is attached hereto as Exhibit 99.1 and is incorporated by reference herein.

Ms. Sriram, age 50, joins the Company from Azenta, Inc. where she served as Senior Vice President - Global Finance and Principal Accounting Officer since November 2022. From September 2021 to October 2022, she served as Azenta’s Vice President of Finance, Corporate Controller and Principal Accounting Officer. Prior to Azenta, she worked at General Electric for over 20 years, most recently as head of financial, planning and analysis of GE Aviation from 2020 to 2021 and previously as Chief Financial Officer for GE Additive from 2018 to 2020, Global Controller for GE Aviation from 2014 to 2018 and in other financial roles. Prior to GE, she spent 5 years at Arthur Andersen in public accounting.

(e) In connection with her employment, Ms. Sriram and the Company have entered into an employment agreement (the “Employment Agreement”) dated August 3, 2023. Under the Employment Agreement, Ms. Sriram’s at-will employment with the Company will commence on August 21, 2023. Her annualized base salary will be \$440,000. Ms. Sriram will be eligible to receive an annual performance bonus with a bonus target of 70% of her annual base salary. Ms. Sriram’s bonus for 2023 will be determined based on her actual base salary earned during the year. Ms. Sriram will also receive a sign-on equity award of \$800,000, consisting of 70% stock options and 30% restricted stock units (the “Sign-On Equity Award”). The Sign-On Equity Award will vest over four years, with one-fourth vesting on the first anniversary of the grant date (the “First Equity Tranche”) and the remainder vesting ratably on a monthly basis over the next three years. Ms. Sriram will also be eligible to receive an annual equity grant as part of the Company’s the Company’s next long-term equity award cycle, which is expected to commence in 2024. The Company will target grant date fair value of the annual equity award of up to \$1,500,000.

If Ms. Sriram’s employment is terminated by the Company without cause or she resigns for good reason, she will receive continued payment of her base salary for six months (the “Severance Period”), payment of an amount equal to her prorated annual target bonus for the year of termination, acceleration of the unvested portion of the First Equity Tranche, if any, and health benefits continuation during the Severance Period. If Ms. Sriram’s employment is terminated by the Company without cause or she resigns for good reason in connection with a change-in-control, all of her outstanding but unvested equity awards will become fully vested. Receipt of the foregoing termination benefits will be subject to Ms. Sriram’s execution of a separation agreement, including certain restrictive covenants and a general release of all claims, in a form acceptable to the Company.

The foregoing description of the principal terms of the Employment Agreement is qualified in its entirety by reference to the Employment Agreement, a copy of which is attached as Exhibit 10.1 and incorporated by reference herein.

Item 9.01 Financial Statements and Exhibits.

(d) Exhibits

Exhibit No.	Description
10.1	Employment Agreement dated August 3, 2023 between Vandana Sriram and the Company.
99.1	Press Release dated August 9, 2023
104	Cover Page Interactive Data File (embedded within the inline XBRL document)

SIGNATURES

Pursuant to the requirements of the Securities Exchange Act of 1934, the Registrant has duly caused this report to be signed on its behalf by the undersigned thereunto duly authorized.

QUANTERIX CORPORATION

By: /s/ Masoud Toloue

Masoud Toloue

Chief Executive Officer

Date: August 9, 2023

The logo for Quanterix, featuring the company name in a bold, sans-serif font against a light green background.

900 Middlesex Turnpike
Building 1
Billerica, MA 01821

tel:617.301.9400
fax:617.301.9401

www.quanterix.com

July 31, 2023

Vandana Sriram
Delivered via Email

Re: **Employment Agreement**

Dear Vandana:

Quanterix Corporation (the "Company") is pleased to offer you the full-time, exempt position of Chief Financial Officer, reporting to Masoud Toloue, Chief Executive Officer. Your start date will be August 21, 2023. This role is required to be on-site 5 days a week in our Billerica, MA office.

Congratulations on this offer and career opportunity with Quanterix!

1. **Base Salary:** The Company will pay you a salary at an annual rate of \$440,000.00, paid at a bi-weekly rate of \$16,923.08 (less all applicable taxes and deductions), subject to periodic review and adjustment at the discretion of the Company. Provided your start date is on or before October 2, 2023, you will be eligible for the 2023 merit increase review process. Merit increases are pro-rated based on start date.
 2. **Annual Bonus:** Your bonus target is 70% of your annual base salary. Provided your start date is on or before October 2, 2023, you will be eligible for a pro-rated bonus in the 2023 annual performance bonus program. The actual bonus award amount is subject to the Company's assessment of your individual performance, as well as existing business conditions of the Company. In order to receive the annual bonus, you must be actively employed on the date the bonus is paid. All bonuses are paid at the discretion of the Company and Company's Board of Directors and are subject to the terms of any applicable bonus plan.
 3. **Benefits:** You will be eligible to participate in the employee benefits and insurance programs generally made available to its full-time employees, including medical insurance, dental insurance, 401K Plan, Flexible Spending Account, term life insurance, and short- and long- term disability insurance. Details of these benefits programs, including mandatory employee contributions, will be made available to you when you start. You will also be eligible to participate under the company's Paid Time Off plan. Details of the policy and specific provisions are set forth in the policy itself.
-

4. **Sign-On Equity Award:** You will be eligible to receive a sign-on equity award having a total value of \$800,000.00, consisting of 70% stock options and 30% restricted stock units (RSUs), as determined on a unit basis. Equity awards are granted on an approximately monthly basis, and we will submit your award for approval as soon as administratively practicable after your start date. Your equity award will vest over four years, with 25% vesting on the first anniversary of the grant date. The remainder will vest in thirty-six (36) equal monthly installments, rounded down to the nearest share, on the 15th day of each month beginning with the first full month following the first anniversary of the Grant Date. The number of RSUs and/or options that you will receive will be determined using the Company's standard valuation methodology as of the date your award is approved, and the exercise price of your stock options will be our closing stock price on the approval date. Your award will also be subject to the terms of our 2017 Employee, Director and Consultant Equity Incentive Plan and the applicable award agreements.
5. **Long-Term Equity Incentive:** You will also be eligible to receive an annual equity grant as part of the Company's next long term equity incentive award cycle, which we expect to commence in the first quarter of 2024. The Company will target a grant date fair value of the annual equity award up to \$1,500,000.00, consisting of 70% stock options and 30% restricted stock units (RSUs), as determined on a unit basis. The value of your award will be discretionary and will be subject to your achievement of the metrics and goals established by the Company. Your equity grant will be subject to valuation methodologies and other terms and conditions applicable to other similarly situated executives of the Company and will be subject to Compensation Committee approval.
6. **At-Will Employment: Accrued Obligations:** Your employment is "at will," meaning you or the Company may terminate your employment at any time for any or no reason. In the event of the termination of your employment for any reason, the Company shall pay you the "Accrued Obligations," defined as (1) your base salary through the date of termination; (2) the amount of any expenses properly incurred by you on behalf of the Company prior to any such termination and not yet reimbursed; and (3) to the extent not theretofore paid or provided, any other amounts or benefits required to be paid or provided or which you have earned under any plan or agreement of or with the Company through the date of termination.
7. **Severance:** Without limiting the at-will nature of your employment relationship, if the Company terminates your employment without Cause, or if you resign for Good Reason, the Company shall provide you with the following termination benefits (the "Termination Benefits").
- (a) **Salary Continuation Payments.** Continuation of your base salary for a period of six (6) months after the date of termination (the "Severance Period") at the salary rate then in effect.
 - (b) **Target Bonus.** An amount equal to your annual target bonus for the year of termination, pro-rated as of the date of termination and paid in one lump sum on the Company's next regularly-scheduled payroll date following the effective date of the separation agreement described below.
-

- (c) Acceleration of Sign-On Equity Award. Notwithstanding anything to the contrary in the applicable equity plan or the award agreement applicable to your Sign-On Equity Award, any outstanding but unvested portion of your Sign-On Equity Award that would have vested between the date of termination and August 21, 2024 had you remained employed during such time shall accelerate and become fully-vested and exercisable as of the later of (A) the termination date, or (B) the effective date of the separation agreement described below.
- (d) Health Benefits Continuation. Continuation of group health plan benefits to the extent authorized by and consistent with 29 U.S.C. § 1161 et seq. (commonly known as "COBRA"), with the cost of the regular premium for such benefits shared in the same relative proportion by the Company and you as in effect on the date of termination until the earlier of (i) the end of the Severance Period; or (ii) the date you become eligible for health benefits through another employer or otherwise become ineligible for COBRA ("Health Benefits Continuation Payments"). Notwithstanding the above, (x) in the event that the Severance Period extends beyond eighteen (18) months following your date of termination, or (y) if the Company otherwise determines in its sole discretion that it cannot provide the foregoing Health Benefits Continuation Payments without potentially violating applicable law (including but not limited to the 2010 Patient Protection and Affordable Care Act, as amended by the 2010 Health Care and Education Reconciliation Act), the Company shall in lieu thereof provide to you a taxable monthly payment in an amount equal to the Company's portion of the monthly COBRA premium (as described above) that you would be required to pay to continue your group health coverage in effect on the date of your termination (which amount shall be based on the premium for the first month of COBRA coverage), which payments shall be made on the last day of each month regardless of whether COBRA continuation coverage remains available (i.e., in the event that the Severance Period extends beyond eighteen (18) months following your date of termination) and shall end on the earlier of (1) the end of the Severance Period, (2) the date you become eligible for health benefits through another employer or otherwise become ineligible for COBRA; or (3) the last day of the twenty-fourth (24th) calendar month following your termination date.

If the Company terminates your employment without Cause, or if you resign for Good Reason, and the effective date of such termination occurs within the 90 day period immediately preceding or the twelve (12) month period immediately following a Change- in-Control (such period the "Change-in-Control Period" and such termination a "Change-in- Control Termination"), then in addition to the Termination Benefits set forth immediately above:

Equity Acceleration. Notwithstanding anything to the contrary in any applicable equity plan or award agreement, all of your outstanding but unvested equity awards shall accelerate and become fully-vested and exercisable as of the later of (A) the termination date, or (B) the effective date of the separation agreement described below, or (C) as of the Change-in-Control.

Notwithstanding anything to the contrary in this Agreement, you shall not be entitled to any Termination Benefits unless (a) within 60 days of your date of termination, you first (i) enter into, do not revoke, and comply with the terms of a separation agreement in a form acceptable to the Company, which shall include a general release in favor of the Company and related persons and entities, and other provisions regarding non-competition, confidentiality, cooperation, non-disparagement and the like as may be included in the Company's then current form of separation agreement (the "Release"); (ii) resign from any and all positions, including, without implication of limitation, as a director, trustee, and officer, that you then hold with the Company and any affiliate of the Company; and (iii) return all Company property and comply with any instructions related to deleting and purging duplicates of such Company property, and (b) you comply with the terms of your Non-Competition, Non-Solicitation, Confidentiality and Assignment Agreement or any other similar agreements with the Company. The Salary Continuation Payments shall commence within 60 days after the date of termination and shall be made on the Company's regular payroll dates; provided, however, that if the 60-day period begins in one calendar year and ends in a second calendar year, the Salary Continuation Payments shall begin to be paid in the second calendar year. In the event you miss a regular payroll period between the date of termination and the first Salary Continuation Payment, the first Salary Continuation Payment shall include a "catch up" payment.

For purposes of this Section:

"Cause" means the occurrence of any of the following (and, if applicable, that the Company has complied with the Cause Process (hereinafter defined) following the occurrence of a circumstance subject to the Cause Process): (i) theft, fraud, embezzlement, misappropriation of assets or property of the Company, or material violation of your Non-Competition, Non-Solicitation, Confidentiality and Assignment Agreement; (ii) dishonesty, gross negligence, misconduct, neglect of duties, or breach of fiduciary duty to the Company; (iii) violation of federal or state securities laws; (iv) breach of an employment, consulting or other agreement with the Company; (v) the conviction of a felony, or any crime involving moral turpitude, including a plea of guilty or *nolo contendere*; or (vi) continued, willful and deliberate non-performance by you of your duties hereunder (other than by reason of your physical or mental illness, incapacity or disability).

"Cause Process" means that (1) the Company has reasonably determined in good faith that a "Cause" condition has occurred; (2) the Company has notified you in writing of the first occurrence of the Cause condition within 60 days of the first occurrence of such condition; (3) you are provided a period of 30 days following such notice (the "Cause Cure Period") to remedy the condition; (4) notwithstanding such efforts, the Company reasonably and in good faith determines at the end of the Cause Cure Period that the Cause condition continues to exist; and (5) the Company terminates your employment within 30 days after the end of the Cause Cure Period. If you cure the Cause condition during the Cause Cure Period, Cause shall be deemed not to have occurred. The Company shall not be required to follow the Cause Process as to those conditions which it reasonably determines in good faith cannot be cured within the Cause Cure Period. For the avoidance of doubt, you and the Company acknowledge and agree that clauses (i), (iii) and (v) cannot be cured and shall not be subject to the requirements of the Cause Process.

"Change-in-Control" means the occurrence of any of the following events: (i) any "Person" (as such term is used in Sections 13(d) and 14(d) of the Securities Exchange Act of 1934, as amended) becomes the "Beneficial Owner" (as defined in Rule 13d-3 under said Act), directly or indirectly, of securities of the Company representing fifty percent (50%) or more of the total voting power represented by the Company's then outstanding voting securities; (ii) a change in the composition of the Company's Board of Directors occurring within a two-year period, as a result of which fewer than a majority of the directors are Incumbent Directors. "Incumbent Directors" will mean directors who either (A) are directors of the Company as of the date hereof, or (B) are elected, or nominated for election, to the Board with the affirmative votes of at least a majority of the Incumbent Directors at the time of such election or nomination (but will not include an individual whose election or nomination is in connection with an actual or threatened proxy contest relating to the election of directors to the Company); (iii) the consummation of a merger or consolidation of the Company, other than a merger or consolidation which would result in the voting securities of the Company outstanding immediately prior thereto continuing to represent (either by remaining outstanding or by being converted into voting securities of the surviving entity or the parent of such corporation) at least fifty percent (50%) of the total voting power represented by the voting securities of the Company or such surviving entity or parent of such corporation, as the case may be, outstanding immediately after such merger or consolidation; or (iv) the consummation of the sale or disposition by the Company of all or substantially all of the Company's assets.

"Good Reason" means that you have complied with the Good Reason Process following the occurrence of any of the following actions undertaken by the Company without your express prior written consent: (i) the material diminution in your authority, duties and responsibilities; (ii) a material reduction in your base salary, provided, however, that Good Reason shall not be deemed to have occurred in the event of a reduction in your base salary that is pursuant to a salary reduction program affecting or a material portion of the similarly situated senior level employees of the Company and that does not adversely affect you to a greater extent than such similarly situated employees; and (iii) a change in the geographic location at which you must regularly report to work and perform services of more than thirty (30) miles, except for required travel on the Company's business; or (v) a material breach by the Company of any of its obligations to you under its employment agreements with you.

"Good Reason Process" means that (1) you have reasonably determined in good faith that a "Good Reason" condition has occurred; (2) you have notified the Company in writing of the first occurrence of the Good Reason condition within 30 days of the first occurrence of such condition; (3) the Company is provided with a period of 30 days following such notice (the "Cure Period") to remedy the condition; (4) notwithstanding such efforts, you reasonably and in good faith determine at the end of the Cure Period that the Good Reason condition continues to exist; and (5) you terminate your employment within 30 days after the end of the Cure Period. If the Company cures the Good Reason condition during the Cure Period, Good Reason shall be deemed not to have occurred.

8. Section 280G:

- (a) If any payment or benefit you would receive under this Agreement, when combined with any other payment or benefit you receive pursuant to a Change-in- Control (for purposes of this Section, a "Payment") would constitute a "parachute payment" within the meaning of Section 280G of the Internal Revenue Code of 1986, as amended (the "Code") and, but for this sentence, be subject to the excise tax imposed by Code Section 4999 (the "Excise Tax"), then such Payment shall be either: (i) the full amount of such Payment; or (ii) such lesser amount (a "Reduced Payment") as would result in no portion of the Payment being subject to the Excise Tax, whichever of the foregoing amounts, taking into account the applicable federal, state and local employment taxes, income taxes and the Excise Tax, results in your receipt, on an after-tax basis, of the greater amount of the Payment notwithstanding that all or some portion of the Payment may be subject to the Excise Tax.
 - (b) With respect to Section 9(a), if there is more than one method of reducing the Reduced Payment amount that would result in no portion of the Payment being subject to the Excise Tax, then the Payment shall be reduced or eliminated in the following order: (i) cash payments; (ii) taxable benefits; (iii) nontaxable benefits; and (iv) accelerated vesting of equity awards in a manner that maximizes the amount to be received by you.
 - (c) The determination of whether Section 9(a)(i) or (ii) applies, and the calculation of the amount of the Reduced Payment if applicable, shall be performed by a nationally recognized certified public accounting firm as may be designated by the Company (the "Accounting Firm"). The Accounting Firm shall provide detailed supporting calculations to both the Company and you within fifteen (15) business days of the receipt of notice from you that there has been a Payment, or such earlier time as is requested by the Company, in a form that can be relied upon for tax filing purposes. All fees and expenses of the Accounting Firm shall be borne solely by the Company.
 - (d) You may receive a Payment that is, in the aggregate, either more or less than the amount described in Section 9(a)(i) or (ii) (as applicable, an "Overpayment" or "Underpayment"). If it is finally determined by a court of competent jurisdiction pursuant to a final non-appealable judgment, or the Internal Revenue Service, or by the Accounting Firm upon request by either the Company or you, that an Overpayment or Underpayment has been made, then: (i) in the event of an Overpayment, you shall promptly repay the Overpayment to the Company, together with interest on the Overpayment at the applicable federal rate from the date of your receipt of such Overpayment until the date of such repayment; and (ii) in the event of an Underpayment, the Company shall promptly pay an amount equal to the Underpayment to you, together with interest on such amount at the applicable federal rate from the date such amount would have been paid to you had the provisions of Section 9(a)(ii) not been applied until the date of payment.
-

9. Section 409A: Anything in this Agreement to the contrary notwithstanding, if at the time of your separation from service within the meaning of Section 409A of the Code, the Company determines that you are a "specified employee" within the meaning of Section 409A(a)(2)(B)(i) of the Code, then to the extent any payment or benefit that you become entitled to under this Agreement on account of your separation from service would be considered deferred compensation subject to the 20 percent additional tax imposed pursuant to Section 409A(a) of the Code, such payment shall not be payable and such benefit shall not be provided until the date that is the earlier of (A) six months and one day after your separation from service, or (B) your death. If any such delayed cash payment is otherwise payable on an installment basis, the first payment shall include a catch-up payment covering amounts that would otherwise have been paid during the six-month period but for the application of this provision, and the balance of the installments shall be payable in accordance with their original schedule. All in-kind benefits provided and expenses eligible for reimbursement under this Agreement shall be provided by the Company or incurred by you during the time periods set forth in this Agreement. All reimbursements shall be paid as soon as administratively practicable, but in no event shall any reimbursement be paid after the last day of the taxable year following the taxable year in which the expense was incurred. The amount of in-kind benefits provided or reimbursable expense incurred in one taxable year shall not affect the in-kind benefits to be provided or the expenses eligible for reimbursement in any other taxable year. Such right to reimbursement or in-kind benefits is not subject to liquidation or exchange for another benefit. To the extent that any payment or benefit described in this Agreement constitutes "non-qualified deferred compensation" under Section 409A of the Code, and to the extent that such payment or benefit is payable upon your termination of employment, then such payments or benefits shall be payable only upon your "separation from service." The Company and you intend that this Agreement will be administered in accordance with Section 409A of the Code. To the extent that any provision of this Agreement is ambiguous as to its compliance with Section 409A of the Code, the provision shall be read in such a manner so that all payments hereunder comply with Section 409A of the Code. It is intended that each installment of the severance payments and benefits provided under this Agreement shall be treated as a separate "payment" for purposes of Section 409A. Neither the Company nor you shall have the right to accelerate or defer the delivery of any such payments or benefits except to the extent specifically permitted or required by Section 409A. The Company makes no representation or warranty and shall have no liability to you or any other person if any provisions of this Agreement are determined to constitute deferred compensation subject to Section 409A of the Code but do not satisfy an exemption from, or the conditions of, such Section.

10. *No Guarantee of Tax Consequences:* The Company makes no guarantee of any tax consequences with respect to any payment hereunder, including, without limitation, under Section 409A of the Code.
11. *No Mitigation:* In no event, except as set forth expressly in this or another agreement signed by you, shall you be obligated to seek other employment or take any other action by way of mitigation of the amounts payable to you under any of the provisions of this Agreement and, subject to the aforesaid exception, such amounts shall not be reduced whether or not you obtain other employment.
12. *Return of Company Property:* Upon termination of employment for any reason, you shall promptly return to the Company any keys, credit cards, passes, confidential documents or material, computer equipment, or other property belonging to the Company, and you shall also return all writings, files, records, correspondence, notebooks, notes and other documents and things (including any copies thereof) containing confidential information or relating to the business or proposed business of the Company or its affiliated entities or containing any trade secrets relating to the Company or its affiliated entities. For purposes of the preceding sentence, the term "trade secrets" shall have the meaning ascribed to it under the Uniform Trade Secrets Act. You agree to represent in writing to the Company upon termination of employment that you have complied with the foregoing provisions of this Section.
13. *Assistance with Claims:* You agree that consistent with your business and personal affairs, during and after your employment by the Company, you will assist the Company and its affiliated entities in the defense of any claims, or potential claims that may be made or are threatened to be made against any of them in any action, suit or proceeding, whether civil, criminal, administrative or investigative (a "*Proceeding*"), and will assist the Company and its affiliated entities in the prosecution of any claims that may be made by the Company or its affiliated entities in any Proceeding, to the extent that such claims may relate to your employment or the period of your employment by the Company. The Company agrees to reimburse you for your reasonable out-of-pocket expenses associated with such assistance, including travel expenses. Any amounts to be paid to you pursuant to this Section 14 shall be paid by the Company within no later than thirty (30) days of the date on which you provide documentation to the Company that such expenses were incurred.
-

14. Representation Regarding Other Obligations: This offer is conditioned on your representation that you are not subject to any confidentiality, non-competition agreement or any other similar type of restriction that may affect your ability to devote full time and attention to your work at the Company. If you have entered into any agreement that may restrict your activities on behalf of the Company, please provide me with a copy of the agreement as soon as possible.

15. Restrictive Covenants: You will be required to sign the Company's standard "Employee Non-Competition, Non-Solicitation, Confidentiality and Assignment Agreement" as a condition of your employment. A copy of that Agreement is enclosed. If the Company elects to enforce a non-competition provision for which post-employment payments are required under applicable law ("Non- Competition Payment"), the Company may apply the amount of any Non- Competition Payment to the Termination Benefits.

16. Other Terms: Your employment with the Company shall be on an at-will basis. In other words, you or the Company may terminate employment for any reason and at any time, with or without notice. In addition, as with all employees, our offer to you is contingent on your submission of satisfactory proof of your identity and your legal authorization to work in the United States.

We are excited about the opportunity to work with you at Quanterix. If you have any questions about this information, please do not hesitate to call. Otherwise, please confirm your acceptance of this offer of employment by signing below and returning a copy to me no later than August 3, 2023.

We are confident that with your background and skills, you will have an immediate positive impact on our organization.

[Signature Page Follows]

Sincerely,

/s/ Masoud Toloue

Masoud Toloue
Chief Executive Officer

Offer accepted:

/s/ Vandana Sriram

Name: Vandana Sriram
Date: August 3, 2023

Quanterix Appoints Vandana Sriram as Chief Financial Officer

Billerica, Mass. – August 9, 2023 — Quanterix Corporation (NASDAQ: QTRX), a company fueling scientific discovery and breakthrough diagnostics through ultrasensitive biomarker detection, today announced that it has appointed Vandana Sriram, CPA, as Chief Financial Officer (CFO), effective August 21, 2023. Sriram replaces Michael Doyle, who has served as the Company’s CFO since July 2021. As previously announced, Mike will transition to the role of Executive Director, Finance, focusing on strategic growth programs until his retirement in March 2024.

“Vandana is experienced in developing and executing strategic roadmaps and financial plans to foster growth opportunities,” said Masoud Toloue, CEO at Quanterix. “The executive team and I are thrilled to welcome her as a partner to further strengthen our current market position and build robust scale. I would like to thank Mike for his guidance and partnership over the last year as we kicked off and continue to successfully execute a major corporate transformation. I expect a smooth transition between Mike and Vandana, and I look forward to them working together into 2024.”

Sriram is a global senior executive with more than 20 years of diversified experience supporting the growth of high-tech manufacturing industries through financial strategy and transformation. Most recently, Sriram served as Senior Vice President of Global Finance at Azenta Life Sciences where she was responsible for the Controllershship, FP&A, and segment CFO activities as a member of the leadership team. She was instrumental in the separation of Azenta and Brooks Automation, and the subsequent deployment of capital to three acquisitions within eight months and management of their integration processes. Prior to Azenta, Sriram served as Senior Executive and Head of FP&A for GE Aviation, a leading provider of commercial and military jet engines and components. While there, she led overall financial planning and drove simplification of financial processes and operating rhythms during a highly dynamic COVID downturn period. Throughout her career with GE spanning over 20 years, and at Azenta for two years, she has partnered with operating teams to drive engagement and accountability on financial performance, resulting in more focused operational actions and outcomes.

“I look forward to applying my career experience in financial management and business strategy to a company that is driving innovation in medicine and breaking new ground in the field of scientific research,” said Sriram. “I have great respect for the market momentum Quanterix has created through its Simoa technology for breakthrough discovery in research and clinical applications, and I look forward to building on this foundation for continued growth.”

Sriram brings global experience to Quanterix through her previous roles at GE Additive in Germany, GE Global Growth Organization MENAT (Middle East, North Africa and Turkey) in Dubai, GE Capital India and Arthur Andersen in New Delhi. She is a Chartered Accountant from the Institute of Chartered Accountants of India. Sriram also graduated from the GE Experience Financial Leadership Program and is a registered CPA.

To learn more about Quanterix, visit: www.quanterix.com/about.

About Quanterix

From discovery to diagnostics, Quanterix's ultrasensitive biomarker detection is fueling breakthroughs only made possible through its unparalleled sensitivity and flexibility. The Company's Simoa® technology has delivered the gold standard for earlier biomarker detection in blood, serum or plasma, with the ability to quantify proteins that are far lower than the Limit of Quantification (LoQ) of conventional analog methods. Its industry-leading precision instruments, digital immunoassay technology and CLIA-certified Accelerator laboratory have supported research that advances disease understanding and management in neurology, oncology, immunology, cardiology and infectious disease. Quanterix has been a trusted partner of the scientific community for nearly two decades, powering research published in more than 2,000 peer-reviewed journals. Find additional information about the Billerica, Massachusetts-based company at <https://www.quanterix.com> or follow us on Twitter and LinkedIn.

Forward-Looking Statements

This press release contains forward-looking statements within the meaning of the Private Securities Litigation Reform Act of 1995. Words such as “may,” “will,” “expect,” “plan,” “anticipate,” “estimate,” “intend” and similar expressions (as well as other words or expressions referencing future events, conditions or circumstances) are intended to identify forward-looking statements. Forward-looking statements in this news release are based on Quanterix's expectations and assumptions as of the date of this press release. Each of these forward-looking statements involves risks and uncertainties. Factors that may cause Quanterix's actual results to differ from those expressed or implied in the forward-looking statements in this press release are discussed in Quanterix's filings with the U.S. Securities and Exchange Commission, including the “Risk Factors” sections contained therein. Except as required by law, Quanterix assumes no obligation to update any forward-looking statements contained herein to reflect any change in expectations, even as new information becomes available.

Media Contact:

Maya Nimnicht, PAN Communications
(510) 334-6273
pan.quanterix@pancomm.com

Investor Relations Contact:

Ed Joyce, Quanterix
(610) 306-9917
ejoyce@quanterix.com
