Instruction 1(b)

FORM 4

Check this box if no longer subject to Section 16. Form 4 or Form 5 obligations may continue. See

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL	
OMB Number:	3235-0287
Estimated average burden	
hours per response:	0.5

Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934 or Section 30(h) of the Investment Company Act of 1940 5. Relationship of Reporting Person(s) to Issuer 2. Issuer Name and Ticker or Trading Symbol 1. Name and Address of Reporting Person (Check all applicable) Quanterix Corp [QTRX] DOYLE MICHAEL A Director 10% Owner Officer (give title Other (specify 3. Date of Earliest Transaction (Month/Day/Year) below) below) (First) (Middle) (Last) 03/31/2023 CFO and Treasurer C/O QUANTERIX CORPORATION 4. If Amendment, Date of Original Filed (Month/Day/Year) 6. Individual or Joint/Group Filing (Check Applicable 900 MIDDLESEX TURNPIKE Line) Form filed by One Reporting Person X (Street) Form filed by More than One Reporting 01821 BILLERICA MA Person Rule 10b5-1(c) Transaction Indication (City) (State) (Zip) Check this box to indicate that a transaction was made pursuant to a contract, instruction or written plan that is intended to satisfy the affirmative defense conditions of Rule 10b5-1(c). See Instruction 10. Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned 1. Title of Security (Instr. 3) 2. Transaction 2A. Deemed 4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 and 5. Amount of 6. Ownership 7. Nature Form: Direct (D) or Indirect of Indirect Beneficial Date **Execution Date** Transaction Securities Code (Instr. (Month/Day/Year) if any Beneficially 5) (Month/Day/Year) 8) Owned Following (I) (Instr. 4) Ownership Reported Transaction(s) (Instr. 4) (A) or (D) ν Price Code Amount (Instr. 3 and 4) 03/31/2023 35⁽¹⁾ 90.042(2) Common Stock F D \$11.27 D Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities) 3. Transaction 1. Title of 3A. Deemed 5. Number 6. Date Exercisable and 7. Title and 8. Price of 9. Number of 10. 11. Nature Date (Month/Day/Year) Expiration Date (Month/Day/Year) Conversion Execution Date, Transaction Derivative Security or Exercise if any Code (Instr. Securities Security Securities Form: Beneficial Direct (D) or Indirect Price of Derivative Securities Acquired Underlying Derivative Beneficially Owned Ownership (Instr. 4) (Instr. 3) (Month/Day/Year) 8) (Instr. 5) (A) or Disposed of (D) Security (Instr. 3 and 4) (I) (Instr. 4) Security Following Reported Transaction(s) (Instr. 3, 4 and 5) (Instr. 4) Amount

Date

Exercisable

(A) (D)

Explanation of Responses:

- 1. Shares withheld by Quanterix Corporation solely to cover tax obligations upon the vesting of 115 RSUs.
- 2. Includes 58,360 restricted stock units

Remarks:

/s/ Brian Keane, as Attorney-

Number

Shares

04/03/2023

in-Fact

Title

Expiration

Date

** Signature of Reporting Person

Reminder: Report on a separate line for each class of securities beneficially owned directly or indirectly.

- * If the form is filed by more than one reporting person, see Instruction 4 (b)(v).
- ** Intentional misstatements or omissions of facts constitute Federal Criminal Violations See 18 U.S.C. 1001 and 15 U.S.C. 78ff(a).

Code

Note: File three copies of this Form, one of which must be manually signed. If space is insufficient, see Instruction 6 for procedure.

Persons who respond to the collection of information contained in this form are not required to respond unless the form displays a currently valid OMB Number.

Know all by these presents, that the undersigned hereby constitutes and appoints John Fry and Brian Keane, each of Quanterix Corporation, and Megan Gates, John Condon, Keunjung Cho, Amanda Mei, Anne Leland and Brenda Meyette, each of Mintz, Levin, Cohn, Ferris, Glovsky and Popeo, P.C., signing singly, with full power of substitution, the undersigned's true and lawful attorney-in-fact to:

- (1) execute for and on behalf of the undersigned, forms and authentication documents for EDGAR Filing Access;
- (2) do and perform any and all acts for and on behalf of the undersigned which may be necessary or desirable to complete and execute any such forms and authentication documents;
- (3) execute for and on behalf of the undersigned, in the undersigned's capacity as an officer, director and/or 10% shareholder of the Company, Forms 3, 4 and 5 in accordance with Section 16(a) of the Securities Exchange Act of 1934 and the rules thereunder;
- (4) do and perform any and all acts for and on behalf of the undersigned which may be necessary or desirable to complete and execute any such Form 3, 4 or 5 and timely file such form with the United States Securities and Exchange Commission and any stock exchange or similar authority; and
- (5) take any other action of any type whatsoever in connection with the foregoing which, in the opinion of such attorney-in-fact, may be of benefit to, in the best interest of, or legally required by, the undersigned, it being understood that the documents executed by such attorney-in-fact, on behalf of the undersigned pursuant to this Power of Attorney, shall be in such form and shall contain such terms and conditions as such attorney-in-fact may approve in such attorney-in-fact's discretion.

The undersigned hereby grants to each such attorney-in-fact full power and authority to do and perform any and every act and thing whatsoever requisite, necessary, or proper to be done in the exercise of any of the rights and powers herein granted, as fully to all intents and purposes as the undersigned might or could do if personally present, with full power of substitution or revocation, hereby ratifying and confirming all that such attorney-in-fact, or such attorney-in-fact's substitute or substitutes, shall lawfully do or cause to be done by virtue of this power of attorney and the rights and powers herein granted. The undersigned acknowledges that the foregoing attorneys-in-fact, in serving in such capacity at the request of the undersigned, is not assuming, nor is the Company assuming, any of the undersigned's responsibilities to comply with Section 16 of the Securities Exchange Act of 1934.

This Power of Attorney shall remain in full force and effect until the undersigned is no longer required to file Forms 3, 4 and 5 with respect to the undersigned's holdings of and transactions in securities issued by the Company, unless earlier revoked by the undersigned in a signed writing delivered to the foregoing attorneysin-fact.

IN WITNESS WHEREOF, the undersigned has caused this Power of Attorney to be executed this 18th day of June 2021. /s/ Michael Doyle Michael Doyle