FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, I	D.C.	20549
---------------	------	-------

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP Check this box if no longer subject to Section 16. Form 4 or Form 5 obligations may continue. See Instruction 1(b).

## **OMB APPROVAL** 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

1. Name and Address of Reporting Person*  DOYLE MICHAEL A					2. Issuer Name <b>and</b> Ticker or Trading Symbol Quanterix Corp [ QTRX ]										all app	licable)	g Person(s) to Issi 10% Owr Other (sp		wner
(Last) (First) (Middle) C/O QUANTERIX CORPORATION 900 MIDDLESEX TURNPIKE					3. Date of Earliest Transaction (Month/Day/Year) 07/13/2022									Α	below) below)  CFO and Treasurer				
(Street) BILLER (City)	Street) BILLERICA MA 01821					4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Indiv Line) X					
	Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																		
1. Title of Security (Instr. 3)  2. Transaction Date (Month/Day/Yes				Execution Da		n Date,		Transaction Code (Instr.		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 ar			nd 5) Secu Bene		cially d Following	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)		7. Nature of Indirect Beneficial Ownership (Instr. 4)	
							Co	de	v	Amount	(A) or (D)	Price		Transa	action(s) 3 and 4)	(5	´	(111341.4)	
Common Stock 07/13/20				07/13/202	22				5		434(1)	D	\$18.1	554 18		,651(2)	Г		
Common Stock 08/01/202			22				5	$\neg$	37(3)	D	\$15.61	47(4)	18,614 <sup>(5)</sup>		Г				
	Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities)																		
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	ear)   Execution Date, if any (Month/Day/Year)   Transaction Code (Instr. Sec Acc (A) Disport (Instr. Sec C)   Code (Instr. Sec C)				5. Numl of Derivat Securit Acquire (A) or Dispos of (D) (Instr. 3 and 5)	ive ies ed	Expi	ate Exe ration I nth/Day		Amo Secu Unde Deriv	le and unt of irities erlying vative irity (Instr. 14)	Der Sec (Ins	Price of rivative curity str. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	Ownersh Form: Direct (D or Indirec (I) (Instr.	nership m: ect (D) ndirect	Beneficial Ownership t (Instr. 4)
					Code V (A)		(A) (		Date ) Exercisab		Expiration Date	ı Title	Amount or Number of Shares						

## **Explanation of Responses:**

- $1. \ The sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 1,382 RSUs.$
- 2. Includes 17,703 restricted stock units.
- 3. The sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 115 RSUs.
- 4. This transaction was executed in multiple trades at prices ranging from \$15.58 to \$15.6412. The price reported reflects the weighted average sale price. The reporting person undertakes to provide upon request to the SEC staff, the issuer or any security holder of the issuer full information regarding the number of shares and prices at which the transaction was effected.
- 5. Includes 17,588 restricted stock units.

## Remarks:

/s/ Brian Keane, as Attorney-

in-Fact

08/03/2022

\*\* Signature of Reporting Person

Date

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).

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