FORM 4

## 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									

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

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  (Last) (First) (Middle)  C/O QUANTERIX CORPORATION  900 MIDDLESEX TURNPIKE  (Street)  BILLERICA MA 01821						Issuer Name and Ticker or Trading Symbol Quanterix Corp [ QTRX ]      Date of Earliest Transaction (Month/Day/Year)      17/12/2021  4. If Amendment, Date of Original Filed (Month/Day/Year)							6. I Lin	5. Relationship of Reporting Person(s) to Issuer (Check all applicable)  Director 10% Owner  X Officer (give title Other (specify below)  CFO and Treasurer  6. Individual or Joint/Group Filing (Check Applicable Line)  X Form filed by One Reporting Person				
(City)	(S	tate)	(Zip)											Persoi	filed by More than One Reporting n			lung
Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																		
1. Title of Security (Instr. 3)  2. Transac Date (Month/Da				Execution Date,		3. Transac Code (II					Benefici Owned	es Formally (D) (Following (I) (II)		wnership n: Direct or Indirect nstr. 4)	7. Nature of Indirect Beneficial Ownership			
								Code	v	Amount	(A) o (D)	Price	Reporte Transac (Instr. 3	tion(s)			(Instr. 4)	
Common Stock 07/12/					/2021		A		5,530	5,530 <sup>(1)</sup> A		5,530 <sup>(2)</sup>			D			
	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)	3A. Deemed Execution Date, if any (Month/Day/Year)		ansacti ode (Ins	ion of str. Do Se Ac (A Di of	of		6. Date Exercisable a Expiration Date (Month/Day/Year)			d 7. Title and Am of Securities Underlying Derivative Secu (Instr. 3 and 4)		8. Price of Derivative Security (Instr. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	e S Illy	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership t (Instr. 4)
				Co	ode V	/ (A	A)		Date Exercisabl		expiration Date	Title	Amount or Number of Shares	ber				
Stock Option (right to buy)	\$59.67	07/12/2021			A	12	2,904		(3)	0	7/12/2031	Common Stock	12,904	\$0.00	12,90	4	D	

## **Explanation of Responses:**

- 1. Represents shares of stock issuable on vesting of restricted stock units. The restricted stock units vest as to 25% of the shares on July 12, 2022, and the remaining 75% vest in 36 equal monthly installments at the end of each month beginning on July 31, 2022.
- 3. The option vests as to 25% of the shares on July 12, 2022, and the remaining 75% vests in 36 equal monthly installments at the end of each month beginning on July 31, 2022.

## Remarks:

/s/ Brian Keane, as Attorney-in-07/13/2021 **Fact** 

\*\* Signature of Reporting Person Date

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

- $^{\star}$  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