SEC Form 4																				
FORM 4 UNITE				ED STATES SECURITIES AND EXCHANGE COMMISSION Washington, D.C. 20549														OMB APPROVAL		
Check this box if no longer subject to Section 16. Form 4 or Form 5 obligations may continue. See Instruction 1(b).							to Sectior	n 16(a	a) of the	e Secu	iritie	s Exchan	-	SHIP OMB Number: 3235-0 Estimated average burden hours per response:			0.5			
1. Name and Address of Reporting Person* DOYLE MICHAEL A						or Section 30(h) of the Investment Company Act of 1940 2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX] 3. Date of Earliest Transaction (Month/Day/Year)										telationship eck all applie Directo X Officer below)	cable) or (give title	ig Pers	ier ner pecify	
(Last) (First) (Middle) C/O QUANTERIX CORPORATION 900 MIDDLESEX TURNPIKE						02/02/2023												and Treasurer		
(Street) BILLERICA MA 01821					4.1											Individual or Joint/Group Filing (Check Applicable le) X Form filed by One Reporting Person Form filed by More than One Reporting Person				
(City)	(S	tate)	(Zip)																	
Table I - Non-Deriva 1. Title of Security (Instr. 3) 2. Transa Date (Month/D)					action	ction 2A. Deemed Execution Date			3. Transactio Code (Insti		on tr.	4. Securities Acquired (A) Disposed Of (D) (Instr. 3, 5)		d (A) or	5. Amou Securitie Benefici	nt of 6. Ov es Forn ally (D) o Following d tion(s)		r Indirect str. 4)	'. Nature of Indirect Beneficial Dwnership Instr. 4)	
Common Stock 02/02/					/202	3			A		╡	41,692	(1)	A	\$0.0) 90,118 ⁽²⁾			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. Deeme Execution if any (Month/Day	Date, T	Transa Code (action Instr.	5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		6. Date Exercisa Expiration Date (Month/Day/Yea			r) of Secu		curitie rlying ative \$	Security	8. Price of Derivative Security (Instr. 5)	9. Numbe derivative Securities Beneficia Owned Following Reported Transacti (Instr. 4)	e s Illy J	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)
				c	Code	v	(A)	(D)	Date Exerci	isable	Ex Da	piration ate	Title		Amount or Number of Shares					
Stock Option (right to buy)	\$14.82	02/02/2023			A		97,281		(3	3)	02	2/02/2033	Comn Stoc		97,281	\$0.00	97,281		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 February 2, 2024, and the remaining 75% vest in 36 equal monthly installments at the end of each month beginning on March 15, 2024.

2. Includes 58,590 restricted stock units.

3. The option vests as to 25% of the shares on February 2, 2024, and the remaining 75% vests in 36 equal monthly installments at the end of each month beginning on March 15, 2024.

Remarks:

/s/ Brian Keane, as Attorney-in-Fact 02/06/2023

** Signature of Reporting Person Date

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

 \ast 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 a stock exclant of the form with the formation of the securities and execute any stock exclant of any type whatsoever in a stock exclant of the formation of any type whatsoever in a stock exclant of the formation of any type whatsoever in a stock exclant of any type whatsoever in a stock exclant of the formation of any type whatsoever in a stock exclant of the formation of any type whatsoever in a stock exclant of the formation of any type whatsoever in a stock exclant of the formation of any type whatsoever in a stock exclant of the formation of any type whatsoever in a stock exclant of the formation of any type whatsoever in a stock exclant of the formation of any type whatsoever in a stock exclant of the formation of any type whatsoever in a stock exclant of the formation of any type whatsoever in a stock exclant of the formation of the stock exclant of the stock exclant of the formation of the stock exclant of the stock exclant of the formation of the stock exclant of the stock ex

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