FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Vashington,	D.C. 2	20549
-------------	--------	-------

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

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

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

Name and Address of Reporting Person* DOYLE MICHAEL A						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]									k all app Direc	licable) tor	ng Pe	rson(s) to Is	vner	
(Last)	,	irst) (I	Middle)		3. Date of Earliest Transaction (Month/Day/Year) 04/07/2023										below	er (give title /) CFO and	Trea	Other (s below) asurer	specify	
900 MIDDLESEX TURNPIKE					4. If Amendment, Date of Original Filed (Month/Day/Year)							6. Individual or Joint/Group Filing (Check Applicable Line)								
(Street) BILLER	ICA M	A 0	1821												X Form filed by One Reporting Person Form filed by More than One Reporting Person					
(City)	(S	tate) (2	Zip)		<u> </u>	Check tl	0b5-1(c) Transaction Indication this box to indicate that a transaction was made pursuant to a contract, instruction or written plan that is intended to the affirmative defense conditions of Rule 10b5-1(c). See Instruction 10.										nded to			
		Table	I - No	n-Deriva	tive S	Secui	rities	Acq	uired,	Dis	posed of	, or E	Benef	iciall	y Own	ed				
1. Title of Security (Instr. 3) 2. Transact Date (Month/Day				y/Year) Execu		Deemed cution Date, y hth/Day/Year)					es Acquired (A) Of (D) (Instr. 3,		4 and Secur Benef		cially Following	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)		7. Nature of Indirect Beneficial Ownership (Instr. 4)		
						Code	v	Amount	(A) (D)	or Pi	ice	Transa	action(s) . 3 and 4)			(111341. 4)				
Common Stock 04/07/2						2023			F		1,006(1)	006 ⁽¹⁾ D S		12.63	.63 89,036 ⁽²⁾			D		
Common Stock 04/30/2					2023				F		119(3)	Г	\$	12.64	88	88,917(4)		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)	Derivative Conversion Date Execution Descurity or Exercise (Month/Day/Year) if any			ion Date,	4. Transaction Code (Instr. 8)		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		6. Date Exercisable and Expiration Date (Month/Day/Year)			7. Title and Amount of Securities Underlying Derivative Security (In: 3 and 4)		De Se (In	. Price of lerivative ecurity nstr. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	y	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership (Instr. 4)	
					Code	Code V (A) (D)		(D)	Date Exercisable		Expiration Date	Title	of Share	s						

Explanation of Responses:

- 1. Shares withheld by Quanterix Corporation solely to cover tax obligations upon the vesting of 3,388 RSUs.
- 2. Includes 54,972 restricted stock units.
- 3. Shares withheld by Quanterix Corporation solely to cover tax obligations upon the vesting of 397 RSUs.
- 4. Includes 54,575 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorney-

05/02/2023

in-Fact

** 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