FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

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

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

MENT 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

1. Name and Address of Reporting Person* DOYLE MICHAEL A						Quanterix Corp [QTRX]									ck all app Direc Office	tor er (give title	ng Pe	10% O	wner	
`		st) (M CORPORATION TURNPIKE	Middle)			te of E 8/202		Transa	action (M	/lonth/	Day/Year)			X	belov	v) CFO and	l Trea	below) asurer		
(Street) BILLER (City)	ICA M	A 0	1821 Zip)		4. If Amendment, Date of Orig						f Original Filed (Month/Day/Year)					rividual or Joint/Group Filing (Check Applicable Form filed by One Reporting Person Form filed by More than One Reporting Person				
		Table	I - Non	-Deriva	tive S	Secu	rities	Acq	uired,	Dis	osed of	, or E	Bene	ficiall	y Own	ed				
Date			2. Transac Date (Month/Da	Execution D		Date,	Code (In					s, 4 and Securi Benefi		ies cially Following	Forn (D) c	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)			
									Code	v	Amount	(A) (D)	or F	rice	Transa	ction(s) 3 and 4)			(11150.4)	
Common Stock 02/2					/2023		F		41(1)	I)	\$11	90,077 ⁽²⁾			D				
		Tal									osed of, onvertib				Owne	d				
1. Title of Derivative Security (Instr. 3) 2. Conversion or Exercise (Month/Day/Year) Price of Derivative Security 3. Transaction Date (Month/Day/Year) (Month/Day/Year) (Month/Day/Year)		on Date,	4. Transaction Code (Instr. 8)		of	r osed (. 3, 4	6. Date Expirati (Month/	on Da	Securities Underlying Derivative Security (In 3 and 4) Amo		unt of rities rlying ative rity (Ins	Str.	Price of erivative ecurity estr. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	у	Ownershi Form: Direct (D) or Indirect (I) (Instr. 4	Beneficial Ownership t (Instr. 4)			
				Code	v	(A)	(D)	Date Exercisa	able	Expiration Date	Title	of								

Explanation of Responses:

- $1. \ Shares \ withheld \ by \ Quanterix \ Corporation \ solely \ to \ cover \ tax \ obligations \ upon \ the \ vesting \ of \ 115 \ RSUs.$
- 2. Includes 58,475 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorneyin-Fact

03/01/2023

** 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