Instruction 1(b).

FORM 4

Check this box if no longer subject to Section 16. Form 4 or Form 5 obligations may continue. See

## **UNITED STATES SECURITIES AND EXCHANGE COMMISSION**

Washington, D.C. 20549	9
------------------------	---

STATEMENT	OF CHANGES	IN BENEFICIAL	OWNERSHIP

OMB APPROVAL										
OMB Number:	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*     Toloue Masoud					2. Issuer Name <b>and</b> Ticker or Trading Symbol Quanterix Corp [ QTRX ]										lationship of Reporting ck all applicable) Director		ng Person(s) to Is 10% Ov		
(Last)	(Fi	rst) (M	∕liddle)		3. Date of Earliest Transaction (Month/Day/Year) 05/15/2024									X	Office below	cer (give title ow) President &		Other (specify below) CEO	
900 MIDDLESEX TURNPIKE				4. If A	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	X Form filed by One Reporting Person  Form filed by More than One Reporting  Person					
(City)	(St	ate) (Z	Zip)		Rul	le 10	)b5-	1(c)	Tran	sac	tion Indi	catio	n						
					Check this box to indicate that a transaction was made pursuant to a contract, instruction or written plan that is intended to satisfy the affirmative defense conditions of Rule 10b5-1(c). See Instruction 10.														
		Table	I - No	n-Deriva	tive S	Secui	rities	Acq	uired,	Dis	posed of	, or B	enefi	cially	Own	ed			
1. Title of Security (Instr. 3)  2. Transac Date (Month/Date)			Execution Date,		Date,	Transaction Disposed (Code (Instr. 5)		ies Acquired (A) Of (D) (Instr. 3, 4		4 and Secur Benef Owne		cially Following	6. Own Form: I (D) or I (I) (Inst	Direct of Indirect Etr. 4)	7. Nature of Indirect Beneficial Ownership				
							Code V		Amount	(A) (D)	Prio	e	Reported Transaction(s) (Instr. 3 and 4)				Instr. 4)		
Common	Stock			05/15/2	2024		F		1,120(1)	D	D \$17.88		339,149(2)		I				
		Tal									osed of, convertib				Owned	d			
1. Title of Derivative Security (Instr. 3)  2. Conversion or Exercise Price of Derivative Security  Output  3. Transaction Date Execution Date, if any (Month/Day/Year)			nsaction de (Instr.  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 (Instr. 3 and 4)		8. Price of Derivative Security (Instr. 5)		9. Number derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	Ownershi Form: Direct (D) or Indirect (I) (Instr. 4		Beneficial Ownership (Instr. 4)					
					Code V (A) (D)		Date Expiration Date		Title	Amoun or Numbe of Shares	r								

## **Explanation of Responses:**

- 1. Shares withheld by Quanterix Corporation solely to cover tax obligations upon the vesting of 2,316 RSUs.
- 2. Includes 162,915 restricted stock units.

## Remarks:

/s/ Brian Keane, as Attorneyin-Fact

05/15/2024

\*\* 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.

## POWER OF ATTORNEY

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

in-fact.
IN WITNESS WHEREOF, the undersigned has caused this Power of Attorney to be executed this 8th day of June 2021.
/s/ Masaoud Toloue

undersigned in a signed writing delivered to the foregoing attorneys-

Masoud Toloue