Instruction 1(b).

FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934

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

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP Check this box if no longer subject to Section 16. Form 4 or Form 5 obligations may continue. See

OMB APPROVAL										
OMB Number: 3235-028										
Estimated average burden										
hours per response	: 0.5									

					or Sec	ction 3	0(h) of the I	ivestme	nt Cor	mpany Act o	f 1940							
Name and Address of Reporting Person* The second sec				2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]							5. Relationship of Reporting Person(s) to Issuer (Check all applicable)							
<u>Toloue Masoud</u>											X	Direc	tor	10% Owner				
(Last)	(Fir	rst) (N	Middle)		Date of Earliest Transaction (Month/Day/Year)						\dashv	X	Office	er (give title v)		Other (specify below)		
C/O QUANTERIX CORPORATION					01/31/2023							President & CEO						
900 MIDDLESEX TURNPIKE																		
(Street)					4. If Amendment, Date of Original Filed (Month/Day/Year)							6. Individual or Joint/Group Filing (Check Applicable Line)						
BILLER	CIA M.	Α 0	1821									X	Form filed by One Reporting Person					
,													Form filed by More than One Reporting Person					
(City)	(St	ate) (Z	Zip)															
		Table	I - No	n-Deriva	tive S	ecur	ities Acq	uired,	Dis	posed of	, or Be	nefici	ially	Own	ed			
1. Title of Security (Instr. 3) 2. Transact Date (Month/Day					Execution Date,		3. Transaction Code (Instr. 8) 4. Securities Acquir Disposed Of (D) (Instr. 5)			nd	5. Amount of Securities Beneficially Owned Following Reported		6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)		7. Nature of Indirect Beneficial Ownership (Instr. 4)			
								Code	v	Amount	(A) or (D)	Price		Transaction(s) (Instr. 3 and 4)				(11150.4)
Common Stock 01/31/2				2023			F		493(1)	D	\$14.	14.13 185,330 ⁽²⁾			D			
		Tal					ies Acqu varrants,							Owne	t			
1. Title of Derivative Security (Instr. 3)	erivative Conversion Date Execution Date, ecurity or Exercise (Month/Day/Year) if any		4. Transac Code (li 8)					ite	7. Title and Amount of Securities Underlying Derivative Security (Ins 3 and 4)		Deri Sec	8. Price of Derivative Security Instr. 5) 9. Number derivative Securities Beneficial Owned Following Reported Transactic (Instr. 4)		y [C	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)		

Date Exercisable

Expiration Date

Explanation of Responses:

- 1. Shares withheld by Quanterix Corporation solely to cover tax obligations upon the vesting of 1,408 RSUs.
- 2. Includes 59,100 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorneyin-Fact

Amount Number

of Shares

Title

02/02/2023

** 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).

Code

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.

(A) (D)

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 undersigned in a signed writing delivered to the foregoing attorneys-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

Masoud Toloue