FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

STATEMENT	OF	<b>CHANGES</b>	IN	<b>BENEFICIAL</b>	OWNERSHIP

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

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

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* <u>Toloue Masoud</u>			2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [ QTRX ]								(Che	Relationship of Reporting Person(s) to Issuer (Check all applicable)     X Director 10% Owner						
(Last) (First) (Middle) C/O QUANTERIX CORPORATION 900 MIDDLESEX TURNPIKE				3. Date of Earliest Transaction (Month/Day/Year) 02/02/2023								_ x	Officer (give title below)  President &			Other (s below) CEO	specify	
(Street) BILLER(		IA tate)	01821 (Zip)		4. If An	nendment,	Date of	f Original I	Filed	(Month/Da	ıy/Year	·)	6. Inc Line)		ed by One	Repo	(Check App rting Persor One Repor	1
Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																		
Date			2. Transa Date (Month/Da	Execution Date,		Date,	, Transaction Dispose Code (Instr.			ities Acquired (A) or d Of (D) (Instr. 3, 4 a			5. Amoun Securities Beneficia Owned Fo	s illy ollowing	Form (D) or	: Direct Indirect str. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
								Code	v	Amount		(A) or (D)	Price	Transacti (Instr. 3 a	on(s)			(1115tr. 4)
Common Stock			02/02/	2/2023		A		111,178 <sup>(1)</sup> A		A	\$0.00	296,5	296,508(2)		D			
Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities)																		
Derivative Conversion Date Execution E Security or Exercise (Month/Day/Year) if any			3A. Deemed Execution Da if any (Month/Day/Y	Transaction Code (Instr.		Derivative		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 of derivative Securities Beneficially Owned Following Reported Transaction(s)		10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership (Instr. 4)		
				Co	de V	(A)		Date Exercisab		Expiration Date	Title	OI N	mount r umber f Shares		(Instr. 4)	ion(s)		
Stock Option (right to buy)	\$14.82	02/02/2023		А		259,416		(3)	0	)2/02/2033	Comr		59,416	\$0.00	259,41	16	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 170,278 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-02/06/2023

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.

## 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:

- 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;
- 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;
- do and perform any and all acts for and on behalf of the (4) 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
- take any other action of any type whatsoever in connection with (5) 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-infact'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 8th day of June 2021. /s/ Masaoud Toloue

Masoud Toloue