FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington D.C. 20549

wasnington,	D.C.	20549	

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

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

1. Name and Address of Reporting Person* <u>Toloue Masoud</u>					2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX] 5. Relationship of Reporting Person(s) to Issuer (Check all applicable)															
					~ "	unitor	<u> </u>	<u>01 p</u>	L ×	1101				X	Direc	tor		10% O	wner	
(Last)	(Fir	rst) (N	Middl	e)	Date of Earliest Transaction (Month/Day/Year)									X	Office below	er (give title v)		Other (s	specify	
C/O QUANTERIX CORPORATION					07/05/2022								President & CEO							
		TURNPIKE																		
					4. If Amendment, Date of Original Filed (Month/Day/Year)									6. Individual or Joint/Group Filing (Check Applicable						
(Street)	CIA M		103												ne) X Form filed by One Reporting Person					
BILLER	CIA M.	A 0	182	I 										Form filed by More than One Reporting						
(City)	(St	ate) (Ž	Zip)												Perso	on				
, ,,			.,	Jan Bankar	41		.141	A -		I D:			D 6' -			1				
		labie	1 - 1	Non-Deriva	tive	Secui	rities	AC	quir	ea, Di	sposea c	or, or	Benetic	lally	Own	ea				
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Ye.				Execution Date		, T	3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 at			nd 5) Sec Ber Ow		ount of ities icially d Following	6. Ownership Form: Direct (D) or Indirect (I)		7. Nature of Indirect Beneficial Ownership			
							-	Code V		Amount	(A) or (D)	A) or D) Price		Reported Transaction(s) (Instr. 3 and 4)		(Instr. 4)		(Instr. 4)		
Common	Stock			07/05/202	22			s 4329		432(1)	D	\$16.56	6.5689(2)		81,269(3)		D			
		Tal	ble l	II - Derivati	ve S	ecurit	ties A	Acqu	uire	d, Dis	posed of	or B	eneficia	ılly (Owne	d		<u> </u>		
				(e.g., pu	ıts, c	alls, v	warra	nts	, op	tions,	converti	ble s	curitie	s) ¯						
1. Title of Derivative Security (Instr. 3) 2. Conversion of Exercise Price of Derivative Security (Instr. 3)				Transaction of Code (Instr. Derivativ		rative rities ired r osed)	Expiration Date (Month/Day/Year)			Amo Secu Und Deri Secu	7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		rice of ivative urity tr. 5)	ative derivative rity Securities		0. Ownership Form: Direct (D) or Indirect I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)			
		Code	v	(A)	(D)	Date Exercisab		Expiration Date	Title	Amount or Number of Shares	er									

Explanation of Responses:

- 1. The sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 1,408 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$16.18 to \$17.10. The price reported reflects the weighted average sale price. The reporting person undertakes to provide upon request to the SEC staff, the issuer or any security holder of the issuer full information regarding the number of shares and prices at which the transaction was effected.
- 3. Includes 68,956 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorneyin-Fact

07/06/2022

** 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:

- (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