SEC Form 4	
------------	--

## 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 STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL

OMB Number:	3235-0287
Estimated average b	urden
hours per response:	0.5

Instruction 1(b).	tinue. See	Filed p	oursuant to Section 16(a) of the Securities Exchange Act of 1934 or Section 30(h) of the Investment Company Act of 1940		l	hours per resp	onse: C	0.5
1. Name and Addre Chaubal Ame	ss of Reporting Pers <u>ol</u>	on*	2. Issuer Name <b>and</b> Ticker or Trading Symbol <u>Quanterix Corp</u> [ QTRX ]	(Check a	onship of F II applicab Director Officer (giv	,	on(s) to Issuer 10% Owner Other (specify	
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE		,	3. Date of Earliest Transaction (Month/Day/Year) 08/03/2020	Λ	below) Chief	below)		
(Street) BILLERCIA MA 01821		4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Individ Line)	ual or Joir	Joint/Group Filing (Check Applicable iled by One Reporting Person			
	01821		X	Form filed				
		01021			Form filed by More than O Person		One Reporting	
(City)	(State)	(Zip)						

#### Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned

1. Title of Security (Instr. 3)	2. Transaction Date (Month/Day/Year)	2A. Deemed Execution Date, if any (Month/Day/Year)	3. Transa Code ( 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 and 5)		5. Amount of Securities Beneficially Owned Following Reported	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership	
			Code	v	Amount	Int (A) or Price Transa		Transaction(s) (Instr. 3 and 4)	(1150.4)	(Instr. 4)
Common Stock	08/03/2020		S		111(1)	D	\$32.2295 <sup>(2)</sup>	25,616 <sup>(3)</sup>	D	
Common Stock	08/03/2020		S		3(1)	D	\$33.33	25,613 <sup>(3)</sup>	D	

### Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities)

			( 371	,	,		,	• •				,			
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	3A. Deemed Execution Date, if any (Month/Day/Year)	4. Transa Code ( 8)		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 of derivative Securities Beneficially Owned Following Reported Transaction(s) (Instr. 4)	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)
				Code	v	(A)	(D)	Date Exercisable	Expiration Date	Title	Amount or Number of Shares				

#### Explanation of Responses:

1. The sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 380 RSUs.

2. This transaction was executed in multiple trades at prices ranging from \$32.2444 to \$33.18. 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 21,461 restricted stock units.

#### **Remarks:**

<u>/s/ Brian Keane, as Attorney-</u> <u>in-Fact</u> <u>08/05/2020</u>

\*\* 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 present, that the undersigned hereby constitutes and appoints each of, Megan Gates, John Condon, Brian Shea, Daniel Marden, Jacqueline A. Cannata, Anne T. Leland and Brenda L. Meyette of Mintz, Levin, Cohn, Ferris, Glovsky and Popeo, P.C., and Brian P. Keane and Joseph Driscoll signing singly, with full power of substitution, the undersigned's true and lawful attorney-infact 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 attorneyin-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 2nd day of April, 2019.

/s/ Amol Chaubal Amol Chaubal