SEC Form 3

FORM 3

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

INITIAL STATEMENT OF BENEFICIAL OWNERSHIP OF SECURITIES

OMB APPROVAL

OMB Number: 3235-0104 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

3. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]				
. Relationship of Reporting Persor Check all applicable) Director	10% Owner	(Mon	5. If Amendment, Date of Original Filed (Month/Day/Year)	
X Officer (give title below)	below)	0.110	6. Individual or Joint/Group Filing (Check Applicable Line)	
Chief Financial O	fficer	X		y One Reporting Person y More than One erson
(City) (State) (Zip) Table I - Non-Derivative Securities Beneficially Owned				
2. Amount of Securities 3. Ownership			4. Nature of Indirect Beneficial Ownership (Instr. 5)	
Table II - Derivative Securities Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities)				
		onversion r Exercise	5. Ownership Form: Direct (D)	6. Nature of Indirect Beneficial Ownership (Instr. 5)
Title	Amount D	Derivative Security	or Indirect (I) (Instr. 5)	
	heck all applicable) Director X Officer (give title below) Chief Financial O Chief Financial O e Securities Beneficially mount of Securities eficially Owned (Instr. 4) Securities Beneficially C s, options, convertible 3. Title and Amount of Security	heck all applicable) Director Chief Cive title below) Chief Financial Officer Chief Financial Officer Chief Financial Officer Chief Financial Officer Chief Financial Officer Chief Financial Officer Securities Beneficially Owned s, options, convertible securities Underlying Derivative Security (Instr. 4) Amount or Number of	heck all applicable) Image: Constraint of the securities of the security of the securit	heck all applicable) Image: Constraint of the securities is the security is the secure securits is the security is the security is the secur

Remarks:

No securities are beneficially owned.

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

04/15/2019

Date

** Signature of Reporting Person

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 5 (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