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

Hlavinka Sarah E. (M			Date of Event Requiring Staten Month/Day/Year 17/25/2019	nent	3. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]							
(Last) (First) (Middle) C/O QUANTERIX CORPORATION,					lationship of Reporting Person(s) to Issuer ck all applicable) Director 10% Owner		(M	5. If Amendment, Date of Original Filed (Month/Day/Year)				
900 MIDDLESEX TURNPIKE					Officer (give title Other (specify below) below)			6. Individual or Joint/Group Filing (Check Applicable Line) X Form filed by One Reporting Person				
(Street) BILLERICA	MA	01821								y More than One		
(City)	(State)	(Zip)										
Table I - Non-Derivative Securities Beneficially Owned												
1. Title of Security (Instr. 4)					int of Securities ially Owned (Instr. 4)			4. Nature of Indirect Beneficial Ownership (Instr. 5)				
Table II - Derivative Securities Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities)												
Exp			2. Date Exercisable and Expiration Date (Month/Day/Year)			3. Title and Amount of Securities Underlying Derivative Security (Inst		4. Conversion	Form:	6. Nature of Indirect Beneficial Ownership (Instr. 5)		
		Date Exercisable	Expiration Date	n Title	3	Amount or Number of Shares	Price of Derivative Security	Direct (D) or Indirect (I) (Instr. 5)				

Explanation of Responses:

Remarks:

Exhibit 24.1 - Power of Attorney

No securities are beneficially owned.

/s/ Brian Keane, Attorney-in-

07/25/2019

<u>fact</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 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 presents, that the undersigned hereby constitutes and appoints Brian Keane and John Fry, each of Quanterix Corporation, and Megan Gates, John Condon, Brian Shea and Keunjung Cho, 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
- 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 25th day of July 2019.

/s/ Sarah E. Hlavinka Sarah E. Hlavinka