FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934 or Section 30(h) of the Investment Company Act of 1940

OMB APPROVAL 3235-0287 OMB Number: Estimated average burden hours per response 0.5

1. Name and Addre	1 0		2. Issuer Name and Ticker or Trading Symbol <u>Quanterix Corp</u> [QTRX]		5. Relationship of Reporting Person(s) to Issuer (Check all applicable)					
HRUSOVSK	YEKEVIN		Semicim Oork [41111]	X	Director	10% Owner				
(Last)	(First)	(Middle)	3. Date of Earliest Transaction (Month/Day/Year)	X	Officer (give title below)	Other (specify below)				
C/O QUANTE	()	(<i>)</i>	03/12/2019		Chairman, Presic	,				
113 HARTWEI	L AVENUE									
(Street)			4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Indiv Line) X	/idual or Joint/Group Filir Form filed by One Re					
LEXINGTON	MA	02421			Form filed by More that Person	0				
(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)			4. Securities A Disposed Of (I			Securities Beneficially	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
			Code	v	Amount	(A) or (D)	Price	Transaction(s) (Instr. 3 and 4)		(11311.4)	
Common Stock	03/12/2019		A		31,732 ⁽¹⁾	Α	\$0.00	1,189,227 ⁽²⁾	D		

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

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)	Transaction Code (Instr. 8)		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	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		Transaction(s) (Instr. 4)			
Stock Option (right to buy)	\$22.69	03/12/2019		A		108,177		(3)	03/12/2029	Common Stock	108,177	\$0.00	108,177	D		

Explanation of Responses:

1. Represents shares of common stock issuable upon vesting of restricted stock units. The restricted stock units vest as to 1,983 of the shares on March 31, 2019 and vest as to an additional 661 shares at the end of each month thereafter for 44 months and as to 665 shares at the end of the 45th month thereafter.

2. Includes 39,815 shares of common stock that are subject to performance based vesting and 244,232 restricted stock units.

3. The option vests as to 25% of the shares on March 12, 2020 and vests as to an additional 2.083% of the shares per month for 36 months thereafter.

Remarks:

/s/ Brian Keane, as Attorney-in-03/18/2019 Fact 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 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.

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