FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

	OMB APPROVAL								
	OMB Number:	3235-0287							
	Estimated average burd	den							
1	hours per response:	0.5							

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

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

1. Name and Address of Reporting Person* <u>CRANDELL KEITH</u>						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]									5. Relationship of Reporting Person(s) to Issuer (Check all applicable) X Director X 10% Owner					
(Last) (First) (Middle) 8755 WEST HIGGINS ROAD, SUITE 1025					3. Date of Earliest Transaction (Month/Day/Year) 01/02/2019									Officer below)	(give title		Other (below)	specify		
(Street) CHICAGO IL 60631				_ 4.	If Ame	endment, D	ate of	Original F	Filed ((Month/Day	Individual or Joint/Group Filing (Check Applicable Line) X Form filed by One Reporting Person Form filed by More than One Reporting Person									
(City)	(S		(Zip)																	
1. Title of Security (Instr. 3) 2. Transa						action 2A. Deemed			quired, Disposed of, or Benefic 3. 4. Securities Acquired (A)) or 5. Amount of 6. Ownership 7. Nature					
Date (Month/Da				/Day/Y	ear)	Execution Date, if any (Month/Day/Year)		Transaction Code (Instr. 8)				4 anu	Beneficia Owned F	Beneficially Owned Following Reported		r Indirect str. 4)	Beneficial Ownership (Instr. 4)			
									Code	v	Amount	(A) (D)	or F	rice	Transaction(s) (Instr. 3 and 4)				(III3ti. 4)	
Common	Stock			01/0	2/201	2/2019		A		2,270 ⁽¹⁾ A		4	\$0.00	0 4,540 ⁽²⁾) D				
Common Stock													2,412,001		I		Held by ARCH Venture Fund VI, L.P. ⁽³⁾			
Common Stock														1,53	7,684		I	Held by ARCH Venture Fund VIII Overage, L.P. ⁽⁴⁾		
			Table II -								sed of, onvertib				Owned					
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	3A. Deemed Execution E if any (Month/Day	Date,	Code (Instr.		Derivative I		6. Date Exercis Expiration Date (Month/Day/Yea			7. Title and Amount of Securities Underlying Derivative Se (Instr. 3 and			8. Price of Derivative Security (Instr. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s	e s Illy	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership t (Instr. 4)	
					Code	v	(A)		Date Exercisab		Expiration Date	Title	or Nu of	nount mber ares		(Instr. 4)	Sin(3)			
Stock Option (Right to Buy)	\$18.24	01/02/2019			A		7,900 ⁽⁵⁾		12/31/201	19 (01/02/2029	Commo Stock	n 7,	900	\$0.00	7,900)	D		

- 1. Represents restricted stock units that vest as to 100% of the shares on December 31, 2019 granted pursuant to the Quanterix Corporation 2018 Non-Employee Director Compensation Policy. Each restricted stock unit represents the right to receive one share of the Issuer's common stock.
- 2. Includes 2,270 restricted stock units.
- 3. The shares are held of record by ARCH Venture Fund VI, L.P. ("ARCH Fund VI"). The sole general partner of ARCH Fund VI is ARCH Venture Partners VI, L.P. ("ARCH Partners VI"), which may be deemed to beneficially own the shares held by ARCH Fund VI. The sole general partner of ARCH Partners VI is ARCH Venture Partners VI, LLC ("ARCH VI LLC"), which may be deemed to beneficially own the shares held by ARCH Fund VI. ARCH Partners VI and ARCH VI LLC disclaim beneficial ownership of such shares, except to the extent of any pecuniary interest therein. As a managing director of ARCH VI LLC, the Reporting Person may be deemed to beneficially own the shares held by ARCH Fund VI. The Reporting Person disclaims beneficial ownership of such shares, except to the extent of any pecuniary interest therein.
- 4. The shares are held of record by ARCH Venture Fund VIII Overage, L.P. ("ARCH Fund Overage"). The sole general partner of ARCH Fund Overage is ARCH Venture Partners VIII, LLC ("ARCH VIII LLC"), which may be deemed to beneficially own the shares held by ARCH Fund Overage. ARCH VIII LLC disclaims beneficial ownership of such shares, except to the extent of any pecuniary interest therein. As a managing director of ARCH VIII LLC, the Reporting Person may be deemed to beneficially own the shares held by ARCH Fund Overage. The Reporting Person disclaims beneficial ownership of such shares, except to the extent of any pecuniary interest therein.
- 5. Represents non-qualified stock option to purchase 7,900 shares of common stock granted pursuant to the Quanterix Corporation 2018 Non-Employee Director Compensation Policy.

Remarks:

/s/ Brian Keane, as Attorney-in-Fact for Keith L. Crandell

** 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 I	Number.