FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

l	OMB APPRO	VAL				
	OMB Number:	3235-0287				
l	Estimated average burde	en				
l	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(b)

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 an	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) 07/01/2019									Offic belov	er (give title v)		Other below)	specify
(Street) CHICAGO IL 60631				4. If <i>A</i>	4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Indiv Line) X	vidual or Joint/Group Filing (Check Applicable Form filed by One Reporting Person Form filed by More than One Reporting Person				on	
(City)	(St		Zip)																
	Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																		
1. Title of Security (Instr. 3) 2. Transact Date (Month/Day					y/Year) Ex		2A. Deemed Execution Date, if any (Month/Day/Year)		3. Transaction Code (Instr. 8)		4. Securities Acquired (A) of Disposed Of (D) (Instr. 3, 4 a			and 5) Securities Beneficiall		ities cially d Following	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)		7. Nature of Indirect Beneficial Ownership
							·		Code	v	Amount	(A) oi (D)	Pric	Transaction(s)				(Instr. 4)	
Common	Stock			07/01/2	.019				A		332(1)	A	\$3	5.38 ⁽²⁾	5,374(3)		D		
Common Stock														2,412,001		I		Held by ARCH Venture Fund VI, L.P. ⁽⁴⁾	
Common Stock															1,5	537,684]	Ī	Held by ARCH Venture Fund VIII Overage, L.P. ⁽⁵⁾
		Та	ıble II -								osed of, o				wned				
1. Title of Derivative Security (Instr. 3)	of 2. ve Conversion Date Execution Date, T or Exercise (Month/Day/Year)			ransaction		5. Number of		6. Date Exerci Expiration Dat (Month/Day/Ye		isable and te	7. Title and Amount of Securities Underlying Derivative Security (Instr. and 4)		8. Price of Derivative Security (Instr. 5)		9. Number of derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	Owi Fori Dire or Ii (I) (I	nership m: ect (D) ndirect nstr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)	
Evaluation				c	Code	v	(A)	(D)	Date Exercis	able	Expiration Date	Title	Amour or Number of Shares	r					

- 1. Consists of stock granted in lieu of cash fees for service on the Company's Board of Directors and committees thereof for the second quarter of 2019, pursuant to an election made in December 2018.
- 2. Closing price of the Company's common stock on the Nasdaq Global Market on July 1, 2019.
- 3. Includes 2,270 restricted stock units.

4. 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.

5. 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.

Remarks:

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

07/03/2019

** 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.