FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

	Washington, D.C. 20549						
Check this box if no longer subject to Section 16. Form 4 or Form 5 obligations may continue. See Instruction 1(b).	STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP						
	Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934						

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

5. Relationship of Reporting Person(s) to Issuer

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

2. Issuer Name and Ticker or Trading Symbol

1. Name and Address of Reporting Person* CRANDELL KEITH					2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]								Relationship of Reporting Person(s) to Issuer (Check all applicable) X Director X 10% Owner Officer (give title Other (specify					
(Last) (First) (Middle) 8755 WEST HIGGINS ROAD, SUITE 1025				3. Date of Earliest Transaction (Month/Day/Year) 12/11/2017								below)	IVC UUC		below)	роспу		
(Street)	GO I	L	60631		4. If Amendment, Date of Original Filed (Month/Day/Year)								Individual or Joint/Group Filing (Check Applicable Line) X Form filed by One Reporting Person Form filed by More than One Reporting Person					
(City)	(State)	(Zip)															
			Table I - Nor	n-Deriv	ative \$	Secur	ities Acc	uired,	Dis	posed of, o	or Bene	ficially C	Owned					
1. Title of Security (Instr. 3)		2. Transaction Date (Month/Day/Year)		2A. Deemed Execution Date, if any (Month/Day/Year)	3. Transaction Code (Instr. 8)		4. Securities Disposed Of			Securities Beneficially Following F	5. Amount of Securities Beneficially Owned Following Reported Fransaction(s)	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Direct ndirect r. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)				
								Code	v	Amount	(A) or (D)	Price	(Instr. 3 and	i 4)				
Common	Stock			12/11/	/2017			С		398,257	A	(1)	398,2	257		I :	Held by ARCH Venture Fund VI, L.P. ⁽²⁾	
Common	Stock			12/11/	/2017			С		1,045,426	A	(1)	1,443,	683		I :	Held by ARCH Venture Fund VI, L.P. ⁽²⁾	
Common	Stock			12/11/	/2017			С		477,042	A	(1)	1,920,	725		I :	Held by ARCH Venture Fund VI, L.P. ⁽²⁾	
Common	Stock			12/11/	/2017			С		109,769	A	(1)	2,030,	494		I :	Held by ARCH Venture Fund VI, L.P. ⁽²⁾	
Common	Stock			12/11/	/2017			С		381,507	A	(1)	2,412,	001		I :	Held by ARCH Venture Fund VI, L.P. ⁽²⁾	
Common	Stock			12/11/	/2017			С		1,271,684	A	(1)	1,271,	684		I :	Held by ARCH Venture Fund VIII Overage, L.P. ⁽³⁾	
Common	mmon Stock			12/11/	1/2017			P		266,000(4)	A	\$15	1,537,684			I ;	Held by ARCH Venture Fund VIII Overage, L.P. ⁽³⁾	
										osed of, or onvertible			vned					
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)	4. Transa Code (action (Instr.	5. Numb Derivati Securiti Acquire Dispose	per of ve		Exercis	sable and 7. e So ar) D	Title and	Amount of Inderlying Security	8. Price of Derivative Security (Instr. 5)	9. Numb derivativ Securiti Benefici Owned Followir	ve es ially ng	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership (Instr. 4)	
				Code	v	(A)	(D)	Date Exercisa		Expiration Date Ti		Amount or Number of Shares	or Rep		tion(s)			

Derivative C Security (Instr. 3) F	2. Conversion or Exercise Price of	3. Transaction Date	3A. Deemed					Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities)													
	Derivative Security			ive ies ed (A) or ed of (D)	Expiration Date (Month/Day/Year) (A) or of (D)			d Amount of Underlying Security nd 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)								
Series A-1 Preferred Stock	(1)	12/11/2017		С			1,280,000	(5)	(6)	Common Stock	398,257	(1)	0	I	Held by ARCH Venture Fund VI, L.P. ⁽²⁾						
Series A-2 Preferred Stock	(1)	12/11/2017		С			3,360,000	(5)	(6)	Common Stock	1,045,426	(1)	0	I	Held by ARCH Venture Fund VI, L.P. ⁽²⁾						
Series B Preferred Stock	(1)	12/11/2017		С			1,533,214	(5)	(6)	Common Stock	477,042	(1)	0	I	Held by ARCH Venture Fund VI, L.P. ⁽²⁾						
Warrant to Purchase Series C Preferred Stock (right to buy) ⁽⁷⁾	\$3.3299	11/27/2017		X ⁽⁸⁾			18,450	03/01/2012	11/30/2017	Series C Preferred Stock	18,450	(8)	14,761	I	Held by ARCH Venture Fund VI, L.P. ⁽²⁾						
Series C Preferred Stock ⁽⁷⁾	(1)	11/27/2017		X ⁽⁸⁾		18,450		(5)	(6)	Common Stock	5,740	(1)	363,869	I	Held by ARCH Venture Fund VI, L.P. ⁽²⁾						
Series C Preferred Stock ⁽⁷⁾	(1)	11/27/2017		S ⁽⁸⁾			14,350	(5)	(6)	Common Stock	4,464	\$4.2815 ⁽⁸⁾	349,519	I	Held by ARCH Venture Fund VI, L.P. ⁽²⁾						
Warrant to Purchase Series C Preferred Stock (right to buy) ⁽⁷⁾	\$3.3299	11/27/2017		X ⁽⁹⁾			11,993	07/30/2012	11/30/2017	Series C Preferred Stock	11,993	(9)	2,768	I	Held by ARCH Venture Fund VI, L.P. ⁽²⁾						
Series C Preferred Stock ⁽⁷⁾	(1)	11/27/2017		X ⁽⁹⁾		11,993		(5)	(6)	Common Stock	3,731	(1)	361,512	I	Held by ARCH Venture Fund VI, L.P. ⁽²⁾						
Series C Preferred Stock ⁽⁷⁾	(1)	11/27/2017		S ⁽⁹⁾			9,328	(5)	(6)	Common Stock	2,902	\$4.2815 ⁽⁹⁾	352,184	I	Held by ARCH Venture Fund VI, L.P. ⁽²⁾						
Warrant to Purchase Series C Preferred Stock (right to buy) ⁽⁷⁾	\$3.3299	11/27/2017		X ⁽¹⁰⁾			2,768	11/05/2012	11/30/2017	Series C Preferred Stock	2,768	(10)	0	I	Held by ARCH Venture Fund VI, L.P. ⁽²⁾						
Series C Preferred Stock ⁽⁷⁾	(1)	11/27/2017		X ⁽¹⁰⁾		2,768		(5)	(6)	Common Stock	861	(1)	354,952	I	Held by ARCH Venture Fund VI, L.P. ⁽²⁾						
Series C Preferred Stock ⁽⁷⁾	(1)	11/27/2017		S ⁽¹⁰⁾			2,152	(5)	(6)	Common Stock	669	\$4.2815 ⁽¹⁰⁾	352,800	I	Held by ARCH Venture Fund VI, L.P. ⁽²⁾						
Series C Preferred Stock	(1)	12/11/2017		С			352,800	(5)	(6)	Common Stock	109,769	(1)	0	I	Held by ARCH Venture Fund VI, L.P. ⁽²⁾						
Series D Preferred Stock	(1)	12/11/2017		С			1,226,158	(5)	(6)	Common Stock	381,507	(1)	0	I	Held by ARCH Venture Fund VI, L.P. ⁽²⁾						
Series D Preferred Stock	(1)	12/11/2017		С			4,087,193	(5)	(6)	Common Stock	1,271,684	(1)	0	I	Held by ARCH Venture Fund VIII Overage, L.P. ⁽³⁾						

Explanation of Responses:

^{1.} The shares of preferred stock automatically converted into common stock immediately prior to the closing of the Issuer's initial public offering using a calculation defined in the Issuer's Restated Certificate of Incorporation, as amended, based on the initial purchase price and the conversion rate in effect at the time of conversion. The shares of preferred stock automatically converted on a 1-for-3.214 basis and such conversion rate is reflected in the amount of common stock underlying the security.

^{2.} 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.

- 3. 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.
- 4. Reflects shares purchased in the Issuer's initial public offering.
- 5. These shares were convertible into the Issuer's common stock in accordance with the Issuer's Restated Certificate of Incorporation, as amended, at any time after the issuance of such shares, at the holder's election.
- Not applicable
- 7. This transaction, which occurred prior to the Issuer's initial public offering of Common Stock, is being reported as required by Rule 16a-2(a) under the Securities Exchange Act of 1934, as amended.
- 8. These warrants to purchase Series C preferred stock were exercised on a cashless basis for shares of Series C preferred stock, resulting in the Issuer's withholding of 14,350 shares of Series C preferred stock to pay the exercise price and issuing to the reporting person the remaining 4,100 shares of Series C preferred stock. The Issuer's withholding of Series C preferred stock was at a price per share of Series C preferred stock of \$4.2815, which is equal to approximately \$13.76 per share on an as-converted-to-common stock basis.
- 9. These warrants to purchase Series C preferred stock were exercised on a cashless basis for shares of Series C preferred stock, resulting in the Issuer's withholding of 9,328 shares of Series C preferred stock to pay the exercise price and issuing to the reporting person the remaining 2,665 shares of Series C preferred stock. The Issuer's withholding of Series C preferred stock was at a price per share of Series C preferred stock of \$4.2815, which is equal to approximately \$13.76 per share on an as-converted-to-common stock basis.
- 10. These warrants to purchase Series C preferred stock were exercised on a cashless basis for shares of Series C preferred stock, resulting in the Issuer's withholding of 2,152 shares of Series C preferred stock to pay the exercise price and issuing to the reporting person the remaining 616 shares of Series C preferred stock. The Issuer's withholding of Series C preferred stock was at a price per share of Series C preferred stock of \$4.2815, which is equal to approximately \$13.76 per share on an as-converted-to-common stock basis.

/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 Number.