FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

## STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

l	OMB APPROVAL								
	OMB Number:	3235-0287							
	Estimated average burde	en							
	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

	nd Address of David C.		2. Issuer Name <b>and</b> Ticker or Trading Symbol  Quanterix Corp [ QTRX ]									eck all appli Directo	*		on(s) to Issuer  10% Owner Other (specify				
•	t) (First) (Middle)  QUANTERIX CORPORATION,  HARTWELL AVENUE						3. Date of Earliest Transaction (Month/Day/Year) 12/16/2019									SVP R&D and CTO			
(Street) LEXINGTON MA 02421						4. If Amendment, Date of Original Filed (Month/Day/Year)									e) X Form f Form f	ual or Joint/Group Filing (Check Applicable Form filed by One Reporting Person Form filed by More than One Reporting Person			
(City)	(S		(Zip)																
Table I - Non-Deriva  1. Title of Security (Instr. 3)  2. Transac Date (Month/Da					ction	tion 2A. Deen Executio			3. Trans Code	, Dis	4. Securities Acquired (A) Disposed Of (D) (Instr. 3, 4			(A) or	5. Amou Securitie Benefici	int of es ally Following	Form (D) o	: Direct r Indirect str. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
						Code	v	Amount	(1	A) or D)	Price	Transac (Instr. 3	tion(s)			(111341. 4)			
Common Stock 12/16/					/2019	2019		M		1,667	7	A	\$1.93	3 46	46,427		D		
Common Stock 12/				12/16	5/2019				S		1,667	(1)	D	\$24.1	5 44,	44,760(2)		D	
		Т	able II -								osed of converti				Owned				
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security		3A. Deem Executior if any (Month/Da	Date, Transac Code (I					6. Date E Expiratio (Month/D	n Date	•	7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)			8. Price of Derivative Security (Instr. 5)	9. Number derivative Securities Beneficial Owned Following Reported Transactio (Instr. 4)	Owners Form: Direct (I or Indire (I) (Instr	Ownership	Beneficial Ownership (Instr. 4)
					Code	ode V		(D)	Date Exercisa		expiration Date	Title	O N O	Amount or Number of Shares					
Stock Option (Right to Buy)	\$1.93	12/16/2019			M			1,667	(3)	(	07/06/2021	Comm		1,667	\$0.00	1,941		D	

## **Explanation of Responses:**

- 1. The sale of shares reported on this Form 4 was made pursuant to a Rule 10b5-1 trading plan adopted on May 16, 2019.
- 2. Includes 8,949 restricted stock units.
- 3. Option is fully vested.

## Remarks:

/s/ Brian Keane, Attorney-in-

12/16/2019

fact

\*\* 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.