FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION Washington, D.C. 20549

OMB APPROVAL

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

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

0									
OMB Number:	3235-0287								
Estimated average burden									
hours per response:	0.5								

1 Indiric and Address of Reporting Leson			2. Issuer Name and Ticker or Trading Symbol <u>Quanterix Corp</u> [QTRX]	5. Relationship of Reporting Person(s) to Issuer (Check all applicable) Director 10% Owner Officer (give title Other (specify
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 113 HARTWELL AVENUE		,	3. Date of Earliest Transaction (Month/Day/Year) 03/01/2019	X below) below) Sr VP, Stgc Mrktg & Assay Tech
(Street) LEXINGTON (City)	MA (State)	02421 (Zip)	4. If Amendment, Date of Original Filed (Month/Day/Year)	 6. Individual or Joint/Group Filing (Check Applicable Line) X Form filed by One Reporting Person Form filed by More than One Reporting Person

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)	Transaction Code (Instr.		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 and 5)			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)		(1150.4)
Common Stock	03/01/2019		S		75(1)	D	\$25.0043 ⁽²⁾	16,715 ⁽³⁾	D	

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

				1		1	_	1		1		1	1	r	
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 (8)		5. Nu of Deriv Secu Acqu (A) or Dispo of (D) (Instr and 5	ative rities ired osed	Expiration Date (Month/Day/Year) (Month/Day/Year) (Month/Day/Year)		and 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 Transaction(s) (Instr. 4)		11. Nature of Indirect Beneficial Ownership (Instr. 4)
				Code	v	(A)	(D)	Date Exercisable	Expiration Date	Title	Amount or Number of Shares				

Explanation of Responses:

1. The sale of shares reported on this Form 4 was made solely to cover tax obligations upon the vesting of 208 RSUs and was effected pursuant to a Rule 10b5-1 trading plan adopted on December 14, 2018. 2. This transaction was executed in multiple trades at prices ranging from \$24.85 to \$25.0738. The price reported reflects the weighted average sale price. The reporting person undertakes to provide upon request to the SEC staff, the issuer or any security holder of the issuer full information regarding the number of shares and prices at which the transaction was effected. 3. Includes 14,972 shares of restricted stock units.

Remarks:

<u>/s/ Brian Keane, as Attorney-</u> <u>in-Fact</u>

03/01/2019

** 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.