SEC Form 4	
------------	--

 \square

FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

	Check this box if no longer subject to
1	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 or Section 30(h) of the Investment Company Act of 1940

OMB Number:	3235-0287										
Estimated average b	urden										

Estimated average burden	
hours per response:	0.5

1 I Marile and Address of Reporting Leson			2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]	5. Relationship of Reporting Person(s) to Issuer (Check all applicable)					
				X	Director	10% Owner			
 (Last) (First) (Middle)		(Middle)	3. Date of Earliest Transaction (Month/Day/Year)	X	Officer (give title below)	Other (specify below)			
C/O QUANTERIX CORPORATION,			06/04/2019		Chairman, President & CEO				
900 MIDDLESEX TURNPIKE									
(Street)			4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Individual or Joint/Group Filing (Check App Line)					
BILLERICA	МА	01821		X	Form filed by One Report	rting Person			
					Form filed by More than Person	One Reporting			
(City)	(State)	(Zip)							

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)	3. Transa Code (8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 and 5)		5. Amount of Securities Beneficially Owned Following Reported	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)		(1130.4)
Common Stock	06/04/2019		S		203(1)	D	\$25.81	1,180,065	D	
Common Stock	06/04/2019		S		2,870 ⁽¹⁾	D	\$26.0093 ⁽²⁾	1,177,195 ⁽³⁾	D	

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

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)		n of		6. Date Exerc Expiration Da (Month/Day/Y	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)	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)
				Code	v			Date Exercisable	Expiration Date	Title	Amount or Number of Shares			

Explanation of Responses:

1. The sales of shares reported on this Form 4 was made solely to cover tax obligations upon the vesting of 661 RSUs and 6,250 RSUs, respectively.

2. This transaction was executed in multiple trades at prices ranging from \$26.00 to \$26.05. 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 39,815 shares of common stock that are subject to performance based vesting and 222,177 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorney-

06/05/2019

** Signature of Reporting Person

in-Fact

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.

Date