SEC Form 4

П

FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

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 or Section 30(h) of the Investment Company Act of 1940

OMB APPROVAL OMB Number: 3235-0287 Estimated average burden

Estimated average burden	
hours per response:	0.5

I I Name and Address of Reporting Feison			2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]	5. Relationship of Reporting Person(s) to Issuer (Check all applicable) Director 10% Owner					
				x	Officer (give title	Other (specify			
(Last) (First) (Middle)		(Middle)	3. Date of Earliest Transaction (Month/Day/Year)		below)	below)			
C/O QUANTERIX CORPORATION,			06/05/2018		SVP Finance and Corporate Dev				
113 HARTWELL AVENUE									
,			4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Indiv Line)	idual or Joint/Group Filin	g (Check Applicable			
(Street)		02.421		X	Form filed by One Rep	orting Person			
LEXINGTON	MA	02421			Form filed by More tha Person	n One Reporting			
(City)	(State)	(Zip)			reison				

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.					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 Pri		Price	Transaction(s) (Instr. 3 and 4)		(1150. 4)
Common Stock	06/05/2018		М		25,000	A	\$2.28	78,802(1)	D	
Common Stock	06/05/2018		S		25,000	D	\$17.1453	53,802 ⁽¹⁾	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)		of		6. Date Exercisable and Expiration Date (Month/Day/Year)		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	(A)	(D)	Date Exercisable	Expiration Date	Title	Amount or Number of Shares				
Stock Option (Right to Buy)	\$2.28	06/05/2018		М			25,000	(2)	02/03/2022	Common Stock	25,000	\$0.00	21,670	D	

Explanation of Responses:

1. Includes 7,131 restricted stock units that vest as to 25% of the shares on May 10, 2019 and as to an additional 2.083% of the shares per month for 36 months thereafter.

2. Vests 23,335 share(s) on 02-Jan-2013, 1,944 share(s) on 31-Jan-2013, 1,944 share(s) on 28-Feb-2013, 1,944 share(s) on 31-Mar-2013, 1,944 share(s) on 30-Jun-2013, 1,944 share(s) on 31-Juh-2013, 1,944 share(s) on 31-Juh-2014, 1,944 share(s) on 31-Juh-2015, 1,944 share(s) on 31-

/s/ Brian Keane, as Attorney-in-06/05/2018

<u>Fact</u>

** Signature of Reporting Person Date

Reminder: Report on a separate line for each class of securities beneficially owned directly or indirectly.

 \ast 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.