SEC F	orm 4
-------	-------

 \square

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

1. Name and Address of Reporting Person* <u>Duffy David C.</u>			2. Issuer Name and Ticker or Trading Symbol <u>Quanterix Corp</u> [QTRX]	(Check	tionship of Reporting Per all applicable) Director Officer (give title	son(s) to Issuer 10% Owner Other (specify
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 113 HARTWELL AVENUE		()	3. Date of Earliest Transaction (Month/Day/Year) 10/31/2018	X	below) SVP R&D and	below)
(Street) LEXINGTON (City)	MA (State)	02421 (Zip)	4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Indiv Line) X	idual or Joint/Group Filin Form filed by One Rep Form filed by More tha Person	orting 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)	3. Transaction Code (Instr. 8)					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)		(1150.4)
Common Stock	10/31/2018		М		1,354	Α	\$1.93	37 , 952 ⁽¹⁾	D	
Common Stock	10/31/2018		М		1,860	Α	\$2.28	39,8 12 ⁽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 Deri Sec Acq (A) o Disp of (I	oosed D) tr. 3, 4	Expiration Date (Month/Day/Year) ed ed		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)	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	10/31/2018		М			1,860	(2)	02/03/2022	Common Stock	1,860	\$0.00	29,253	D	
Stock Option (Right to Buy)	\$1.93	10/31/2018		М			1,354	(3)	07/06/2021	Common Stock	1,354	\$0.00	14,202	D	

Explanation of Responses:

1. Includes 5,485 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. 25% @ 1 year, monthly on last day of month thereafter for 3 years (starting in 12th month) beginning on 01-Nov-2011

3. 25% @ 1 year, monthly on last day of month thereafter for 3 years (starting in 12th month) beginning on 06-Jul-2011

Remarks:

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

<u>fact</u>

10/31/2018

** 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.