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

П

FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

Check this box if no longer subject	ct 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

OND NUMBER.	0200 0201
Estimated average burd	len
hours per response:	0.5

1. Name and Address of Reporting Person* Duffy David C. (Last) (First) (Middle) C/O QUANTERIX CORPORATION, 113 HARTWELL AVENUE		n*	2. Issuer Name and Ticker or Trading Symbol <u>Quanterix Corp</u> [QTRX]		tionship of Reporting Perso all applicable) Director	on(s) to Issuer 10% Owner
			3. Date of Earliest Transaction (Month/Day/Year) 11/18/2019	- X	Officer (give title below) SVP R&D and	Other (specify below) CTO
(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 Filing Form filed by One Repor Form filed by More than Person	ting 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. 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)		(1150.4)
Common Stock	11/18/2019		М		1,667	A	\$1.93	46,427	D	
Common Stock	11/18/2019		S		1,667(1)	D	\$22.5643 ⁽²⁾	44,760 ⁽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)		of Deri Secu Acq (A) o Disp of (E (Inst	5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		ate	Amount of		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	4			
Stock Option (Right to Buy)	\$1.93	11/18/2019		М			1,667	(4)	07/06/2021	Common Stock	1,667	\$0.00	3,608	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. This transaction was executed in multiple trades at prices ranging from \$22.53 to \$22.58. 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 9,063 restricted stock units.

4. Option is fully vested.

Remarks:

/s/ Brian Keane, Attorney-in-

<u>fact</u>

11/19/2019

** 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.