\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

OMB APPROVAL OMB Number: 3235-0287 Estimated average burden 0.5

			or Section 30(h) of the Investment Company Act of 1940						
1	dress of Reporting	Person [*]	2. Issuer Name and Ticker or Trading Symbol <u>Quanterix Corp</u> [QTRX]		5. Relationship of Reporting Person(s) to Issuer (Check all applicable)				
Duffy David C.					Director	10% Owner			
			Dete of Sectors they (Marth Dev/Mart)	x	Officer (give title below)	Other (specify below)			
(Last) C/O QUANT	(First) ERIX CORPO	(Middle) RATION,	3. Date of Earliest Transaction (Month/Day/Year) 12/18/2019		SVP R&D and CTO				
113 HARTW	ELL AVENUE								
(Stroot)			4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Indiv Line)	/idual or Joint/Group Fil	ing (Check Applicable			
(Street)	N MA	02421		X	Form filed by One Re	eporting Person			
	N IV17 1	02421			Form filed by More the Person	nan 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)	if any 🧴						5. Amount of Securities Beneficially Owned Following Reported		7. Nature of Indirect Beneficial Ownership (Instr. 4)
			Code	v	Amount	(A) or (D)	Price	Transaction(s) (Instr. 3 and 4)		(iiisti: 4)
Common Stock	12/18/2019		М		6,600	Α	\$2.28	51,360 ⁽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	12/18/2019		М			6,600	(2)	02/03/2022	Common Stock	6,600	\$0.00	24,513	D	

Explanation of Responses:

1. Includes 8,949 restricted stock units.

2. Option is fully vested.

Remarks:

/s/ Brian Keane, Attorney-in-

12/20/2019

** Signature of Reporting Person Date

<u>fact</u>

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.

hours per response: