FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL									
OMB Number: 3235-028									
Estimated average burden									
hours per response	e: 0.5								

Check this box if no longer subject to Section 16. Form 4 or Form 5 obligations may continue. See Instruction 1(b).

Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934 or Section 30(h) of the Investment Company Act of 1940

Name and Address of Reporting Person* Duffy David C.						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]								(Chec	k all app Direc	,	ng Per	rson(s) to Is 10% O Other (wner
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						3. Date of Earliest Transaction (Month/Day/Year) 05/08/2020								X		below) SVP R&D		below)	
(Street) BILLERICA MA 01821 (City) (State) (Zip)					4. If A	4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Indi	·				
		Table	I - No	n-Deriva	tive S	Secu	rities	Acq	juired,	, Dis	posed of	, or E	Benef	icially	y Own	ed			
1. Title of Security (Instr. 3) 2. Transacti Date (Month/Day)					Execution Date,				es Acquired (A Of (D) (Instr. 3,			Securi Benefi	urities eficially ned Following		vnership n: Direct r Indirect nstr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)			
									Code	v	Amount	(A) (D)	or Pi	rice	Transa	action(s) 3 and 4)			(311 4)
Common Stock 05/08/2					2020			S		1,667(1)	1) \$	26.52	51,192(2)			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) 3. Deemed Execution Date, if any (Month/Day/Year)		Code (8)	ransaction of Code (Instr. Derivative		6. Date Exercisable and Expiration Date (Month/Day/Year) Date Expiration Exercisable Date		7. Title and Amount of Securities Underlying Derivative Security (Inst 3 and 4) Amount or Number of Title Shares		De Se (In	Price of rivative curity str. 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)				

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 December 3, 2019.
- 2. Includes 13,237 restricted stock units.

Remarks:

/s/ Brian Keane, Attorney-in-

05/12/2020

fact

** 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.