FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL											
OMB Number: 3235-0287											
Estimated average burden											
hours per response:	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

1. Name and Address of Reporting Person* Duffy David C.						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]								(Checl	all app Direc	ionship of Reporti all applicable) Director Officer (give title		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) 03/19/2021								X	below				,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,
(Street) BILLER (City)	4. If A	4. If Amendment, Date of Original Filed (Month/Day/Year)							6. Indir Line) X	·									
		Table	I - No	n-Deriva	tive S	Secu	rities	Acq	uired,	Dis	posed of	, or E	Benefi	cially	Own	ed			
1. Title of Security (Instr. 3) 2. Transacti Date (Month/Day)					Execution Date,				es Acquired (A Of (D) (Instr. 3,		4 and Secu Bene		cially Following	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)		7. Nature of Indirect Beneficial Ownership (Instr. 4)			
				Code	v	Amount	(A) (D)	or Prid	се	Transa	nsaction(s) str. 3 and 4)			, , ,					
Common Stock 03/19/20					2021			S		1,667(1)	7 ⁽¹⁾ D §		8.36	38,285(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)		4. Transaction Code (Instr. 8) Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5) Code V (A) (D)		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) Amour or Numbe of Title Shares		Der Sec (Ins	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 May 15, 2020.
- 2. Includes 11,313 restricted stock units.

Remarks:

/s/ Brian Keane, Attorney-in-

fact

** Signature of Reporting Person

Date

03/19/2021

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.