FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington,	D.C.	20549
vvasimigton,	D.C.	20040

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

											_										
Name and Address of Reporting Person* Duffy David C.						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]										eck all applic Directo	,		son(s) to Iss 10% Ov Other (s	vner	
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						3. Date of Earliest Transaction (Month/Day/Year) 01/03/2022										below)		D and	below)		
(Street) BILLER			01821		4. If Amendment, Date of Original Filed (Month/Day/Year)								ar)	Line	Individual or Joint/Group Filing (Check Applicable ine) X Form filed by One Reporting Person Form filed by More than One Reporting Person						
(City)	(S	tate)	(Zip)																		
			ole I - Nor			_				i, D	isp								1		
Date				Date	saction n/Day/Year)		2A. Deemed Execution Date, if any (Month/Day/Year)		e, Trar Cod	Code (Instr						Securitie Benefici Owned F	5. Amount of Securities Beneficially Owned Following		n: Direct or Indirect nstr. 4)	7. Nature of Indirect Beneficial Ownership	
									Cod	e v		Amount (A		(A) or (D)	Price	Reported Transact (Instr. 3	tion(s)			(Instr. 4)	
Common Stock					3/2022			М			4,667		A	\$2.9	5 27	,949		D			
Common	on Stock			01/03	01/03/2022				М			15,556		A	\$2.8	9 43	43,505		D		
Common	Common Stock 01/			01/03	3/2022		М			21,023 A		Α	\$2.2	64,5	64,528(1)		D				
		-	Table II - I									sed of, onvertil				Owned					
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	3A. Deemed Execution D if any (Month/Day/	Date, T	ransaction Code (Instr.		5. Number of		6. Date Exercis Expiration Date (Month/Day/Yea		ate	e oʻ ar) U D		7. Title and Amor of Securities Underlying Derivative Secur (Instr. 3 and 4)		8. Price of Derivative Security (Instr. 5)	9. Number derivative Securities Beneficial Owned Following Reported Transactio (Instr. 4)	Owners Form: Direct (or Indir	Ownership	Beneficial Ownership (Instr. 4)	
					Code	v	(A)	(D)	Date Exercis	able		kpiration ate	Title		Amount or Number of Shares						
Stock Option (right to buy)	\$2.96	01/03/2022			М			4,667	(2)		07	7/08/2024		nmon ock	4,667	\$0.00	0.00		D		
Stock Option (right to buy)	\$2.89	01/03/2022			М			15,556	(2)		05	5/21/2023		nmon ock	15,556	\$0.00	0.00		D		
Stock Option (Right to buy)	\$2.28	01/03/2022		T	M			21,023	(2)		02	2/03/2022		nmon ock	21,023	\$0.00	0.00		D		

Explanation of Responses:

- 1. Includes 7,653 restricted stock units.
- 2. All shares underlying this option have vested.

Remarks:

/s/ Brian Keane, Attorney-in-

01/05/2022

lact

** 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.