FORM 4

Check this box if no longer subject to Section 16. Form 4 or Form 5 UNITED STATES SECURITIES AND EXCHANGE COMMISSION Washington, D.C. 20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL

-	-								
OMB Number:	3235-0287								
Estimated average burden									
hours per response:	0.5								

Instruction 1(b).	munue. See		Filed pursuant to Section	led pursuant to Section 16(a) of the Securities Exchange Act of 1934								
				f the Investment Company Act of 1940	1004							
1. Name and Address Duffy David C	1 0	rson*	2. Issuer Name ar Quanterix Co	nd Ticker or Trading Symbol orp [QTRX]		tionship of R all applicabl Director Officer (giv	10% Owner		er			
(Last) C/O QUANTERI 900 MIDDLESE)			3. Date of Earliest 07/14/2021	Transaction (Month/Day/Year)		below) be SVP R&D and CT(below) CTO				
			4. If Amendment,	Date of Original Filed (Month/Day/Year)		idual or Join	t/Group Filing	(Check App	licable			
(Street) BILLERICA	MA	01821			Line) X		by One Repo by More than	0				
(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)	2A. Deemed Execution Date, if any (Month/Day/Year)						Securities Beneficially	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)		(1130.4)
Common Stock	07/14/2021		S		1,667(1)	D	\$57.77	31,6 17 ⁽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)	3A. Deemed Execution Date, if any (Month/Day/Year)	4. Transa Code (8)		5. Number n of		6. Date Exercisable and Expiration Date (Month/Day/Year)				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				

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 March 8, 2021.

2. Includes 9,849 restricted stock units.

Remarks:

/s/ Brian Keane, Attorney-in-

fact

07/14/2021

** 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.