Instruction 1(b).

FORM 4

## **UNITED STATES SECURITIES AND EXCHANGE COMMISSION**

Washington, D.	C. 20549
----------------	----------

Check this box if no longer subject	STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP
to Section 16. Form 4 or Form 5	
obligations may continue. See	

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

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 ]									(Che	ck all app Direc	,		10% O	
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE					3. Date of Earliest Transaction (Month/Day/Year) 05/28/2021									X	belov			below)	
(Street) BILLER (City)	ICA M.	A 0	1821 Zip)		4. If A	4. If Amendment, Date of Original Filed (Month/Day/Year)							6. Inc Line) X						
		Table	I - Nor	n-Deriva	tive S	Secu	rities	Acq	uired,	Dis	posed of	, or E	3ene	ficiall	y Own	ed			
1. Title of Security (Instr. 3)  2. Transact Date (Month/Day				Executacy/Year) if any		Deemed ecution Date, ny onth/Day/Year)				ies Acquired (A Of (D) (Instr. 3			Securit Benefic Owned	5. Amount of Securities Beneficially Owned Following Reported		vnership n: Direct r Indirect nstr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)		
								Code	v	Amount	(A) (D)	or F	rice	Transa	ction(s) 3 and 4)				
Common Stock 05/28/.				/2021			S		1,667(1)	1	D	\$51	34,951(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)	ve Conversion Date Execution Date, or Exercise (Month/Day/Year) if any			Code ( 8)	Transaction of Code (Instr. Derivative		rative rities ired r osed ) : 3, 4	6. Date Exercisable and Expiration Date (Month/Day/Year)  Date Expiration Exercisable Date			7. Title and Amount of Securities Underlying Derivative Security (Ins 3 and 4)		nt er		9. Number derivative Securities Beneficiall Owned Following Reported Transactio (Instr. 4)	у	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 March 8, 2021.
- 2. Includes 10,581 restricted stock units.

## Remarks:

/s/ Brian Keane, Attorney-in-

fact

\*\* Signature of Reporting Person

Date

06/02/2021

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.