FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

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

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

MENT OF CHANGES IN BENEFICIAL OWNERSHIP

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 ar	2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]									(Checl	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) 10/15/2021								X		SVP R&D and CTO			
(Street) BILLER (City)	ICA M	A 0	1821 Zip)		4. If A	4. If Amendment, Date of Original Filed (Month/Day/Year)							6. Indi Line) X	Form filed by One Reporting Person Form filed by More than One Reporting Person					
		Table	I - No	n-Deriva	tive S	Secu	rities	Acq	uired	, Dis	posed of	, or E	Benef	ficially	Own	ed			
1. Title of Security (Instr. 3) 2. Transacti Date (Month/Day					Execution Date,		Date,	3. Transaction Code (Instr. 8) 4. Securitie Disposed C				4 and Securi		cially I Following	Form (D) o	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)		
									Code	v	Amount	(A) or (D)		rice	Transaction(s) (Instr. 3 and 4)				(111501.4)
Common Stock 10/15/2				2021			S		1,667(1)	I) \$	51.47	.47 26,616(2)			D			
		Tal							,		osed of, convertib			•	Owne	d	,		
1. Title of Derivative Security (Instr. 3)	Derivative Conversion Date Execution Date, Gecurity or Exercise (Month/Day/Year) if any					ransaction of Code (Instr. Derivativ		rative rities iired r osed)	6. Date Exercisable and Expiration Date (Month/Day/Year)			7. Title and Amount of Securities Underlying Derivative Security (Ins 3 and 4)		Sec (Ins	Price of ivative curity str. 5)		Ownershi Form: Direct (D) or Indirect (I) (Instr. 4	Ownership	Beneficial Ownership t (Instr. 4)
					Code V		(A)	(D)	Date Exercisable		Expiration Date	Title	Numb of Share						

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 9,117 restricted stock units.

Remarks:

/s/ Brian Keane, Attorney-in-

10/19/2021

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.