FORM 4

Check this box if no longer subject

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

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

STATEMENT	OF CHANGES	IN BENEFICIAL	OWNERSHIP
O :/ \: = :\:	0. 0.17		O 1111E1 (O1 111

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

to Section 16. Form 4 or Form 5 obligations may continue. See Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934 or Section 30(h) of the Investment Company Act of 1940 Instruction 1(b).

1. Name and Address of Reporting Person* Roskey Mark T.						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]								Check	all app	ship of Reporting I applicable) irector fficer (give title		n(s) to Is 10% Ov Other (s	ner
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						3. Date of Earliest Transaction (Month/Day/Year) 03/13/2020								X	tor				
(Street) BILLERICA MA 01821 (City) (State) (Zip)						4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Individual or Joint/Group Filing (Check Applicable Line) X Form filed by One Reporting Person Form filed by More than One Reporting Person					
			I - I	Non-Deriva				_	<u>. </u>		•	<u> </u>		ially					
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Ye			Execution Date,		, Ti	3. Transaction Code (Instr. 8)					nd 5) See Be Ow		Securities Beneficially Dwned Following		Direct	7. Nature of Indirect Beneficial Ownership (Instr. 4)			
								С	ode	V A	Amount	(A) or (D)	Price		Reported Transaction(s) (Instr. 3 and 4)		(Instr. 4)		(111501.4)
Common	Stock			03/13/202	20			S		566(1)	D	D \$16.04		34 ⁽²⁾ 21,929 ⁽³⁾		D			
		Tal	ole	II - Derivati (e.g., pu							oosed of, convertib				Owne	t			
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	Exe if ar	Deemed cution Date, iy nth/Day/Year)	4. Transa Code 8)	(Instr.	5. Numof of Deriv. Securi Acqui (A) or Dispo of (D) (Instr. and 5	ative rities ired osed	Exp (Mo	oiration E	Year)	Amount of Securities Underlying Derivative Security (Ins 3 and 4)		nt er		9. Number derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	Ow Foi Dir or (I)	nership	11. Nature of Indirect Beneficial Ownership (Instr. 4)

Explanation of Responses:

- $1. \ The sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 1,807 RSUs.$
- 2. This transaction was executed in multiple trades at prices ranging from \$15.88 to \$16.25. The price reported reflects the weighted average sale price. The reporting person undertakes to provide upon request to the SEC staff, the issuer or any security holder of the issuer full information regarding the number of shares and prices at which the transaction was effected.
- 3. Includes 17,647 restricted stock units.

Remarks:

/s/ Brian Keane, Attorney-in-

fact

** Signature of Reporting Person Date

03/17/2020

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.