## FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

	9	,		

l	OMB APPROVAL									
l	OMB Number:	3235-0287								
l	Estimated average burden									
l	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

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

Name and Address of Reporting Person*     Roskey Mark T.						2. Issuer Name <b>and</b> Ticker or Trading Symbol Quanterix Corp [ QTRX ]								Check	all app	olicable)	g Person(s) to I: 10% ( Other	
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						3. Date of Earliest Transaction (Month/Day/Year) 12/02/2019								X	below) SVP, Commercial		below)	
(Street) BILLERICA MA 01821					4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)								Individual or Joint/Group Filing (Check Applicable Line)     X Form filed by One Reporting Person     Form filed by More than One Reporting Person				
(City)	(St	ate) (	Zip)												Peis	OH		
		Tabl	e I - N	Non-Deriv	ative	Seci	uritie	s Ac	quire	ed, Di	isposed o	f, or E	Benefici	ally (	Owne	ed		
1. Title of Security (Instr. 3)  2. Transaction Date (Month/Day/Ye					Execution Date,		·	3. Transaction Code (Instr. 8)  4. Securities Acq Disposed Of (D) (			Acquired (A) or (D) (Instr. 3, 4 and 5		I 5) Secu Bene		ficially d Following	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
										v	Amount	(A) or (D)	Price		Trans	action(s) 3 and 4)		(111501.4)
Common Stock 12/02/201					19			S		47(1)	D	\$25.06	29 <sup>(2)</sup>	<sup>2)</sup> 14,285 <sup>(3)</sup>		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)  3. Transaction Date (Month/Day/Year)  3. Transaction Date (Month/Day/Year)		4. Transa Code ( 8)		Str. Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		Expiration Date		7. Title and Amount of Securities Underlying Derivative Security (Instr. and 4)  Amour or Numbe of Title Shares		t r		9. Number of derivative Securities Beneficially Owned Following Reported Transaction ((Instr. 4)	Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)				

## **Explanation of Responses:**

- 1. The sales of shares reported on this Form 4 was made solely to cover tax obligations upon the vesting of 148 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$24.99 to \$25.13. 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 11,541 restricted stock units.

## Remarks:

/s/ Brian Keane, Attorney-infact 12/03/2019

\*\* 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.