Instruction 1(b).

FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

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

**OMB APPROVAL** OMB Number: 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*  Roskey Mark T.						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [ QTRX ]									eck all applic Directo	cable) or (give title	g Pers	10% Ov Other (s below)	/ner
•	Last) (First) (Middle) C/O QUANTERIX CORPORATION, 13 HARTWELL AVENUE					3. Date of Earliest Transaction (Month/Day/Year) 05/10/2018									,	SVP and GM A		,	nts
(Street) LEXINGTON MA 02421				4. 1	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					
(City)	(\$	State)	(Zip)												Persor	1			
		Tak	ole I - Nor	n-Deri	vativ	e Se	curities	s Acc	quired, [	Disp	osed o	f, or B	ene	ficiall	y Owned				
Dat			Date	. Transaction Date Month/Day/Year)		2A. Deemed Execution Date, if any (Month/Day/Year)		Transaction Disposed Code (Instr. 5)			ities Acquired (A) or d Of (D) (Instr. 3, 4 and			5. Amour Securitie Beneficia Owned F	Form (D) (collowing (I) (I)		n: Direct or Indirect nstr. 4)	7. Nature of Indirect Beneficial Ownership	
								Code	v	Amount	(A) or (D) Pi		Price	Reported Transact (Instr. 3 a	ion(s)			(Instr. 4)	
Common	Stock			05/1	0/201	.8			A		7,131	(1)		\$ <mark>0</mark>	7,131 <sup>(2)</sup> D				
		•	Table II -						ired, Di options						Owned				
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	e (Month/Day/Year)	3A. Deemee Execution I if any (Month/Day	Date,	4. Transaction Code (Instr 8)		n of		6. Date Exercisable and Expiration Date (Month/Day/Year)			7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)			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)	Beneficial Ownership (Instr. 4)
					Code	v	(A)		Date Exercisable		xpiration ate	Title	OI N Of	umber					
Stock Option (right to	\$19.12	05/10/2018			A		24,311		(3)	0	5/10/2028	Commo: Stock	2	4,311	\$0	24,311	1	D	

## **Explanation of Responses:**

- 1. Represents shares of common stock issuable upon vesting of restricted stock units. The restricted stock units vest as to 25% of the shares on May 10, 2019 and vest as to an additional 2.083% of the shares per month for 36 months thereafter.
- 2. Represents 7,131 restricted stock units.
- 3. The option vests as to 25% of the shares on May 10, 2019 and vests as to an additional 2.083% of the shares per month for 36 months thereafter.

/s/ Brian Keane, as Attorney-in-05/14/2018 **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.