FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

## STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL										
OMB Number:	3235-028									
Estimated average burden										

0.5

hours per response:

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

1. Name and Address of Reporting Person* <u>Roskey Mark T.</u>						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [ QTRX ]									eck all appli Directo	cable) or	g Pers	son(s) to Iss	vner	
_	O QUANTERIX CORPORATION,						3. Date of Earliest Transaction (Month/Day/Year) 03/12/2019									ficer (give title low) P and GM Ap		Other (s below) and Reager		
113 HARTWELL AVENUE						4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Ir	Individual or Joint/Group Filing (Check Applicable					
(Street) LEXINGTON MA 02421													- 1	X Form filed by One Reporting Person Form filed by More than One Reporting Person						
(City)	(5)	State)	(Zip)												Person					
		Tak	ole I - Noi	n-Deri	vativ	e Se	curities	s Acc	quired, [	Disp	osed o	f, or E	ene	eficiall	y Owned					
1. Title of Security (Instr. 3)  2. Transa Date (Month/L						2A. Deemed Execution Date, if any (Month/Day/Year)		Transaction Dispose Code (Instr. 5)		4. Securi Disposed 5)				Securitie Benefici Owned I	Securities For Beneficially (D)		n: Direct or Indirect nstr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)		
									Code	v	Amount	unt (A) or (D)		Price			Transac			
Common Stock 03/12					2/201	/2019		A		7,228	7,228 <sup>(1)</sup> A :		\$0.00	14,3	14,359 <sup>(2)</sup>		D			
			Table II -						uired, Di , options						Owned					
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	3A. Deeme Execution if any (Month/Day	Date,	4. Transa Code ( 8)		of		6. Date Exe Expiration (Month/Day	Date	of Securi (r) Underlyii		ying tive Security		8. Price of Derivative Security (Instr. 5)	9. Numbe derivative Securities Beneficial Owned Following Reported Transactio (Instr. 4)	e S Illy	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership t (Instr. 4)	
					Code	v	(A)		Date Exercisable		expiration pate	Title		mount or lumber of shares						
Stock Option (right to buy)	\$22.69	03/12/2019			A		24,640		(3)	0	3/12/2029	Commo Stock	n 2	24,640	\$0.00	24,64	0	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 March 12, 2020 and vest as to an additional 2.083% of the shares per month for 36 months thereafter.
- 2. Includes 14,359 restricted stock units.
- 3. The option vests as to 25% of the shares on March 12, 2020 and vests as to an additional 2.083% of the shares per month for 36 months thereafter.

## Remarks:

/s/ Brian Keane, Attorney-in-

03/14/2019

<u>fact</u>

\*\* 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.