FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

	9	

STATEMENT	OF CHANGES	S IN BENEFICIAL	OWNERSHIP

OMB APPROVAL										
OMB Number: 3235-02										
Estimated average burden										
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

Name and Address of Reporting Person*     Roskey Mark T.						2. Issuer Name <b>and</b> Ticker or Trading Symbol Quanterix Corp [ QTRX ]								(Check	all appli Directo	or		10% Ov	n(s) to Issuer  10% Owner  Other (specify	
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						3. Date of Earliest Transaction (Month/Day/Year) 06/15/2020								X	below)				·	
(Street) BILLER	ICA M	[A	01821		4. 1	4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Indiv Line) X	Form f	al or Joint/Group Filing (Ch orm filed by One Reporting orm filed by More than On erson		orting Perso	ng Person	
(City)	(S	tate)	(Zip)																	
		Tab	le I - I	Non-Deriv	<i>r</i> ative	e Sec	uriti	ies A	cquir	ed, D	isposed o	of, or B	enefic	ially	Owned	I				
1. Title of Security (Instr. 3)  2. Transaction Date (Month/Day/Ye					.	Execution Date,		3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 and			Benefic Owned		ies Fo cially (D) Following (I)		n: Direct or Indirect nstr. 4)	7. Nature of Indirect Beneficial Ownership		
									Code	v	Amount	(A) or (D)	Price			ed ction(s) 3 and 4)			(Instr. 4)	
Common Stock 06/15/202						0			M		5,000	A	\$2.9	96 26		,648 <sup>(1)</sup>		D		
Common Stock 06/15/202					020	20		S		5,000(2)	D	\$24.80	.8049 <sup>(3)</sup>		21,648 <sup>(1)</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	Date (Month/Day/Year) if	Execution Date, if any			ransaction code (Instr.		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		e Exer ation D h/Day/		7. Title and Amount of Securities Underlying Derivative Secur (Instr. 3 and 4)		De Se (Ir	Price of Perivative Security Instr. 5)	9. Number derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	у	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)	
					Code	v	(A)	(D)	Date Exerc	isable	Expiration Date	Title	Amou or Numb of Share	er						
Stock Option (Right to Buy)	\$2.96	06/15/2020			M			5,000	(4	4)	09/23/2024	Common Stock	5,00	0	\$0.00	94,455		D		

## **Explanation of Responses:**

- 1. Includes 16,753 restricted stock units.
- 2. The sale of shares reported on this Form 4 was made pursuant to a Rule 10b5-1 trading plan adopted on December 13, 2019.
- 3. This transaction was executed in multiple trades at prices ranging from \$24.57 to \$25.10. 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.
- 4. Option is fully vested.

## Remarks:

/s/ Brian Keane, Attorney-in-

06/15/2020

**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.