Che

FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

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

Check this box if no longer subject
to Section 16. Form 4 or Form 5
obligations may continue. See
Instruction 1(h)

## STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

**OMB APPROVAL** 3235-0287 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

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 Direc			.0% O	
_	(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE					3. Date of Earliest Transaction (Month/Day/Year) 07/02/2021								X	below			below)	
(Street) BILLER (City)	(Street) BILLERICA MA 01821				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 Person					
		Table	1 - 1	Non-Deriva	itive	Secui	rities A	cqu	uirec	d, Di	sposed o	of, or I	3enefic	ially	Own	ed			
1. Title of Security (Instr. 3)			2. Transaction Date (Month/Day/Ye	ear)	2A. Deen Executio if any (Month/D	n Date,		Transaction Code (Instr.		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 a			nd 5) Secu Bene Own		cially I Following	6. Ownership Form: Direct (D) or Indirect (I)		7. Nature of Indirect Beneficial Ownership	
								Cod	de V	<i>,</i>	Amount	(A) or (D)	Price		Reported Transaction(s) (Instr. 3 and 4)		(Instr. 4)		(Instr. 4)
Common Stock				07/02/2021				S			70(1)	D	\$57.63	96(2) 24		,066 <sup>(3)</sup>	D		
Common Stock				07/02/202	1			S			22(1)	D	\$58.38	381(4) 2		24,044 <sup>(3)</sup>			
Common Stock				07/02/202	1			S			54 <sup>(1)</sup>	D	\$59.34	32 <sup>(5)</sup>	23,990(3)		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)	Derivative Conversion Date Execution Date Courity or Exercise (Month/Day/Year) if any			cution Date, ny	4. Transaction Code (Instr. 8)		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		Expiration (Month/Da			Amo Secu Unde Deriv	,	Der Sec (Ins	rice of ivative urity tr. 5)	9. Number derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	Owr Forr Dire or Ir (I) (I	ership n: ct (D) direct nstr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)
					Cod	e V	(A) (I		Date Exerc	isable	Expiration Date	n Title	Amount or Number of Shares						

## **Explanation of Responses:**

- 1. The sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 472 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$57.13 to \$58.11. 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 13,293 restricted stock units
- 4. This transaction was executed in multiple trades at prices ranging from \$58.17 to \$59.03. 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
- 5. This transaction was executed in multiple trades at prices ranging from \$59.17 to \$59.51. 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.

## Remarks:

/s/ Brian Keane, Attorney-in-

07/06/2021

**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.