FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

D.C. 20549
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											
	OMB Number: 3235-028											
	Estimated average burden											
ı	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 <b>and</b> Ticker or Trading Symbol  Quanterix Corp [ QTRX ]								Check	all app	o of Reportir licable) tor er (give title	ng Pei	rson(s) to Is 10% O Other (	wner
(Last) C/O QU.		3. Date of Earliest Transaction (Month/Day/Year) 06/01/2021								X	below)  SVP, Strategic F		ic Pai	below)					
900 MIDDLESEX TURNPIKE  (Street)  BILLERICA MA 01821  (City) (State) (Zip)						4. If Amendment, Date of Original Filed (Month/Day/Year)								. Indiv ine) X	′				
		Table	I - N	Non-Deriva	tive	Secui	rities	Ac	quir	ed, Di	isposed c	f, or I	<b>Benefic</b>	ially	Own	ed			
1. Title of Security (Instr. 3)  2. Transaction Date (Month/Day/Yea					ear) if	2A. Deemed Execution Date, if any (Month/Day/Year)			3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 au			nd 5) Secu Bend		rities eficially ed Following		n: Direct or rect (I)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
		(	Code	v	Amount			(A) or (D)	Price		Transa	r. 3 and 4)		.,	()				
Common Stock 06/01/2021						21			S		153 <sup>(1)</sup>	D	\$48.3181(2)		24,136 <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)	rivative conversion or Exercise str. 3) str. 3) Price of Derivative Security  Date (Month/Day/Year) (Month/Day/Year)  Execution Date, if any (Month/Day/Year)			Code 8)	Transaction of Code (Instr. Derivativ			Expiration Date (Month/Day/Year)			Amo Secu Unde Deriv Secu 3 and	Amount or Number of	ınt per				10. 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 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 \$48.00 to \$48.75. 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 13,765 restricted stock units.

## Remarks:

/s/ Brian Keane, Attorney-in-

<u>fact</u>

\*\* Signature of Reporting Person Date

06/02/2021

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.