SEC Form 4

FORM 4

Check this box if no longer subject

UNITED STATES SECURITIES AND EXCHANGE COMMISSION Washington, D.C. 20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL

OMB Number:	3235-0287								
Estimated average burden									
hours per response:	0.5								

obligations may continue. See Instruction 1(b).	Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934	hours per response: 0.5
1. Name and Address of Reporting Person* <u>Roskey Mark T.</u>	or Section 30(h) of the Investment Company Act of 1940 2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]	5. Relationship of Reporting Person(s) to Issuer (Check all applicable)
(Last) (First) (Middle) C/O QUANTERIX CORPORATION,	3. Date of Earliest Transaction (Month/Day/Year) 04/01/2021	Director 10% Owner X Officer (give title below) Other (specify below) SVP, Strategic Partnerships
900 MIDDLESEX TURNPIKE	4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Individual or Joint/Group Filing (Check Applicable
(Street) BILLERICA MA 01821		Line) X Form filed by One Reporting Person Form filed by More than One Reporting Person
(City) (State) (Zip)		

Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned

1. Title of Security (Instr. 3)	2. Transaction Date (Month/Day/Year)	2A. Deemed Execution Date, if any (Month/Day/Year)	3. Transa Code (8)		4. Securities Disposed Of			5. Amount of Securities Beneficially Owned Following	6. Ownership Form: Direct (D) or Indirect (I)	7. Nature of Indirect Beneficial Ownership
			Code	v	Amount	(A) or (D)	Price	Reported Transaction(s) (Instr. 3 and 4)	(Instr. 4)	(Instr. 4)
Common Stock	04/01/2021		S		41 ⁽¹⁾	D	\$61.9196 ⁽²⁾	24,531 ⁽³⁾	D	
Common Stock	04/01/2021		S		85(1)	D	\$ 62.6984 ⁽⁴⁾	24,446 ⁽³⁾	D	
Common Stock	04/01/2021		S		10(1)	D	\$63.36	24,436 ⁽³⁾	D	

Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities)

L	(e.g., puis, cans, warrants, options, conventible securities)															
	1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	3A. Deemed Execution Date, if any (Month/Day/Year)	4. Transa Code (8)		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)	11. Nature of Indirect Beneficial Ownership (Instr. 4)
					Code	v	(A)	(D)	Date Exercisable	Expiration Date	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 \$61.30 to \$62.27. 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 14,709 restricted stock units.

4. This transaction was executed in multiple trades at prices ranging from \$62.30 to \$63.21. 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-

fact

04/05/2021

** 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.