## FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPRO	DVAL
OMB Number:	3235-0287
Estimated average burd	en
hours per response:	0.5
	OMB Number: Estimated average burd

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 ]									all app	olicable)		Owner (specify
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						3. Date of Earliest Transaction (Month/Day/Year) 11/01/2019								Λ	SVP,	ow) below) P, Commercial & Accelerator		
(Street) BILLERICA MA 01821  (City) (State) (Zip)						4. If Amendment, Date of Original Filed (Month/Day/Year)								Individual or Joint/Group Filing (Check Applicable Line)     X Form filed by One Reporting Person     Form filed by More than One Reporting Person				
Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																		
1. Title of Security (Instr. 3)  2. Transaction Date (Month/Day/Ye						Execution Date,		,	3. 4. Securities Acquire Disposed Of (D) (Instance)						Securities Beneficially		6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
						v	Amount	(A) or (D)	Price		Trans	action(s) 3 and 4)		()				
Common	19			S		46(1)	D	D \$20.859		4 <sup>(2)</sup> 14,397 <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)	2. Conversion or Exercise Price of Derivative Security	onversion Exercise (ice of erivative ecurity    Continue   Continu			saction le (Instr. Scurities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)			Date Exercisable and Expiration Date (Month/Day/Year)  Date Exercisable Expiration Date			7. Title and Amount of Securities Underlying Derivative Security (Instr. and 4)  Amount or Number of Shares				9. Number of derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	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 was made solely to cover tax obligations upon the vesting of 148 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$20.79 to \$20.92. 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 11,689 restricted stock units.

## Remarks:

/s/ Brian Keane, Attorney-in-11/04/2019 fact

\*\* Signature of Reporting Person

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.