FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington,	D.C.	20549	

wasnington,	D.C.	20548

Estimated average burden	OVAL	
STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP	OMB Number:	3235-0287
	OMB Number: Estimated average bure	rden
	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(b) of the Investment Company Act of 1940

1. Name and Address of Reporting Person* Roskey Mark T. (Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE (Street)						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX] 3. Date of Earliest Transaction (Month/Day/Year) 05/03/2021 4. If Amendment, Date of Original Filed (Month/Day/Year)						r) 6	S. Relationship of Reporting Person(s) to Issuer (Check all applicable) Director 10% Owner X Officer (give title below) SVP, Strategic Partnerships 6. Individual or Joint/Group Filing (Check Applic Line)				wner specify s		
(City))182 Zip)	1									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. Transac Date			2. Transaction Date (Month/Day/Ye	Execution ear) if any		2A. Deemed Execution Date, if any (Month/Day/Year)		3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or			5. A Sec Ber Ow		ount of ties cially d Following	6. Ownership Form: Direct (D) or Indirect (I)		7. Nature of Indirect Beneficial Ownership	
								С		v	Amount	(A) or (D)	Price			orted (Instr. 4) saction(s) tr. 3 and 4)		4)	(Instr. 4)
Common	Common Stock 05/03/202							\top	S	$\dagger \dagger$	75 ⁽¹⁾	D	\$57.24	71(2)	24	24,361 ⁽³⁾		D	
Common	Stock			05/03/202	(2021 S 58 ⁽¹⁾ D \$58.1591 ⁽⁴⁾ 24,303 ⁽³⁾]	D						
Common	Stock			05/03/202	1				S		9(1)	D	\$59.01	29 ⁽⁵⁾	24	,294 ⁽³⁾		D	
Common	Stock			05/03/202	1				S		5 ⁽¹⁾	D	\$61.	36	6 24,289 ⁽³⁾ D				
	Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities)																		
Derivative Conversion Date		3. Transaction Date (Month/Day/Year)	Exe if ar	Deemed cution Date, ny inth/Day/Year)		action (Instr.	5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		Expiration D (Month/Day/		ay/Year) Securities Underlying Derivative Security (In 3 and 4) Amo or Num of		unt of unities erlying vative unity (Instr. d 4) Amount or Number of	Deri Sec (Ins	rice of ivative urity tr. 5)	9. Number derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	y Di Oi (I)	D. wnership orm: irect (D) r Indirect ((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 \$56.67 to \$57.63. 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 14,237 restricted stock units
- 4. This transaction was executed in multiple trades at prices ranging from \$58.75 to \$59.45. 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 \$57.67 to \$58.64. 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-

05/05/2021

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.