FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

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

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

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 and Ticker or Trading Symbol Quanterix Corp [QTRX]										all app Direc	ionship of Reportir all applicable) Director Officer (give title		erson(s) to Is 10% Ov Other (s	wner
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						3. Date of Earliest Transaction (Month/Day/Year) 03/01/2022										below) SVP, Strategio		c Pa	below)	· /
(Street) BILLERICA MA 01821 (City) (State) (Zip)					4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)									i. Indiv ine) X	I				
		Table	I - N	Non-Deriva	tive	Secui	rities	Ac	qui	red, D)is	posed o	f, or I	3enefic	ially	Own	ed			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Yea					ear) it	2A. Deemed Execution D if any (Month/Day		, 	3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 an			d (A) or r. 3, 4 and	5)	Securi Benefi Owned			m: Direct or irect (I)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
						Code V		An	nount	(A) or (D)	Price		Reported (Ir Transaction(s) (Instr. 3 and 4)		(1115	4)	(111511.4)			
Common	03/01/202	22				S			122(1)	D	\$31.20	66 ⁽²⁾	22,540 ⁽³⁾			D				
Common Stock 03/01/20					2				S			44(1)	D	\$32.15		22,496 ⁽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, Gecurity or Exercise (Month/Day/Year) if any					5. Numb of Operivati Securiti Acquire (A) or Dispose of (D) (Instr. 3, and 5)		vative rities iired r osed) r. 3, 4	Expiration Dat (Month/Day/Ye			n Date Amount of Securities Underlying Derivative Security (Inst 3 and 4)		unt of rities erlying rative rity (Instr.	Deri Sec (Ins	rice of ivative urity tr. 5)	tive derivative ity Securities		10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)
					Code	Code V (A) (E		(D)		Date Exercisab		Expiration Date	Title	Number of					L	

Explanation of Responses:

- 1. The sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 542 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$30.84 to \$31.83. 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 8,604 restricted stock units.

Remarks:

/s/ Brian Keane, Attorney-in-

fact

03/03/2022

** 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.