FORM 4

Check this box if no longer subject to Section 16. Form 4 or Form 5 obligations may continue. See

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

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

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

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

Instruc	tion 1(b).			Filed							ities Exchar ompany Act						,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,,		
Name and Address of Reporting Person* Roskey Mark T.				2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]									all app	nship of Reportin applicable) Director Officer (give title		rson(s) to Is 10% Ov Other (s	ner		
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE				3. Date of Earliest Transaction (Month/Day/Year) 01/03/2022									Λ	below SV	below) P, Strategic Partnerships				
(Street) BILLER (City)	ICA M	A 0	1821 'ip)	1	4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)								i. Indiv ine) X	vidual or Joint/Group Filing (Check Applicable Form filed by One Reporting Person Form filed by More than One Reporting Person				
		Table	I - N	Non-Deriva	tive	Secui	rities	Acc	quire	ed, Di	sposed c	f, or I	Benefic	ially	Own	ed			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Yes				Execution ear) if any		on Date, Day/Year)		3. Transaction Code (Instr. 8)					ind 5) S B		5. Amount of Securities Beneficially Owned Following Reported		n: Direct or rect (I)	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
								ode	V A	mount	(A) or (D)	Price		Transaction(s) (Instr. 3 and 4)		(Instr. 4)		(111501.4)	
Common Stock 01/03/2023					22			S		149(1)	D	\$41.90	1.9066(2)		23,120(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)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	Exe if an	Deemed cution Date, ny nth/Day/Year)	4. Transaction Code (Instr. 8) 5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)				Expiration Date (Month/Day/Year) Securi Underl Deriva Securi 3 and 4			unt of crities crlying vative crity (Instr. d 4)	Deri Sec (Ins	s. Price of Derivative Security Instr. 5)	9. Number derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	Owner Form: Direct or Indi (I) (Ins	Ownership	11. Nature of Indirect Beneficial Ownership (Instr. 4)	
					Code V (A) (D)					e rcisable	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 \$41.71 to \$42.16. 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 10,461 restricted stock units.

Remarks:

/s/ Brian Keane, Attorney-in-

fact

** Signature of Reporting Person Date

01/05/2022

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.