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:	0.5							

Instruc	tion 1(b).			Filed								es Exchan pany Act					liouis	perit		0.5
Name and Address of Reporting Person* Roskey Mark T.				2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]									5. Relationship of Reporting Person(s) to Issuer (Check all applicable) Director 10% Owner X Officer (give title Other (specific					wner		
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE					3. Date of Earliest Transaction (Month/Day/Year) 12/01/2021									SVP, Strategic Partnerships						
(Street) BILLER (City)			182 Zip)	1	4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)								Indivi ne) X	iividual or Joint/Group Filing (Check Applicable Form filed by One Reporting Person Form filed by More than One Reporting Person					
		Table	I - I	Non-Deriva	ative	Secu	rities	s Ac	cquir	ed, D	isp	osed o	f, or I	3enefici	ally	Own	ed			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Ye				Execution Da		on Date, Tr			Transaction Disposed Of (Code (Instr.				nd 5) Sec Ben Owi		Amount of curities neficially ned Following ported		m: Direct or rect (I)	7. Nature of Indirect Beneficial Ownership (Instr. 4)		
								Code	v	Amo	ount	(A) or (D)	Price	Transaci (Instr. 3		ction(s)	L.	,	, ,	
Common	Common Stock 12/01/202				1			S		142(1)		D	\$40.2643(2)		23,274 ⁽³⁾			D		
Common Stock 12/01/202					1			S			5(1)	D	\$41		23,269			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)	rivative Conversion Date Execution Date, curity or Exercise (Month/Day/Year) if any			action (Instr.			Expiration Date (Month/Day/Year)			7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)				9. Number derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	,	10. 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 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 \$40.00 to \$40.92. 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 10,933 restricted stock units.

Remarks:

/s/ Brian Keane, Attorney-in-

fact

12/02/2021

** 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.