FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

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

theck this box if no longer subject
Section 16. Form 4 or Form 5
bligations may continue. See
etruction 1(h)

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL							
OMB Number: 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		ng Pers	son(s) to Is 10% O Other (wner
(Last) (First) (Middle) C/O QUANTERIX CORPORATION,						3. Date of Earliest Transaction (Month/Day/Year) 12/01/2020									SVP,			below) Accelerator	
900 MIDDLESEX TURNPIKE (Street) BILLERICA MA 01821					4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Indiv Line)	-7				
(City)	(St		Zip)	Jan Barina	<u> </u>	tive Securities Acquired, Disposed of, or Beneficially Owned													
		Table	I - I	Non-Deriva	itive	Secu	rities A	4CC	quire	ea, D	isposed (ot, or	Benet	icially	Own	ed			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Yell)				ear) i	2A. Deemed Execution I if any (Month/Day		Ti C	3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 an			d 5) Secur Benef		icially d Following	Form (D) o	n: Direct r ect (I)	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
								С	ode	v	Amount	(A) or (D)	Price		Transa	saction(s) r. 3 and 4)		4)	(111341. 4)
Common Stock				12/01/202	20				S		42(1)	D	\$43.3	8058 ⁽²⁾	21	21,984 ⁽³⁾		D	
Common Stock 12/01/2				12/01/202	20				s 49 ⁽¹⁾		D	\$44.1	.341(4)	21	1,935 ⁽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)	perivative Conversion Date Execution Date, if any					ransaction of code (Instr. Deriv		Expiration (Month/Date ed				Amo Seci Und Deri Seci	7. Title and Amount of Securities Underlying Derivative Security (Inst 3 and 4)		rice of ivative urity itr. 5)	rative derivative rity Securities		10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)
					Code V (A) (D)				Date Exe	ate Expiration tercisable Date		n Title	Amou or Numb of Share	er					

Explanation of Responses:

- 1. The sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 298 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$43.00 to \$43.94. 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 14,965 restricted stock units. Also includes 838 shares acquired under the Quanterix Corporation Employee Stock Purchase Plan.
- 4. This transaction was executed in multiple trades at prices ranging from \$44.01 to \$44.39. 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-

fact

12/03/2020

** 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.