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, 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							urities Exc Company			of 1934			lilouis	perio		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 Issu (Check all applicable) Director Officer (give title Other (spe					wner		
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE				3. Date of Earliest Transaction (Month/Day/Year) 11/02/2021										below) below) SVP, Strategic Partnerships						
(Street) BILLER (City)			1821 Zip)		4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)								Indivi ne) X	lividual or Joint/Group Filing (Check Applicable Form filed by One Reporting Person Form filed by More than One Reporting Person					
		Table	I - N	on-Deriva	tive	Secui	rities	Ac	quir	ed, D	ispose	d of	f, or E	Benefici	ally	Own	ed			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Ye			Execution Date,		٠,	Code (Instr.					i (A) or :. 3, 4 and 5	nd 5) Se Be Ov		i. Amount of Securities Beneficially Dwned Following Reported		m: Direct or irect (I)	7. Nature of Indirect Beneficial Ownership (Instr. 4)			
									Code	v	Amount		(A) or (D)	Price	Transaction(s) (Instr. 3 and 4)		ction(s)	, , ,		(
Common	Stock			11/02/202	1				s 100 ⁽¹⁾ D \$52.3974 ⁽²⁾ 23,455 ⁽³⁾ D			D								
Common	ommon Stock 11/02/202			1	s 39 ⁽¹⁾ D \$53.		\$53.144	23,416 ⁽³⁾		,416 ⁽³⁾		D								
		Tal	ble II	- Derivati (e.g., pu												wne	t			
1. Title of Derivative Security (Instr. 3)	ative Conversion Date Execution Date, if any		Code 8)	Transaction of Code (Instr. Derivative		Exp (Mo	Expiration Date (Month/Day/Year) Date Expiration			7. Title and Amount of Securities Underlying Derivative Security (Inst 3 and 4) Amour or Numbe of Title Shares		nt er		9. Number of 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 \$51.86 to \$52.85. 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.

(D)

- 4. This transaction was executed in multiple trades at prices ranging from \$52.96 to \$53.78. 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

11/04/2021

** 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.