FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

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

STATEMENT	OF CHANGES	IN BENEFICIAL	OWNERSHIP

OMB APPROVAL 3235-0287 Estimated average burden hours per response: 0.5

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

Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934 or Section 30(h) of the Investment Company Act of 1940

(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE Street BILLERICA MA 01821	1. Name and Address of Reporting Person* Roskey Mark T.						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]								Check	all app	o of Reportir licable) tor er (give title	ng Per	rson(s) to Is 10% O Other (vner
City (Street) BILLERICA MA O1821 City (State) (Zip) City (Month/DaylYear) (Month	C/O QUANTERIX CORPORATION,															belov	v) ``	below)		
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Year) (Month/Day/Year) 2. A. Deemed Execution Date, if any (Month/Day/Year) (Month/Day/Year) (Month/Day/Year) 2. Transaction Date (Month/Day/Year) 3. Transaction Date (Month/Day/Year) 3. Transaction Date (Month/Day/Year) 3. Transaction Date (Month/Day/Year) 3. Transaction Date (Month/Day/Year) 4. Transaction Date (Month/Day/Year) 3. Transaction Date (Month/Day/Year) 3. Transaction Date (Month/Day/Year) 4. Transaction Date (Month/Day/Year) 3. Transaction Date (Month/Day/Year) 4. Transaction Date (Month/Day/Year) 5. Amount of Securities Peneficially Owned (Instr. 4) 5. Amount (A) or Price (D) or Indirect (I) (Instr. 4) 5. Amount of Securities Peneficially Owned (Instr. 8) 7. Tritle and Amount of Securities Peneficially Owned (Month/Day/Year) 8. Price of Derivative Securities Peneficially Owned (Month/Day/Year) 8. Price of Derivative Securities Peneficially Owned (Month/Day/Year) 9. Number of Derivative Securities Peneficially Owned (Month/Day/Year) 10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4) 11. Nature Price of Derivative Securities Peneficially Owned (Month/Day/Year) 12. Nature Price of Derivative Securities Peneficially Owned (Month/Day/Year) 13. Nature Price of Derivative Securities Peneficially Owned (Instr. 4) 14. Securities Acquired (A) or Price or Disposed of (D) (Instr. 3, 4 and 5) 15. Amount of Securities Peneficially Owned (Instr. 4) 16. Ownership Form: Direct (D) Owned (Instr. 4) 17. Title and Amount of Securities Peneficially Owned (Instr. 4) 18. Price of Derivative Securities Peneficially Owned (Instr. 4) 19. Number of Derivative Securities	(Street) BILLERICA MA 01821					4. If Amendment, Date of Original Filed (Month/Day/Year)								ine)	X Form filed by One Reporting Person Form filed by More than One Reporting					
Date (Month/Day/Year) Execution Date, if any (Month/Day/Year) Execution Date (Month/Day/Year) Execut			Table	I - I	Non-Deriva	tive	Secui	rities	Ac	quire	ed, Di	isposed o	f, or I	3enefic	ially	Own	ed			
Common Stock 11/02/2020 S 91(1) D \$36.9687(2) 21,188(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 (Month/Day/Year) (Month/Day/Year) 3A. Deemed Execution Date (Month/Day/Year) 4. Transaction Code (Instr. 8) 5. Number of Exercisable and Exercisable and Execution Date (Month/Day/Year) 6. Date Exercisable and Execution Date (Month/Day/Year) 7. Title and Amount of Securities (Month/Day/Year) 8. Price of Derivative Securities (Month/Day/Year) 9. Number of derivative Securities (Month/Day/Year) 11. Nad Security (Instr. 5) 12. Ownership Form: Derivative Securities (Month/Day/Year) 13. Transaction Date (Month/Day/Year) 14. Nad Security (Instr. 5) 15. Number of Ownership Securities (Month/Day/Year) 16. Date Exercisable and Expiration Date (Month/Day/Year) 17. Title and Amount of Securities (Month/Day/Year) 18. Price of Derivative Securities (Month/Day/Year) 19. Number of derivative Securities (Month/Day/Year) 19. Number of derivative Securities (Month/Day/Year) 10. Ownership Form: Direct (D) Owned Following Reported Transaction(s) (Instr. 4) 11. Nad Transaction (Securities (Month/Day/Year) 12. Nad Transaction (Securities (Month/Day/Year) 13. Amount of Date (Month/Day/Year) 14. Amount of Derivative Securities (Month/Day/Year) 15. Number of Owner (Month/Day/Year) 16. Date Exercisable and Ex	Date			Execution Date, ar) if any			Transaction Code (Instr.					nd 5) Secur Benef Owne		rities Fe ficially (E ed Following In		n: Direct or rect (I)	7. Nature of Indirect Beneficial Ownership			
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) Price of Derivative Security Secu								-	Code	v .	Amount	(A) or (D)	Price		Transa	action(s)	(11130	4)	(1130.4)	
1. Title of Derivative Security (Instr. 3) In the of Derivative Security (Instr. 4) In the of Derivative Security (Instr. 5) In the of Derivative Security (Instr. 4) In the of Derivative Security (Instr. 4) In the of Derivative Security (Instr. 5) In the of Derivative Security (Instr. 4) In the of Derivative Security (Instr. 5) In the of Derivative Security (Instr. 4) In the of Derivative Security (Instr. 4) In the of Derivative Security (Instr. 4) In the of Derivative Security (Instr. 5) In the of Derivative Security (Instr. 4) In the of Derivative Securit	Common Stock 11/02/2020			.0			S		91(1)	D	D \$36.9687 ⁽²⁾		21,188 ⁽³⁾			D				
Derivative Security (Instr. 3) Price of Derivative Security Security Date (Month/Day/Year) Execution Date, if any (Month/Day/Year) Execution Date, if any (Month/Day/Year) Execution Date, if any (Month/Day/Year) Expiration Date (Month/Day/Year) Expi																				
	Derivative Security	Conversion or Exercise Price of Derivative	Date	Exe if ar	cution Date, ny	Transa Code	nsaction de (Instr. Securitie Acquired (A) or Disposed of (D) (Instr. 3,		ative ities red sed 3, 4	Expiration (Month/Da		Date //Year)	Amo Secu Unde Deriv Secu 3 and	unt of rities erlying vative rity (Instr. i 4) Amount or Number	Derivative Security (Instr. 5)		derivative Securities Beneficially Owned Following Reported Transaction	y	Ownership Form: Direct (D) or Indirect	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 298 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$36.52 to \$37.33. 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 15,263 restricted stock units.

Remarks:

/s/ Brian Keane, Attorney-in-

fact

** Signature of Reporting Person Date

11/03/2020

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.