SEC Form 4	
------------	--

FORM 4

Check this box if no longer subject

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							

to Section 16. Fe obligations may 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			Estimated aver hours per resp	0	0.5
HRUSOVSK (Last)	ss of Reporting Perso <u>Y E KEVIN</u> (First) RIX CORPORATI	(Middle)	2. Issuer Name and Ticker or Trading Symbol <u>Quanterix Corp</u> [QTRX] 3. Date of Earliest Transaction (Month/Day/Year) 07/13/2021		all applicab Director Officer (giv below)	,	10% Owner Other (speci below)	
900 MIDDLESI	EX TURNPIKE							
(Street) BILLERICA	МА	01821	4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Indivi Line) X	Form filed	loint/Group Filing (Check Applicabl led by One Reporting Person led by More than One Reporting		
(City)	(State)	(Zip)		Person			ł	

Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned

1. Title of Security (Instr. 3)	2. Transaction Date (Month/Day/Year)	2A. Deemed Execution Date, if any (Month/Day/Year)	3. Transa Code (8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 and 5)		5. Amount of Securities Beneficially Owned Following Reported	Form: Direct (D) or Indirect (I)	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
	Code		v	Amount	(A) or (D)	Price	Transaction(s) (Instr. 3 and 4)			
Common Stock	07/13/2021		S		3,249 ⁽¹⁾	D	\$57.221 ⁽²⁾	944,018 ⁽³⁾	D	
Common Stock	07/13/2021		S		1,751 ⁽¹⁾	D	\$ 57.8834 ⁽⁴⁾	942,267 ⁽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)	3A. Deemed Execution Date, if any (Month/Day/Year)	4. Transa Code (8)		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)			of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4		of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4		Expiration Date		and 7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		nount of curities derlying rivative curity (Instr. 5)		9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s) (Instr. 4)	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)
				Code	v	(A)	(D)	Date Exercisable	Expiration Date	Title	Amount or Number of Shares										

Explanation of Responses:

1. The sale of shares was effected pursuant to a 10b5-1 trading plan adopted on December 14, 2020.

2. This transaction was executed in multiple trades at prices ranging from \$56.58 to \$57.55. 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 97,244 restricted stock units and 39,815 shares of common stock that are subject to performance based vesting.

4. This transaction was executed in multiple trades at prices ranging from \$57.70 to \$58.24. 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:



in-Fact

07/14/2021

** 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.