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

## 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							

Instruction 1(b).		Filed pursuant to Section 16(a) of the Securities Exchange Act of 19	hours per	hours per response: 0.5			
		or Section 30(h) of the Investment Company Act of 1940					
1. Name and Address of Rep HRUSOVSKY E K	0	2. Issuer Name <b>and</b> Ticker or Trading Symbol <u>Quanterix Corp</u> [ QTRX ]	(Check all appl	icable)	,		
<u></u>			X Direct	or	10% Owner		
(Last) (First)	(Middle)	3. Date of Earliest Transaction (Month/Day/Year)	X Office below	r (give title )	Other (specify below)		
C/O QUANTERIX CO	RPORATION,	10/22/2021		Chairman an	airman and CEO		
900 MIDDLESEX TUP	NPIKE						
(Street)		4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Individual or Line)	Joint/Group Fil	ing (Check Applicable		
BILLERICA MA	01821		X Form	filed by One Re	porting Person		
	01021		Form Perso		an One Reporting		
(City) (State)	(Zip)						

#### 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	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership	
		Code		v	Amount	(A) or (D)	Price	Transaction(s) (Instr. 3 and 4)	(1150.4)	(Instr. 4)
Common Stock	10/22/2021		S		1,900 <sup>(1)</sup>	D	<b>\$50.1891</b> <sup>(2)</sup>	909,416 <sup>(3)</sup>	D	
Common Stock	10/22/2021		S		3,000 <sup>(1)</sup>	D	<b>\$</b> 50.9992 <sup>(4)</sup>	906,416 <sup>(3)</sup>	D	
Common Stock	10/22/2021		S		100(1)	D	\$51.63	906,316 <sup>(3)</sup>	D	

## Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned

	(e.g., puts, calls, warrants, options, convertible securities)														
1. Title of Derivativ Security (Instr. 3)		3. Transaction Date (Month/Day/Year)	3A. Deemed Execution Date, if any (Month/Day/Year)	4. Transa Code ( 8)		of		6. Date Exerc Expiration Da (Month/Day/Y	7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		8. Price of Derivative Security (Instr. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s) (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 \$49.56 to \$50.54. 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 73,439 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 \$50.63 to \$51.50. 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:** 

# <u>/s/ Brian Keane, as Attorney-</u>

in-Fact

<u>10/22/202</u>1

\*\* Signature of Reporting Person Date

Reminder: Report on a separate line for each class of securities beneficially owned directly or indirectly.

 $^{\star}$  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.