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

## FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

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

# STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

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

OMB APPROVAL	
--------------	--

OMB Number:	3235-0287
Estimated average b	ourden
hours per response:	0.5

1. Name and Address of Reporting Person*			2. Issuer Name <b>and</b> Ticker or Trading Symbol <u>Quanterix Corp</u> [ QTRX ]	5. Relationship of Reporting Person(s) to Issuer (Check all applicable)					
<u>HRUSOVSKY E KEVIN</u>			Canar 1	X	Director	10% Owner			
					Officer (give title	Other (specify			
(Last)	(First)	(Middle)	3. Date of Earliest Transaction (Month/Day/Year)	X	below)	below)			
C/O QUANTERIX CORPORATION,			01/05/2021		Chairman, President	: & CEO			
900 MIDDLESEX TURNPIKE									
			4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Individual or Joint/Group Filing (Check Applicat					
(Street)				Line)					
BILLERICA	МА	01821		X	Form filed by One Repor	ting Person			
					Form filed by More than	One Reporting			
					Person				
(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. Transaction Code (Instr. 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)	(1130.4)	(Instr. 4)
Common Stock	01/05/2021		S		3,510 <sup>(1)</sup>	D	<b>\$44.1317</b> <sup>(2)</sup>	1,015,788 <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 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)		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)	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 sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 7,123 RSUs.

2. This transaction was executed in multiple trades at prices ranging from \$43.91 to \$44.62. 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 135,394 restricted stock units and 39,815 shares of common stock that are subject to performance based vesting.

#### Remarks:

#### /s/ Brian Keane, as Attorney-01/06/2021

<u>in-Fact</u>

\*\* 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.