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

			()					
	ss of Reporting Perso	on <sup>*</sup>	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>				X	Director	10% Owner		
				x	Officer (give title	Other (specify below)		
(Last)	(First)	(Middle)	3. Date of Earliest Transaction (Month/Day/Year)		below)	Delow)		
C/O QUANTERIX CORPORATION,			04/01/2022	Chairman and CEO				
900 MIDDLESEX TURNPIKE								
			4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Indiv	idual or Joint/Group Filing	(Check Applicable		
(Street)				Line)				
BILLERICA	МА	01821		X	Form filed by One Repor	rting Person		
DILLENICA	MA	01021			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)	Code (Instr.		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 and 5)			Securities Beneficially	6. Ownership 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)	(Instr. 4)	(11150. 4)
Common Stock	04/01/2022		S		868(1)	D	\$29.7929 <sup>(2)</sup>	886,976 <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)		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		of Expiration Date (Month/Day/Year) Securities Acquired (A) or Disposed of (D) (Instr. 3, 4			nd 7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		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 2,068 RSUs.

2. This transaction was executed in multiple trades at prices ranging from \$29.25 to \$30.00. 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 39,214 restricted stock units and 39,815 shares of common stock that are subject to performance based vesting.

#### **Remarks:**

#### /s/ Brian Keane, as Attorney-04/05/2022

in-Fact

\*\* Signature of Reporting Person Date

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

 $^{\ast}$  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.