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

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

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

1. Name and Address of Reporting Person*  HRUSOVSKY E KEVIN						2. Issuer Name <b>and</b> Ticker or Trading Symbol Quanterix Corp [ QTRX ]										ionship of Reportin all applicable) Director		g Person(s) to Iss 10% Owi		
						3. Date of Earliest Transaction (Month/Day/Year) 03/16/2021									X	below	Officer (give title below)  Chairman, Pres		Other (s below) at & CEC	
(Street) BILLERICA MA 01821 (City) (State) (Zip)						4. If Amendment, Date of Original Filed (Month/Day/Year)									ine) X	vidual or Joint/Group Filing (Check Applicable  Form filed by One Reporting Person  Form filed by More than One Reporting  Person				
Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																				
Date				2. Transaction Date (Month/Day/Ye	Execution Year) if any		med on Date, Day/Year)		3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 an				nd 5) Secu Bene		rities F eficially ( ed Following II		ı: Direct r ect (I)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
								Code	v	Amount		(A) or (D)	Price		Transaction(s) (Instr. 3 and 4)		(111341	(Instr. 4)	(1130.4)	
Common		03/16/202	21				S <sup>(1)</sup>			900	D	\$60.42	23(2) 95		993,100(3)		D			
Common Stock				03/16/202	21				S <sup>(1)</sup>		4	4,100	D	\$61.51	49 <sup>(4)</sup>	989,000(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)	Execu if any	eemed ution Date, , th/Day/Year)		action (Instr.	5. Nu of Deriv Secu Acqu (A) or Dispo of (D) (Instr and 5	vative irities iired r osed ) r. 3, 4	Expiration Date (Month/Day/Year)			te ear)	Amo Secu Unde Deriv Secu 3 and	Amount or Number of	nt er		9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s) (Instr. 4)		LO. Ownership Form: Direct (D) Or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)

## **Explanation of Responses:**

- $1. \ The \ sale \ of \ shares \ was \ effected \ pursuant \ to \ a \ Rule \ 10b5-1 \ trading \ plan \ adopted \ on \ December \ 14, \ 2020.$
- 2. This transaction was executed in multiple trades at prices ranging from \$60.08 to \$60.74. 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 128,984 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 \$61.20 to \$61.72. 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:

/s/ Brian Keane, as Attorney-

03/16/2021

in-Fact

\*\* 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.