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 and Ticker or Trading Symbol Quanterix Corp [QTRX]									all app	onship of Reporting Pe Il applicable) Director		son(s) to Is	
(Last) (First) C/O QUANTERIX CORD 900 MIDDLESEX TURN		dle)	3. Date of Earliest Transaction (Month/Day/Year) 07/01/2020								X	below Cha	airman, Presid				
(Street) BILLERICA MA (City) (State)	0182 (Zip)		4. If Amendment, D			, Date of Original Filed (Month/Day/Year)						ine)	Form	I or Joint/Group Filing (Check Applicable rm filed by One Reporting Person rm filed by More than One Reporting rrson			
	Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																
1. Title of Security (Instr. 3)		2. Transaction Date (Month/Day/Yea	2A. Deem Execution ar) if any (Month/Da		on Date,		3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 an			d 5) Secur Benef		ities For icially (D) d Following Ind		: Direct	7. Nature of Indirect Beneficial Ownership (Instr. 4)
							ode	v	Amount (A		Price	Transa		action(s) 3 and 4)		130.4)	(11150.4)
Common Stock 07/01/2020)				S		3,060(1)	D	\$26.828	85 ⁽²⁾ 1,07		71,577 ⁽³⁾		D		
Common Stock 07/01/202		07/01/2020)				S		237(1)	D	\$27.45	7.45 ⁽⁴⁾		1,071,340(3)		D	
Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities)																	
Derivative Conversion Date	e Exemple Exem		4. Transaction Code (Instr. 8) 5. Nur Code (Instr. 8) 5. Nur Acqui (A) or Dispo of (D) (Instr. and 5) 5. Nur Code V (A)			ative rities ired osed	Expiration I (Month/Day et al., 4		y/Year)	Amo Secu Unde Deri Secu 3 an	Amount or Number of	unt per				.0. Ownership Form: Direct (D) Or Indirect I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)

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 \$26.39 to \$27.22. 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 178,132 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 \$27.39 to \$27.57. 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-

07/06/2020

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.