FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

	OMB APPROVAL									
	OMB Number:	3235-028								
-	Estimated average I	hurden								

0.5

hours per response:

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

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*						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]								5. Relationship of Reporting Person(s) to Issuer (Check all applicable)					
HRUSOVSKY E KEVIN					-									X	Direc	ctor	10%	Owner	
(Last) (First) (Middle)						Date of Earliest Transaction (Month/Day/Year)								X	Offic belov	er (give title w)	Othe belov	r (specify v)	
C/O QUANTERIX CORPORATION,					02/0	02/06/2020									Chairman, President & CEO				
900 MIDDLESEX TURNPIKE																			
(Ctract)						4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Individual or Joint/Group Filing (Check Applicable Line)					
(Street) BILLERICA MA 01821													X	X Form filed by One Reporting Person					
															Form filed by More than One Reporting Person				
(City)	(SI	ate) (Zip)																
		Tabl	e I - No	on-Deriv	ative	Sec	uritie	s Ac	quire	d, Dis	sposed o	f, or E	3enef	icially	Owne	ed			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day)						Execution Date,					es Acquired (A) o Of (D) (Instr. 3, 4 a		and 5) Secu Bene		icially d Following	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)		
									Code	v	Amount	(A) or (D) Pric		се	Transa	action(s) 3 and 4)		(iiisti. 4)	
Common Stock 02/06/20						020		S ⁽¹⁾		6,500	D	\$2	.7.56 ⁽²⁾	1,122,433 ⁽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)	3A. Dee Execution if any (Month/	on Date,	4. Transaction Code (Instr. 8)		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		6. Date Exerc Expiration Da (Month/Day/Y		te	7. Title and Amount of Securities Underlying Derivative Security (Inst and 4)		Der Sed (Ins	rice of ivative urity tr. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership (Instr. 4)	
					Code	v	(A)	(D)	Date Exercis	sable	Expiration Date	Title	Amou or Numb of Share	er					

Explanation of Responses:

- 1. The sale of shares was effected pursuant to a Rule 10b5-1 trading plan adopted on December 5, 2019.
- 2. This transaction was executed in multiple trades at prices ranging from \$27.14 to \$28.12. 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 39,815 shares of common stock that are subject to performance based vesting and 174,750 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorneyin-Fact 02/07/2020

** 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.