FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

STATEMENT	OF CHANGES	S IN BENEFICIA	L OWNERSHIP

l	OMB APPRO	VAL
	OMB Number:	3235-0287
l	Estimated average burd	en
l	hours per response:	0.5

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

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]								5. Relationship of Reporting Person(s) to Issuer (Check all applicable) X Director 10% Owner					
						3. Date of Earliest Transaction (Month/Day/Year) 07/01/2019								X Officer (give title below) Chairman, Pr			belov	,
(Street) BILLERI (City))1821 Zip)		- 4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)							i. Indivi.ine)					
Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																		
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Ye				Execution Date,		·			Acquired (A) or (D) (Instr. 3, 4 and		5) Secur Benef		ficially 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)	Price		Trans	action(s) 3 and 4)		(111511.4)	
Common Stock 07/01/201					019	.9			S	П	301(1)	D	\$34.63	69 ⁽²⁾	1,	176,894	D	
Common Stock 07/01/201				019	.9		S		2,844(1)	D	\$34.63	1.6369 ⁽²⁾		74,050 ⁽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)	Execution Date, if any		4. Transa Code (8)		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		6. Date Exe Expiration I (Month/Day		ate	7. Title and Amount of Securities Underlying Derivative Security (Instr. and 4)			vative ırity r. 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 Exer	cisable	Expiration Date	Title	Amount or Number of Shares					

Explanation of Responses:

- 1. The sales of shares reported on this Form 4 was made solely to cover tax obligations upon the vesting of 661 RSUs and 6,250 RSUs, respectively.
- 2. This transaction was executed in multiple trades at prices ranging from \$34.50 to \$34.8384. 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 215,266 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorneyin-Fact

07/03/2019

** 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.