FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL									
OMB Number:	3235-0287								
Estimated average burden									
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 (Last) (First) (Middle) C/O QUANTERIX CORPORATION, 113 HARTWELL AVENUE (Street)						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX] 3. Date of Earliest Transaction (Month/Day/Year) 10/02/2018 4. If Amendment, Date of Original Filed (Month/Day/Year)								5. Relationship of Reporting Person(s) to Issuer (Check all applicable) X Director 10% Owner X Officer (give title Other (specify below) Chairman, President & CEO 6. Individual or Joint/Group Filing (Check Applicable Line) X Form filed by One Reporting Person				
(City)			02421 (Zip)											71	Form filed by More than One Reporting Person			
		Tabl	le I - N	lon-Deriv	ative	Sec	uritie	s Ac	quire	d, Di	sposed o	f, or E	Benefici	ally	Owne	ed		
1. Title of Security (Instr. 3) 2. Transactio Date (Month/Day/N						Execution Date,		ate,	3. Transaction Code (Instr. 8) 4. Securities Acqui Disposed Of (D) (In						5. Amount of Securities Beneficially Owned Following Reported		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)		(11301. 4)
Common Stock 10/02/20					018)18			S		2,995(1)	D	\$20.9	520.905 ⁽²⁾		73,068 ⁽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)	rivative Conversion Date Execution or Exercise (Month/Day/Year) if any		ion Date,	4. Transaction Code (Instr. 8)		5. Nu of Deriv Secu Acqu (A) or Dispo of (D) (Instr and 5	rative rities ired r osed)	6. Date Exer Expiration D (Month/Day/		ate	7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4) Amount or Number of Shares		8. Price of Derivative Security (Instr. 5)		9. Number of derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	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 reported on this Form 4 was made solely to cover tax obligations upon the vesting of 6,250 RSUs and was effected pursuant to a Rule 10b5-1 trading plan adopted on August 14, 2018.
- 2. This transaction was executed in multiple trades at prices ranging from \$20.63 to \$21.23. 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 243,750 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorneyin-Fact 10/03/2018

** 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.