SEC Form 4	
------------	--

 \square

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

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

OMB APPROVAL								
OMB Number:	3235-0287							
Estimated average burg	len							
hours per response:	0.5							

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)					
HRUSU	<u>VSKIEKEV</u>	<u>11N</u>		X Director 10% Owner					
		EVIN (Middle) RPORATION,		X Officer (give title Other (specify					
(Last)	(First)	(Middle)	3. Date of Earliest Transaction (Month/Day/Year)	below) below)					
C/O QUAN	TERIX CORPO	RATION,	09/04/2019	Chairman, President & CEO					
900 MIDD	LESEX TURNPI	KE							
(Ctract)			4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Individual or Joint/Group Filing (Check Applicable Line)					
(Street)	A MA	01821		X Form filed by One Reporting Person					
				Form filed by More than One Reporting Person					
(City)	(State)	(Zip)		r croui					

Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned

1. Title of Security (Instr. 3)	2. Transaction Date (Month/Day/Year)	2A. Deemed Execution Date, if any (Month/Day/Year)	3. Transa Code (8)			Disposed Of (D) (Instr. 3, 4 and 5)		Securities Beneficially	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	Transaction(s) (Instr. 3 and 4)		(1150.4)
Common Stock	09/04/2019		S		3,362 ⁽¹⁾	D	\$25.971 ⁽²⁾	1,177,192 ⁽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. Deemed Execution Date, if any (Month/Day/Year)	4. Transa Code (8)		of		6. Date Exerc Expiration Da (Month/Day/)	7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		8. Price of Derivative Security (Instr. 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 Exercisable	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 7,123 RSUs.

2. This transaction was executed in multiple trades at prices ranging from \$25.92 to \$26.085. 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 210,365 restricted stock units.

Remarks:

<u>/s/ Brian Keane, as Attorney-</u>

in-Fact

09/05/2019

** 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.