## FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington D.C. 20549

STATEMENT	OF CHANGES	S IN BENEFICIAL	<b>OWNERSHIP</b>

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*					2. Issuer Name <b>and</b> Ticker or Trading Symbol Quanterix Corp [ QTRX ]								5. Relationship of Reporting Person(s) to Issuer (Check all applicable)					
HRUSOVSKY E KEVIN														X	X Director		10%	Owner
(Last)	(Fi	rst) (I	Middle	)	3. D	Date of Earliest Transaction (Month/Day/Year)								X Office below		er (give title w)	Othe belo	r (specify v)
C/O QUANTERIX CORPORATION,						01/15/2020							Chairman, President & CEO					
900 MIDDLESEX TURNPIKE																		
900 WIDDLESEX TORNFIRE				- 1 If	If Amendment, Date of Original Filed (Month/Day/Year)							- 1,	6. Individual or Joint/Group Filing (Check Applicable					
(Street)					7. "	4. II Americinent, Date of Original Filed (Month/Ddy/fedf)								Line)				
BILLERICA MA 01821												X Form filed by One Reporting Person  Form filed by More than One Reporting Person						
											.							
(City)	(St	ate) (2	Zip)												reis	OH		
Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																		
1. Title of Security (Instr. 3)  2. Transaction Date (Month/Day/Ye				ear) Execution Date, if any (Month/Day/Year)					Acquired (A) or (D) (Instr. 3, 4 and		<b>5</b> )	5. Amount of Securities		6. Ownership Form: Direct	7. Nature of Indirect			
			Year)			Code (Instr. 8)			. 5, 4 anu	, Be		ficially d Following	(D) or Indirect (I) (Instr. 4)	Beneficial Ownership				
							Code	v	Amount	(A) or (D)	Price		Repoi		(1) (111341. 4)	(Instr. 4)		
							$\dashv$		+		· /		(0)	`	<u> </u>			
Common Stock 01/15/202				)20	.0		S		6,500(1)	D	\$22.17	<sup>7</sup> 38 <sup>(2)</sup>	1,1	57,342 <sup>(3)</sup>	D			
Common Stock 01/17/202			)20	20		S		6,500(1)	D	\$24.	41	1,1	50,842(3)	D				
		Ta	ble II	I - Derivat	ive S	ecuri	ities	Acq	uired	l, Dist	oosed of,	or Be	neficial	ly Ov	vned		,	, , , , , , , , , , , , , , , , , , ,
											convertib							
1. Title of Derivative Security (Instr. 3)	Derivative Conversion Date Executity or Exercise (Month/Day/Year) if any		Execu if any	eemed Ition Date, Ith/Day/Year)  4. Transa Code (				6. Date Exercisable and Expiration Date (Month/Day/Year)		7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)					Ownership Form: Direct (D) or Indirect (I) (Instr. 4	Beneficial Ownership (Instr. 4)		
					Code	v	(A)	(D)	Date Exer	e rcisable	Expiration Date	Title	Amount or Number of Shares					

## **Explanation of Responses:**

- 1. The sale of shares reported on this Form 4 was made 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 \$22.05 to \$22.25. 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.
- $3. \ Includes \ 39,815 \ shares \ of common \ stock \ that \ are subject \ to \ performance \ based \ vesting \ and \ 181,873 \ restricted \ stock \ units.$

## Remarks:

/s/ Brian Keane, as Attorney-in-Fact

01/17/2020

\*\* Signature of Reporting Person

Date

Reminder: Report on a separate line for each class of securities beneficially owned directly or indirectly.

- $^{\star}$  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.