FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 205	49
----------------------	----

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*					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						Quanterra Corp [Q1KA]								X Direc		ctor		10% O	wner
(Last)	(Fi	rst) (N	Middle))	3. Da	Date of Earliest Transaction (Month/Day/Year)							\dashv	X	X Officer (give title below)		Other (sp below)		specify
C/O QUANTERIX CORPORATION,					05/0	05/01/2020									Chairman, President & CEO				
900 MIDDLESEX TURNPIKE																			
					4. If A	4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Individual or Joint/Group Filing (Check Applicable Line)					
(Street) BILLER	ICA M	A 0	1821											X	Form filed by One Reporting Person				
-															Form filed by More than One Reporting Person				
(City)	(S	ate) (Z	Zip)																
		Table	I - N	on-Deriva	tive S	Secur	rities	Ac	quire	d, Di	sposed of	, or E	enefic	ially	Own	ed			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Y				Year)	Execution Date,		· /	3. Transaction Code (Instr. 8) 4. Securities Acquired (A Disposed Of (D) (Instr. 3,					and 5) Secur Benet		rities Fo ficially (D) ed Following (I)		: Direct	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
									Code	v	Amount	(A) or (D)	Price	Transa		ction(s) 3 and 4)			(mounty)
Common Stock 05/01/202)20				S		3,399(1)	D	\$25.4	.4422 1,1		1,108,468(2)		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)	ative Conversion Date Execution Date, ity or Exercise (Month/Day/Year) if any				ransaction of code (Instr.) Sec (A) Dis of (i		osed) r. 3, 4	Expir	te Exer ation C th/Day/	/Year) Secu Unde Deriv		nt of ities lying ative ity (Instr. 4) Amount or	Deri Sec	rice of ivative urity tr. 5)	9. Number derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	Ownershi Form: Direct (D) or Indirec (I) (Instr. 4		Beneficial Ownership (Instr. 4)	
					Code	v	(A)	(D)	Date Exerc	cisable	Expiration Date	Title	Number of Shares						

Explanation of Responses:

- 1. The sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 7,123 RSUs.
- 2. Includes 192,378 restricted stock units and 39,815 shares of common stock that are subject to performance based vesting.

Remarks:

/s/ Brian Keane, as Attorneyin-Fact

05/04/2020

** 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.