FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

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

## STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

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

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						Quantity Corp [ ditty ]									X	Direc	tor		10% O	wner	
(Last)	(Fir	st) (N	√iddl	e)	3. D	Date of Earliest Transaction (Month/Day/Year)									X		Officer (give title pelow)		Other ( below)	specify	
C/O QUANTERIX CORPORATION,					07/2	07/21/2020									Chairman, President & CEO						
900 MIDDLESEX TURNPIKE																					
					4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)									6. Individual or Joint/Group Filing (Check Applicable						
(Street)															Line)  X Form filed by One Reporting Person						
BILLERICA MA 01821															Form filed by More than One Reporting						
(0)	(0)		·· 、													Perso			01.01.00	5. t.ig	
(City)	(St	ate) (2	Zip)																		
		Table	I - I	Non-Deriva	tive	Secui	rities	Ac	quir	ed, D	ispose	d of	f, or E	Benefic	ially	Own	ed				
1. Title of Security (Instr. 3)  2. Transaction Date (Month/Day/Yea			Execution Date,		,	3. Transaction Code (Instr. 8)  4. Securities A Disposed Of (I					and 5) S		5. Amount of Securities Beneficially Owned Following Reported		n: Direct or rect (I)	7. Nature of Indirect Beneficial Ownership					
								-	Code V		Amount		(A) or (D)	Price		Transa	saction(s) r. 3 and 4)		ıstr. 4)	(Instr. 4)	
Common Stock 07/21/20			07/21/202	20			S <sup>(1)</sup>		2,636		D	\$31.87	724 <sup>(2)</sup> 1,0		1,064,204 <sup>(3)</sup>		D				
	Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned																				
				(e.g., pu																	
1. Title of Derivative Security (Instr. 3)	erivative Conversion Date Execution Date, curity or Exercise (Month/Day/Year) if any			4. Transaction Code (Instr. 8)  5. Nun of Deriva Securi Acquii (A) or Dispos of (D) (Instr. and 5)		rative rities ired r osed ) : 3, 4	Exp (Mo	Date Exe Diration Onth/Day		and	7. Title and Amount of Securities Underlying Derivative Security (Ins 3 and 4)		Deri Sec (Ins	rice of vative urity tr. 5)	9. Number derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	Owner Form: Direct or Indi (I) (Ins	Ownership	11. Nature of Indirect Beneficial Ownership (Instr. 4)			
					Code	v	(A)	(D)	Dat Exe	te ercisable	Expir e Date	ation	Title	Number of Shares							

## **Explanation of Responses:**

- $1. \ The \ sale \ of \ shares \ was \ effected \ pursuant \ to \ a \ Rule \ 10b5-1 \ trading \ plan \ adopted \ on \ December \ 5, \ 2019.$
- 2. This transaction was executed in multiple trades at prices ranging from \$31.85 to \$31.98. 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 178,132 restricted stock units and 39,815 shares of common stock that are subject to performance based vesting.

## Remarks:

/s/ Brian Keane, as Attorney-

07/23/2020

in-Fact \*\* 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.