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(h)

## STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

**OMB APPROVAL** 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 [ Q110x ]										Direc	Director		10% O	wner			
(Last) (First) (Middle)							Date of Earliest Transaction (Month/Day/Year)										Officer (give title below)			Other ( below)				
C/O QUANTERIX CORPORATION,						07/02/2021											Chairman, President & CEO							
900 MIDDLESEX TURNPIKE																								
F							If Amendment, Date of Original Filed (Month/Day/Year)											6. Individual or Joint/Group Filing (Check Applicable						
(Street)						. ,	iaiii	ioni, bi	ato o	JI 011	giriari	nea (me	1101/10	ayi real	, I	Line)					ррисцые			
BILLER	ICA M	A 0	182	1												X	Form	Form filed by One Reporting Person						
,														Form filed by More than One Reporting Person										
(City)	(St	ate) (2	Zip)																					
Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																								
1. Title of Security (Instr. 3)				2. Transaction Date (Month/Day/Ye	ear)	2A. Deemed Execution Date, if any (Month/Day/Year)		Tr	3. Transaction Code (Instr. 8)		4. Securities Acquired (A) of Disposed Of (D) (Instr. 3, 4 a				5)	Securi Benefi Owned	5. Amount of Securities Beneficially Owned Following		wnership m: Direct or rect (I)	7. Nature of Indirect Beneficial Ownership				
								Co	ode	v	Amount		(A) or (D)	Price		Reported Transaction(s) (Instr. 3 and 4)		(Instr. 4)		(Instr. 4)				
Common	Stock		07/02/202	1				S		1,767	(1)	D	\$57.6396(2)		949,158 <sup>(3)</sup>			D						
Common	07/02/202	1					S		549(	1)	D	\$58.38	881 <sup>(4)</sup>	948,609(3)			D							
Common	07/02/202	.1					S		1,342	(1)	D	\$59.3432(5)		947,267 <sup>(3)</sup>			D							
		Tal	ole	II - Derivati													Owne	d						
				(e.g., pu	its, (	calls	, W	arrar	its,	opi	tions	, conv	ertii	ble se	curitie	:S)								
1. Title of Derivative Security (Instr. 3)	rivative Conversion Date Execution Date, Curity or Exercise (Month/Day/Year) if any					4. Transaction Code (Instr. 8)		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		Exp	iration	ercisable and I Date Iy/Year)		Amo Secu Unde Deriv	le and unt of rities rrlying rative rity (Instr 1 4)	Der Sec (Ins	rice of ivative urity tr. 5)	9. Number derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	y	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)			
					Code	Code V		(A) (	D)	Date Exercisab			ration	Title	Amount or 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 7,935 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$57.13 to \$58.11. 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 97,244 restricted stock units and 39,815 shares of common stock that are subject to performance based vesting.
- 4. This transaction was executed in multiple trades at prices ranging from \$58.17 to \$59.03. 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
- 5. This transaction was executed in multiple trades at prices ranging from \$59.17 to \$59.51. 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

## Remarks:

/s/ Brian Keane, as Attorney-

07/06/2021

in-Fact

\*\* Signature of Reporting Person

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.