FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPRO	DVAL
OMB Number:	3235-0287
Estimated average burd	den
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(h)

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* <u>HRUSOVSKY E KEVIN</u>						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]										olicable)	g Person(s) to I		
						3. Date of Earliest Transaction (Month/Day/Year) 01/29/2020									belov	*	Other below esident & CE		
(Street) BILLERICA MA 01821 (City) (State) (Zip)					- 4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)									5. Individual or Joint/Group Filing (Check Applicable Line) X Form filed by One Reporting Person Form filed by More than One Reporting Person				
	Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																		
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Ye						Execution Date,		·	3. Transa Code (8)		4. Securities Disposed Of			d 5) Secu Bene		ficially d Following	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership	
									Code	v	Amount	(A) or (D)	Price		Trans	action(s) 3 and 4)		(Instr. 4)	
Common Stock 01/29				01/29/20	20				S		6,500(1)	D	\$26.98	57 ⁽²⁾	1,1	37,842 ⁽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)	rive Conversion Date Execution Date, or Exercise (Month/Day/Year) if any			4. Transa Code (8)		5. Numof Derive Securion Acquired (A) or Disposof (D) (Instrument)	ative rities ired osed	6. Date Exercisable and Expiration Date (Month/Day/Year)			7. Title and Amount of Securities Underlying Derivative Security (Instr. and 4)				9. Number of derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)		
					Code	v	(A)	(D)	Date Exer	: cisable	Expiration Date	Title	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 \$26.70 to \$27.31. 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 Attorneyin-Fact 01/31/2020

** 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.