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(b).

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* HRUSOVSKY E KEVIN						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]								Check X	all app	tor 10% (on(s) to Is 10% O	wner
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						3. Date of Earliest Transaction (Month/Day/Year) 04/01/2021								X Officer (give the below) Chairman, President & CEO					·
(Street) BILLERICA MA 01821 (City) (State) (Zip)					4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)								i. Indiv ine) X	,				
	Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																		
1. Title of Security (Instr. 3)				2. Transaction Date (Month/Day/Ye	Execution		n Date,	Co	3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 ar			l 5) Securi Benefi		cially d Following		Direct	7. Nature of Indirect Beneficial Ownership (Instr. 4)
							Co	ode	v	Amount	(A) or (D)	Price	Transa		action(s) 3 and 4)	(iiisti.	· - "	(111501.4)	
Common	Stock		04/01/202	1				S		1,024(1)	D	\$61.91	96(2) 98		7,976 ⁽³⁾		D		
Common Stock				04/01/202	1				S		2,148(1)	D	\$62.69	84(4)	985,828(3)			D	
Common Stock				04/01/202	21				S		260(1)	D	\$63.3	\$63.36		985,568(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) 2. Conversion of Exercise Price of Derivative Security (Instr. 3) 3. Transaction Date (Month/Day/Year) (Month/Day/Year) 3A. Deemed Execution Date, if any (Month/Day/Year)				4. Transaction Code (Instr. 8)		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		Exp	Pate Exe piration onth/Day		Amo Secu Unde Deriv	cle and unt of urities erlying vative urity (Instr. d 4)	Deri	vative urity tr. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s (Instr. 4)	/ O F D 0 (I	0. Ownership Form: Direct (D) Ir Indirect I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)	
					Code	Code V (A)		(D)	Date Exercisable		Expiration e Date	n 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 \$61.30 to \$62.27. 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 121,049 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 \$62.30 to \$63.21. 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-

04/05/2021

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.