FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

STATEMENT	ΩF	CHANGES	IN BENEFI	CIAI	OWNERS	HIP
SIAIEMENI	UГ	CHANGES	IIN DEINELI	CIAL	OVVIVERS	HIL

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					Quanterry Corb [Grky]								[`	X Direc		,		wner		
(Last) (First) (Middle)					3. D	Date of Earliest Transaction (Month/Day/Year)								\dashv	X	X Officer (give title below)			Other (specify below)	
C/O QUANTERIX CORPORATION,					01/	01/03/2022								Chairman and CEO						
900 MIDDLESEX TURNPIKE																				
					4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)									6. Individual or Joint/Group Filing (Check Applicable Line)					
(Street) BILLERICA MA 01821														X						
														Form filed by More than One Reporting Person					orting	
(City)	(St	ate) (Z	<u>z</u> ip)																	
		Table	I - I	Non-Deriva	tive	Secui	rities	Acc	quir	ed, D	Disp	osed o	f, or E	Benefic	ially	Own	ed			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Year				ar) E	2A. Deemed Execution Da if any (Month/Day/)		Date, Ti		3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 ar			nd 5) Secur Benet		cially I Following	Form (D) o	n: Direct r ect (I)	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
							C		ode	v	Am	ount	(A) or (D)	Price		Transa	Transaction(s) (Instr. 3 and 4)		1. 4)	(11150.4)
Common Stock 01/			01/03/202	!2				S		3,715 ⁽¹⁾		D	\$41.9066(2)		890,386(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 or Exercise Price of Derivative Security	ercise (Month/Day/Year) if any (Month/Day/Year)				de (Instr. Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)			Expiration Date (Month/Day/Year)			Amou Secul Unde Deriv Secul 3 and	7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4) Amount or Number of Title Shares		rice of vative urity tr. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	,	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)	

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 \$41.71 to \$42.16. 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 49,634 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorneyin-Fact

01/05/2022

** 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.