FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

· · · · · · · · · · · · · · · · · · ·	OMB APPROVAL			
STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP	OMB Number:	3235-0287		
OTATEMENT OF OTTATOLO IN BEINE TOTAL OWNERORIII	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)				
HRUS(OVSKY I	E KEVIN			וַעַנ	ıante	<u> FIX C</u>	<u>югр</u>	Į QT	[QTRX]				X	Direc	,	10% (Owner
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						3. Date of Earliest Transaction (Month/Day/Year) 12/02/2019								X Officer (give titl below) Chairman,			Other below esident & CE	'
(Street) BILLERICA MA 01821				- 4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Individual or Joint/Group Filing (Check Applicable Line) X Form filed by One Reporting Person Form filed by More than One Reporting Person					
(City)	(Si	ate) (Zip)															
		Tabl	e I - I	Non-Deriv	ative	Sec	uritie	s Ad	cquire	ed, Di	isposed o	f, or E	Benefici	ally (Owne	ed		
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Yell)						Execution Date,						Acquired (A) or (D) (Instr. 3, 4 and		5) Secur Benef Owne		ficially d Following	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
									Code	v	Amount	(A) or (D)	Price		Reported Transaction(s) (Instr. 3 and 4)			(iiisii. 4)
Common Stock 12/02/201						19					3,357(1)	D	\$25.06	29 ⁽²⁾	1,1	67,163 ⁽³⁾	D	
		Та	ıble II							, .	oosed of, convertib			•	ned		•	,
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (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)		6. Date Exercisable and Expiration Date (Month/Day/Year)			7. Title and Amount of Securities Underlying Derivative Security (Instr. and 4)		8. Pri Deriv Secu (Instr	rative rity 7. 5)	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	of Shares					

Explanation of Responses:

- $1. \ The sales of shares reported on this Form 4 was made solely to cover tax obligations upon the vesting of 7,123 RSUs.$
- 2. This transaction was executed in multiple trades at prices ranging from \$24.99 to \$25.13. 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 188,996 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorney-12/03/2019

** 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.