FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington,	D.C.	20549
-------------	------	-------

	OMB APPROVAL								
	OMB Number:	3235-0287							
l	Estimated average burden								
l	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* HRUSOVSKY E KEVIN (Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE (Street) BILLERICA MA 01821 (City) (State) (Zip)						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX] 3. Date of Earliest Transaction (Month/Day/Year) 02/03/2020 4. If Amendment, Date of Original Filed (Month/Day/Year)								Check X X	all app Direct Office below Ch	onship of Reporting Person(s) to Issuer Il applicable) Director 10% Owner Officer (give title Other (specify below) below) Chairman, President & CEO ual or Joint/Group Filing (Check Applicable Form filed by One Reporting Person Form filed by More than One Reporting Person		
Table I - Non-Deriva 1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Ye					on	2A. Deemed Execution Date,			3. 4. Securities		4. Securities A	Acquired (A) or (D) (Instr. 3, 4 and 5)			5. Amount of Securities Beneficially Owned Following Reported Transaction(s) (Instr. 3 and 4)		6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
Common Stock 02/03/20)20	0			S		5,900(1)	D	\$27.04	56 ⁽²⁾	<u> </u>	131,942	D	
Common Stock 02/03/20								4	S	\square	600(1)	D	\$27.	_	1,131,342		D	
1. Title of Derivative Conversion Security or Exercise (Month/Day/Year) if any		3A. Do Execu		4. Transa Code (ve Securi Its, calls, 4. Transaction Code (Instr.		Jarrants, op. 5. Number of for Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		s 2,409(3) iired, Disposed of, options, convertib 6. Date Exercisable and Expiration Date (Month/Day/Year)		7. Title Amour Securit Underl Derivat	e securities) 7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4) Amount or Number of		1,128,933 ⁽⁴⁾ Price of derivative surity Securities Beneficiall Owned Following Reported Transactio (Instr. 4)		Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)	

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 \$26.44 to \$27.40. 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. The sales of shares reported on this Form 4 was made solely to cover tax obligations upon the vesting of 7,123 RSUs.
- 4. Includes 39,815 shares of common stock that are subject to performance based vesting and 174,750 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorney-

02/05/2020

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.