FORM 4

obligations may continue. See Instruction 1(b).

Check this box if no longer subject to Section 16. Form 4 or Form 5

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

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

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL OMB Number: 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]									all app Dired	olicable) ctor	10%	Person(s) to Issuer 10% Owner	
•	`	CORPORATION	Middle) N,			3. Date of Earliest Transaction (Month/Day/Year) 11/16/2018									belov	,	belo	Other (specify below) sident & CEO	
(Street) LEXING (City))2421 Zip)		- 4. If	f Amendment, Date of Original Filed (Month/Day/Year)								6. Indi Line) X	ividual or Joint/Group Filing (Check Applicable Form filed by One Reporting Person Form filed by More than One Reporting Person				
		Tab	e I - N	lon-Deri	vative	Sec	uritie	s Ac	quire	d, Di	sposed o	f, or E	Benefic	cially	Own	ed			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/N					Execution Date,			3. Transaction Code (Instr. 8) 4. Securities Disposed Of					nd 5) Secu Bene		icially 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		Trans	action(s) 3 and 4)		(111501.4)	
Common Stock 11/16				11/16/2	2018)18			S		3,940(1)	D	\$16.	\$16.663 ⁽²⁾		66,299(3)	D		
Common	Stock			11/16/2	2018				S		60(1)	D	\$1	17.12 1,166,239 ⁽³⁾ D					
		Ta	ble II								osed of, convertib				wned				
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	se (Month/Day/Year)	Execut if any			ection Instr.	5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		Expira	e Exerc tion D h/Day/		7. Title and Amount of Securities Underlying Derivative Security (Instr. and 4)		Deri Sec (Ins	ivative urity tr. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s (Instr. 4)	Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership (Instr. 4)	
					Code	v	(A)	(D)	Date Exerci	isable	Expiration Date	Title	or Number of Shares	r					

Explanation of Responses:

- 1. The sale of shares was effected pursuant to a Rule 10b5-1 trading plan adopted on June 14, 2018.
- 2. This transaction was executed in multiple trades at prices ranging from \$16.031 to \$17.00. 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 237,500 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorney-

11/20/2018

** 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.