FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Check this box if no longer subject
to Section 16. Form 4 or Form 5
obligations may continue. See
Instruction 1(h)

## 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 <b>and</b> Ticker or Trading Symbol Quanterix Corp [ QTRX ]									all app		ng Per	10% O	wner	
(Last) (First) (Middle)  C/O QUANTERIX CORPORATION,  900 MIDDLESEX TURNPIKE						3. Date of Earliest Transaction (Month/Day/Year) 03/10/2020									belov	er (give title v) airman, Pro	eside	Other ( below) nt & CEC		
(Street) BILLERICA MA 01821 (City) (State) (Zip)					4. If A	4. If Amendment, Date of Original Filed (Month/Day/Year)								. Indivine)	′					
Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																				
1. Title of Security (Instr. 3)  2. Transaction Date (Month/Day/						Execution Date,			3. Transaction Code (Instr. 8)  4. Securities Acquired (A) Disposed Of (D) (Instr. 3, 4)					and 5) Securi Benefi		cially d Following	Form (D) o		7. Nature of Indirect Beneficial Ownership (Instr. 4)	
									Code	v	Amount	(A) o (D)	Price		Transa	action(s) 3 and 4)			(IIISU. 4)	
Common Stock 03/10/20						)20					5,500(1)	D	\$20.5	.53(2) 1,14		40,600(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)	erivative Conversion Date Execution Date, curity or Exercise (Month/Day/Year) if any			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 (Inst 3 and 4)		t		9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s) (Instr. 4)		10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)		
					Code	v	(A)	(D)		isable	Date	Title	Shares							

## **Explanation of Responses:**

- $1.\ The\ sale\ of\ shares\ was\ effected\ pursuant\ to\ a\ Rule\ 10b5-1\ trading\ plan\ adopted\ on\ December\ 5,\ 2019.$
- 2. This transaction was executed in multiple trades at prices ranging from \$20.37 to \$20.54. 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 206,624 restricted stock units and 39,815 shares of common stock that are subject to performance based vesting.

## Remarks:

/s/ Brian Keane, as Attorney-

03/12/2020

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.