FORM 4

## **UNITED STATES SECURITIES AND EXCHANGE COMMISSION**

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

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL									
OMB Number: 3235-0287									
Estimated average burden									
hours per response: 0									

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

Name and Address of Reporting Person*     Hlavinka Sarah E.					2. Issuer Name <b>and</b> Ticker or Trading Symbol  Quanterix Corp [ QTRX ]									all app Direc	tor	ng Pers	10% O	wner	
(Last) (First) (Middle) C/O QUANTERIX CORPORATION					3. Date of Earliest Transaction (Month/Day/Year) 07/01/2020										Officer (give title below)		Other (s below)		specify
900 MIDDLESEX TURNPIKE					4. If Amendment, Date of Original Filed (Month/Day/Year)							6. Individual or Joint/Group Filing (Check Applicable Line)							
(Street) BILLERICA MA 01821														X	<b>,</b>				
(City)	(5	tate) (2	Zip)																
	Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																		
1. Title of Security (Instr. 3)  2. Transact Date (Month/Day				Execution Date,			3. Transaction Code (Instr. 8)  4. Securities Acquired (Disposed Of (D) (Instr. 3)			iired (A) nstr. 3, 4	or I and	Securi Benefi Owned	5. Amount of Securities Beneficially Dwned Following Reported		: Direct	7. Nature of Indirect Beneficial Ownership			
									Code	v	Amount	(A) (D)	(A) or (D) Pri		Transa	Transaction(s) (Instr. 3 and 4)			(Instr. 4)
Common Stock 07/01/2				2020		A		359(1)	A	\$2	7.8(2)	.8 <sup>(2)</sup> 8,441 <sup>(3)</sup>		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)	rivative conversion or Exercise (Month/Day/Year) Execution Date, (month/Day/Year) Frice of Derivative Security			nsaction de (Instr.  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)  Date Expiration Exercisable Date			7. Title and Amount of Securities Underlying Derivative Security (Insi 3 and 4)  Amou or Numbo of Title Share:		Der Sec (Ins	Price of rivative curity str. 5)	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)		

## **Explanation of Responses:**

- 1. Consists of stock granted in lieu of cash fees for service on the Company's Board of Directors and committees thereof for the second quarter of 2020, pursuant to an election made in December 2019.
- 2. Closing price of the Company's common stock on the Nasdaq Global Market on July 1, 2020.
- 3. Includes 6,810 restricted stock units.

## Remarks:

/s/ Brian Keane, as Attorney-

07/06/2020

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.