SEC Form 4

FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

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

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934 or Section 30(h) of the Investment Company Act of 1940 OMB APPROVAL
OMB Number: 3235-0287
Estimated average burden
hours per response: 0.5

			2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]		ationship of Reporting Person(s) to Issuer k all applicable)							
<u> Hidvilika Sal</u>	<u>dll E.</u>			X	Director	10% Owner						
(Last) C/O QUANTEF	(First) RIX CORPORATIO	(Middle) ON,	3. Date of Earliest Transaction (Month/Day/Year) 07/25/2019		Officer (give title below)	Other (specify below)						
900 MIDDLESEX TURNPIKE			4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Individual or Joint/Group Filing (Check Applicable Line)								
(Street)				X	Form filed by One Reporting Person							
BILLERICA	MA	01821			Form filed by More than Person	One Reporting						
(City)	(State)	(Zip)										

Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned

1. Title of Security (Instr. 3)	2. Transaction Date (Month/Day/Year)	2A. Deemed Execution Date, if any (Month/Day/Year)	te, Transaction Code (Instr.		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 and 5)			5. Amount of Securities Beneficially Owned Following Reported	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	Transaction(s) (Instr. 3 and 4)		(1130.4)	
Common Stock	07/25/2019		A		4,540 ⁽¹⁾⁽²⁾	Α	\$ <mark>0</mark>	4,540	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)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	3A. Deemed Execution Date, if any (Month/Day/Year)	4. Transa Code (8)		Derivative		6. Date Exerc Expiration Da (Month/Day/Y	ate	7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		8. Price of Derivative Security (Instr. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s)	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)
				Code	v	(A)	(D)	Date Exercisable	Expiration Date	Title	Amount or Number of Shares		(Instr. 4)		
Stock Option (right to buy)	\$31.91	07/25/2019		A			15,800 ⁽¹⁾	(3)	07/25/2029	Common Stock	15,800	\$0	15,800	D	

Explanation of Responses:

1. Initial equity grant under the Quanterix Corporation 2018 Non-Employee Director Compensation Policy upon Ms. Hlavinka's election to the Quanterix Board of Directors.

2. Represents shares of common stock issuable upon vesting of restricted stock units. The restricted stock units vest as to one-third of the shares on each of July 25, 2020, July 25, 2021 and July 25, 2022.

3. The option vests as to one-third of the shares on July 25, 2020, and the remaining two-thirds of the shares over the following two years in 24 successive equal monthly installments at the end of each month until July 25, 2022.

<u>/s/ Brian Keane, Attorney-in-</u>

fact

** Signature of Reporting Person

Reminder: Report on a separate line for each class of securities beneficially owned directly or indirectly.

 \ast 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.

<u>07/25/2019</u>

Date