## FORM 4

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

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549 STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL

l	OMB Number:	3235-0287									
l	Estimated average burden										
1	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

			or Section 30(n) of the investment Company Act of 1940					
1	dress of Reporting	Person <sup>*</sup>	2. Issuer Name <b>and</b> Ticker or Trading Symbol Quanterix Corp [ QTRX ]	5. Relationship of Reporting Person(s) to Issuer (Check all applicable)				
Madaus M			-<	X Director 10% Owner				
			—	Officer (give title Other (specify				
(Last)	(First)	(Middle)	3. Date of Earliest Transaction (Month/Day/Year) 10/19/2020	below) below)				
C/O QUANT	QUANTERIX CORPORATION,		10/19/2020					
900 MIDDLESEX TURNPIKE				6. Individual or Joint/Group Filing (Check Applicable Line)				
			4. If Amendment, Date of Original Filed (Month/Day/Year)					
(Street)				X Form filed by One Reporting Person				
BILLERICA	MA	01821		Form filed by More than One Reporting Person				
(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)	3. Transa Code ( 8)				(A) or 3, 4 and	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	10/19/2020		М		1,600	A	\$2.28	13,316(1)	D	
Common Stock	10/19/2020		S		1,600 <sup>(2)</sup>	D	\$44.59	11,716 <sup>(1)</sup>	D	
Common Stock	10/20/2020		М		1,600	A	\$2.28	13,316(1)	D	
Common Stock	10/20/2020		S		1,600(2)	D	\$42.57	11,716 <sup>(1)</sup>	D	
Common Stock								164,966	I	See Footnote <sup>(3)</sup>

# Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities)

			(0.9.,	pu.o,	ouno	,	mante	s, options			111100)				
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. 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 (Instr. 3 and 4)		8. Price of Derivative Security (Instr. 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)
				Code	v	(A)	(D)	Date Exercisable	Expiration Date	Title	Amount or Number of Shares				
Stock Option (Right to Buy)	\$2.28	10/19/2020		М			1,600	(4)	10/31/2021	Common Stock	1,600	\$0.00	4,572	D	
Stock Option (Right to Buy)	\$2.28	10/20/2020		М			1,600	(4)	10/31/2021	Common Stock	1,600	\$0.00	2,972	D	

Explanation of Responses:

1. Includes 2,270 Restricted Stock Units.

2. All sales reported in this Form 4 were effected pursuant to a Rule 10b5-1 trading plan adopted by the Reporting Person on September 9, 2019.

3. Held by a trust for which the Reporting Person is a trust advisor who shares voting and investment control.

4. All options are vested and exercisable.

#### Remarks:

/s/ Brian Keane, Attorney-in-Fact

10/20/2020

\*\* 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.