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

Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL							
OMB Number	3235-028						

Estimated average burden hours per response: 0.5

			or Section 30(h) of the Investment Company Act of 1940							
Name and Address of Reporting Person* Madaus Martin D			2. Issuer Name and Ticker or Trading Symbol <u>Quanterix Corp</u> [QTRX]	(Check	5. Relationship of Reporting Person(s) to Issuer (Check all applicable)					
				X	Director	10% Owner				
(Last) C/O QUANTER		(Middle)	3. Date of Earliest Transaction (Month/Day/Year) 12/10/2019		Officer (give title below)	Other (specify below)				
113 HARTWELL AVENUE			4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Indiv	6. Individual or Joint/Group Filing (Check Applicable					
(Street) LEXINGTON	MA	02421		X	Form filed by One Re	. 0				
(City)	(State)	(Zip)			Person					
Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned										

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)				5. Amount of Securities Beneficially Owned Following Reported	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership	
			Code	v	Amount	(A) or (D)	Price	Transaction(s) (Instr. 3 and 4)		(Instr. 4)
Common Stock	12/10/2019		M		1,600	A	\$1.93	8,149	D	
Common Stock	12/10/2019		S ⁽¹⁾		1,600	D	\$24.81	6,549	D	
Common Stock	12/10/2019		S ⁽¹⁾		2,300	D	\$25.1274 ⁽²⁾	213,466	I	See Footnote ⁽³⁾
Common Stock	12/11/2019		M		1,600	A	\$1.93	8,149	D	
Common Stock	12/11/2019		S ⁽¹⁾		1,600	D	\$25.25	6,549	D	
Common Stock	12/11/2019		S ⁽¹⁾		2,300	D	\$25.3571 ⁽⁴⁾	211,166	I	See footnote ⁽³⁾

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)	Code (nsaction of Ex		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)	\$1.93	12/10/2019		M			1,600	(5)	01/05/2021	Common Stock	1,600	\$0.00	63,307	D	
Stock Option (Right to Buy)	\$1.93	12/11/2019		M			1,600	(5)	01/05/2021	Common Stock	1,600	\$0.00	61,707	D	

Explanation of Responses:

- 1. All sales reported in this Form 4 were effected pursuant to Rule 10b5-1 trading plans adopted by the Reporting Person on September 9, 2019.
- 2. This transaction was executed in multiple trades at prices ranging from \$24.93 to \$25.26. 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. Held by a trust for which the Reporting Person is a trust advisor who shares voting and investment control.
- 4. This transaction was executed in multiple trades at prices ranging from \$24.96 to \$25.77. 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
- 5. All options are vested and exercisable.

Remarks:

/s/ Brian Keane, Attorney-in-

12/11/2019

<u>fact</u>

** 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.	