SEC Form 4	
------------	--

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

	OMB Number:	3235-0287
	Estimated average burde	en
I	hours per response:	0.5

or Section 30(h) of the Investment Company Act of 1940			ne Securities Exchange Act of 1 tment Company Act of 1940	934
--	--	--	--	-----

			of Section So(n) of the investment Company Act of 1940						
1. Name and Add	1 0	Person*	2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]	5. Relationship of Reporting Person(s) to Issuer (Check all applicable)					
<u>Madaus Martin D</u>				X Director 10% Owner					
				Officer (give title Other (specify					
			3. Date of Earliest Transaction (Month/Day/Year)	below) below)					
C/O QUANTE	RIX CORPOR	RATION,	05/27/2020						
900 MIDDLES	SEX TURNPIP	Œ							
			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 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)		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	
			Code	v	Amount	(A) or (D)	Price	Transaction(s) (Instr. 3 and 4)		(Instr. 4)
Common Stock	05/27/2020		М		1,600	A	\$1.93	12,145(1)	D	
Common Stock	05/27/2020		S ⁽²⁾		1,600	D	\$26.02	10,545(1)	D	
Common Stock	05/28/2020		М		1,600	A	\$1.93	12,145(1)	D	
Common Stock	05/28/2020		S ⁽²⁾		1,600	D	\$26.51	10,545(1)	D	
Common Stock								164,966	I	See Footnote ⁽³⁾

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

	(eigi, pulo, bulo, bulo, bulois, bonventible sedunices)														
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)		of Deri Sec Acq (A) Disp of (I	rivative (Month/Day/Year) curities quired or sposed (D) str. 3, 4		te Amount of Securities Underlying		Amount of Securities Inderlying Derivative Security		9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s) (Instr. 4)	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	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	05/27/2020		М			1,600	(4)	01/05/2021	Common Stock	1,600	\$0.00	24,907	D	
Stock Option (Right to Buy)	\$1.93	05/28/2020		М			1,600	(4)	01/05/2021	Common Stock	1,600	\$0.00	23,307	D	

Explanation of Responses:

1. Includes 2,270 Restricted Stock Units.

2. 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.

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

05/29/2020

** 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.