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

Vashington,	D.C.	20549
-------------	------	-------

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL OMB Number: 3235-0287 Estimated average burden 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

1. Name and Address of Reporting Person* Madaus Martin D					2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]								5. Relationship of Report (Check all applicable) X Director			10% Own			
(Last) (First) (Middle) C/O QUANTERIX CORPORATION,						3. Date of Earliest Transaction (Month/Day/Year) 06/24/2020								Officer below)	(give title	Э	Other 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) BILLER	ICA M	'A	01821			I '									X Form filed by One Reporting Person Form filed by More than One Reporting Person				
(City)	(S	tate)	(Zip)																
		Tab	le I - No	on-Deriv	/ative	Sec	urit	ies Ac	quired	l, Di	sposed o	of, or Be	neficia	lly Owned	d				
		2. Transaction Date (Month/Day/Year)		2A. Deemed Execution Date, if any (Month/Day/Year)				4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 and 5)			Beneficially Owned Following		6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)		7. Nature of Indirect Beneficial Ownership				
						Code	v	Amount	(A) or (D)	Price	Reported Transacti (Instr. 3 a	on(s)			(Instr. 4)				
Common Stock 06/24			06/24/	2020				М		1,600	A	\$1.93	12,1	45 ⁽¹⁾	D				
Common Stock			06/24/	06/24/2020		0		S ⁽²⁾		1,600	D	\$26.39	10,5	45 ⁽¹⁾	D				
Common Stock			06/25/	06/25/2020)		M		1,600	A	\$1.93	12,1	45 ⁽¹⁾	D				
Common Stock			06/25/	25/2020				S (2)		1,600	D	\$26.08	3 10,5	45(1)	D				
Common Stock													164,	966		I	See Footnote ⁽³⁾		
		Т	able II								posed of converti			/ Owned					
1. Title of Derivative Security (Instr. 3)	titive Conversion Date Execution Date, Tr ity or Exercise (Month/Day/Year) if any C				ransaction code (Instr. 5)		n of		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. Ownersh Form: Direct (D or Indirec (I) (Instr.	Beneficial Ownership t (Instr. 4)		
					Code	v	(A)		Date Exercisa	able	Expiration Date	Title	Amount or Number of Shares						
Stock Option (Right to Buy)	\$1.93	06/24/2020			M			1,600	(4)		01/05/2021	Common Stock	1,600	\$0.00	18,5	07	D		
Stock Option (Right to	\$1.93	06/25/2020			M			1,600	(4)		01/05/2021	Common Stock	1,600	\$0.00	16,9	07	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-

06/26/2020

Fact

** Signature of Reporting Person

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.