FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington,	D.C.	20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL									
OMB Number: 3235-028									
Estimated average burden									
hours	ner response	e: 0.5							

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

Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934 or Section 30(h) of the Investment Company Act of 1940

							00(.	.,		-	ompany 7 tot	0. 20.0								
1. Name and Address of Reporting Person* Madaus Martin D (Last) (First) (Middle) C/O QUANTERIX CORPORATION,						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]								5. Relationship of Reporting Person(s) to Issuer (Check all applicable)						
														X Director		10% Owner		wner		
						3. Date of Earliest Transaction (Month/Day/Year) 05/12/2020									(give title		Other (below)	specify		
900 MIDDLESEX TURNPIKE					4. 11	4. If Amendment, Date of Original Filed (Month/Day/Year)									6. Individual or Joint/Group Filing (Check Applicable					
(Street) BILLER	ICA M	'A	01821											Line) 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	_			quired	l, Di	sposed o	of, or Be	neficia	lly Owne	d					
= This or county (monto)		2. Transaction Date (Month/Day/Year)		2A. Deemed Execution Date, if any (Month/Day/Year)		Transaction Disposed Code (Instr. 5)		Disposed (ies Acquired (A) or Of (D) (Instr. 3, 4 and		5. Amount of Securities Beneficially Owned Following		6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Direct Ir Indirect B tr. 4) C	7. Nature of Indirect Beneficial Ownership					
							Code	v	Amount	(A) or (D)	Price	Reported Transacti (Instr. 3 a	on(s)		(1)	nstr. 4)				
Common Stock 0		05/12/	2/2020				M		1,600	A	\$1.93	3 12,1	45 ⁽¹⁾	D						
Common	Common Stock 05/1			05/12/	2020				S ⁽²⁾		1,600	D	\$27.8	6 10,5	10,545(1)		D			
Common	Stock			05/13/	2020				М		1,600	A	\$1.93	3 12,1	12,145(1)		D			
Common	Stock			05/13/	2020				S (2)		1,600	D	\$26.	5 10,5	10,545(1)		D			
Common	Stock													164,966 I				ee ootnote ⁽³⁾		
		Т	able II								oosed of converti			y Owned		,				
1. Title of Derivative Security (Instr. 3) 2. Conversion or Exercise Price of Derivative Security		3. Transaction Date (Month/Day/Year)	Execution if any	Execution Date, f any		4. Transaction Code (Instr. 8)		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. Numb derivativ Securitie Benefici Owned Followin Reporter Transact (Instr. 4)	e Ownersh Form: Direct (D or Indirect (I) (Instr.		11. Nature of Indirect Beneficial Ownership (Instr. 4)		
					Code	٧	(A)		Date Exercisa	ıble	Expiration Date	Title	Amount or Number of Shares	1						
Stock Option (Right to Buy)	\$1.93	05/12/2020			М			1,600	(4)		01/05/2021	Common Stock	1,600	\$0.00	28,1	07	D			
Stock Option (Right to	\$1.93	05/13/2020			M			1,600	(4)		01/05/2021	Common	1,600	\$0.00	26,5	07	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-

05/14/2020

<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.