FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington,	D.C.	20549	

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL								
OMB Number:	3235-0287							
Estimated average burden								
hours per response.	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

Name and Address of Reporting Person*     Madaus Martin D					2. 19	Section 30(f) of the investment Company Act of 1940  2. Issuer Name and Ticker or Trading Symbol  Quanterix Corp [ QTRX ]								Relationship of Reporting Person(s) to Issuer (Check all applicable)     X Director 10% Owner					
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						3. Date of Earliest Transaction (Month/Day/Year) 07/08/2020									Officer (give title Other (specify below) below)				
(Street) BILLERI	ICA M	A	01821 (Zip)		4. 11	ndmer	nt, Date	e of Original Filed (Month/Day/Year)					6. Individual or Joint/Group Filing (Check Applicable Line)  X Form filed by One Reporting Person  Form filed by More than One Reporting Person						
(City)	(5)	,	,	on-Deriv	/ative	Sec	urit	ies Ac	auired	 I. Di	sposed o	of. or Be	neficia	ally Owne	d				
1. Title of Security (Instr. 3)		2. Transaction Date (Month/Day/Year)		2A. Deemed Execution Date,		3. 4. Securiting Disposed (Code (Instr. 5)		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)		7. Nature of Indirect Beneficial Ownership					
								Code	v	Amount	(A) or (D)	Price	Reported Transacti (Instr. 3 a	on(s)			(Instr. 4)		
Common Stock			07/08/	08/2020		20		M		1,600	A	\$1.9	3 12,7	12,796(1)		D			
Common Stock		07/08/	08/2020		20		<b>S</b> <sup>(2)</sup>		1,600	D	\$27	11,1	11,196(1)		D				
Common	Common Stock		07/09/	07/09/2020				M		1,600	A	\$1.9	3 12,7	12,796(1)		D			
Common	Stock			07/09/	2020	_			S <sup>(2)</sup>		1,600	D	\$26.8	11,1					
Common	Common Stock											164	164,966		I	See Footnote <sup>(3)</sup>			
		Т	able II						,		posed of converti	,		ly Owned		•			
1. Title of Derivative Security (Instr. 3)	e of 2. 3. Transaction Bate Execution Date Execution Date, if any		n Date,	4.		n of		6. Date Exercisa Expiration Date (Month/Day/Year		te	Amount of		8. Price of Derivative Security (Instr. 5)			10. Ownersh Form: Direct (D or Indire (I) (Instr.	Beneficial Ownershi ct (Instr. 4)		
					Code	v	(A)	(D)	Date Exercisa	ıble	Expiration Date	Title	Amoun or Number of Shares						
Stock Option (Right to Buy)	\$1.93	07/08/2020		М				1,600	(4)		01/05/2021	Common Stock	1,600	\$0.00	00 15,307		D		
Stock Option (Right to Buy)	\$1.93	07/09/2020			М			1,600	(4)		01/05/2021	Common Stock	1,600	\$0.00	13,707		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-

07/10/2020

**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.