FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

Washington, D.C. 20040	

OMB APPROVAL OMB Number: Estimated average burden

0.5

hours per response:

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(.				Company 7 tot	. 01 10-10									
Name and Address of Reporting Person*     Madaus Martin D						2. Issuer Name <b>and</b> Ticker or Trading Symbol Quanterix Corp [ QTRX ]									5. Relationship of Reporting Person(s) to Issuer (Check all applicable)						
Maddus Martin D															X Director		10% Owner		wner		
(Last) (First) (Middle) C/O QUANTERIX CORPORATION,					3. Date of Earliest Transaction (Month/Day/Year) 01/22/2020									Officer below)	(give title		Other ( below)	specify			
900 MIDDLESEX TURNPIKE					4. 1	4. If Amendment, Date of Original Filed (Month/Day/Year)									6. Individual or Joint/Group Filing (Check Applicable						
(Street) BILLERICA MA 0182			L		Line) X Fo									Form f	Form filed by One Reporting Person Form filed by More than One Reporting Person						
(City)	(S	state)	(Zip)																		
		Tab	le I -	Non-Deri	vative	Sec	urit	ies A	cquir	ed, C	Disposed (	of, or B	enefic	ially	Owned	i					
1. Title of Security (Instr. 3)				2. Transacti Date (Month/Day		Execut		Date,	3. Transaction Code (Instr. 8)			Acquired (A) or (D) (Instr. 3, 4 and		5)		es ally Following	Form:	Direct I Indirect I str. 4)	7. Nature of ndirect Beneficial Ownership		
									Code	v	Amount	(A) or (D)	Price		Reported Transact (Instr. 3 a	ion(s)			Instr. 4)		
Common Stock				01/22/2020				M		1,600	A	\$1.9	93	11,202			D				
Common Stock				01/22/2020				S <sup>(1)</sup>		1,600	D	\$25.	5.15 9,602		502		D				
Common Stock				01/22/2020				S <sup>(1)</sup>		2,300	D	\$25.42	5.428 <sup>(2)</sup> 199,66		,666			See cootnote <sup>(3)</sup>			
Common Stock				01/23/2020				M		1,600	A	\$1.9	\$1.93 11,20		202		D				
Common Stock				01/23/2020				S <sup>(1)</sup>		1,600	D	\$24.	524.86 9,6		602		D				
Common Stock				01/23/2020				S <sup>(1)</sup>		2,300	D	\$25.59	\$25.5978 <sup>(4)</sup>		197,366			See cootnote <sup>(3)</sup>			
		7	Гable								sposed of				Owned						
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	if any	eemed ution Date,	4. Transa	I. Fransaction Code (Instr.		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)			cisable and Date	7. Title and Amount of Securities Underlying Derivative Secu (Instr. 3 and 4)		8 D S	Price of erivative ecurity nstr. 5)	9. Numbe derivative Securities Beneficia Owned Following Reported Transactie (Instr. 4)	e s I ully I	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership (Instr. 4)		
					Code	v	(A)	(D)	Date Exerc	cisable	Expiration Date	Title	Amou or Numb of Share	oer							
Stock Option (Right to Buy)	\$1.93	01/22/2020			M	<i>A</i> 1,600 (5) 01/05/		01/05/2021	Commo Stock	n 1,60	1,600		53,707		D						
Stock														T							

## **Explanation of Responses:**

\$1.93

- 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 \$24.995 to \$25.850. 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.

1.600

- 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 \$25.400 to \$25.780. 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:

Option

(Right to Buy)

/s/ Brian Keane, Attorney-in-

1,600

\$0.00

01/24/2020

52,107

D

<u>fact</u>

Common

01/05/2021

\*\* 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).

01/23/2020

\*\* 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.	