FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington,	D.C.	20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

l	OWR APPRO	VAL
	OMB Number:	3235-0287
l	Estimated average burd	en
l	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

1. Name and Address of Reporting Person * $\underline{Madaus\ Martin\ D}$						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, 113 HARTWELL AVENUE			11/	3. Date of Earliest Transaction (Month/Day/Year) 11/12/2019 4. If Amendment, Date of Original Filed (Month/Day/Year)									Officer (give title Other (specify below)						
(Street) LEXINGTON MA 02421				- 4. 1 -	t Amen	iament	i, Date	e of Orig	Jinai F	ilea (Montn/D	Individual or Joint/Group Filing (Check Applicable Line) X Form filed by One Reporting Person Form filed by More than One Reporting Person								
(City)	(S	tate)	(Zip)																
		Tab	le I - I	Non-Deri	vative	Sec	uritie	es A	cquir	ed, D	oisposed o	of, or E	Benefic	cially	Owned	1			
1. Title of Security (Instr. 3)			2. Transacti Date (Month/Day		Execut if any	A. Deemed Execution Date, f any Month/Day/Year)		3. Transaction Code (Instr. 8)			Acquired (A) or D) (Instr. 3, 4 and		Benefici Owned F		es ally Following	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Direct Indirect str. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
							. ,		Code	v	Amount	(A) or (D)	Price		Reported Transaction(s) (Instr. 3 and 4)				
Common Stock				11/12/20	019)			M		1,600	A	\$1.9	81.93		8,149		D	
Common Stock				11/12/2019		9			S ⁽¹⁾		1,600	D	\$22.27	771 ⁽²⁾ 6,5		549		D	
Common Stock			11/12/20	019				S ⁽¹⁾		2,300	D	\$22.23	322 ⁽³⁾	222	,666			See footnote ⁽⁴⁾	
Common Stock				11/13/2019					M		1,600	A	\$1.9	.93 8,3		149		D	
Common Stock 11/1:			11/13/20	019	19			S ⁽¹⁾ 1,		1,600	D	\$22.	22.04 6,5		549		D		
Common Stock				11/13/2019)			S ⁽¹⁾		2,300	D	\$21.78	315 ⁽⁵⁾	220,366				See footnote ⁽⁴⁾
		ī	able								sposed of s, converti				Owned				
1. Title of Derivative Security (Instr. 3)	2. 3. Transaction Date (Month/Day/Ye Price of Derivative Security		Execu if any	eemed Ition Date, h/Day/Year)		Transaction Code (Instr.		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		e Exer ation D h/Day/		7. Title and Amount of Securities Underlying Derivative Sect (Instr. 3 and 4)		S (I	. Price of Perivative Pecurity Instr. 5)	9. Numbe derivative Securities Beneficia Owned Following Reported Transacti (Instr. 4)	ve Ces F ally Cong (d tion(s)	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership (Instr. 4)
						v	(A)	(D)	Date Exerc	isable	Expiration Date	Title	Amou or Numl of Share	ber					
Stock Option (Right to Buy)	\$1.93	11/12/2019			M			1,600	((6)	01/05/2021	Commo Stock		00	\$0.00	69,70	7	D	
Stock Option (Right to Buy)	\$1.93	11/13/2019			M			1,600	(1	6)	01/05/2021	Commo Stock		00	\$0.00	68,10	7	D	
vnlanation	of Respons	2001																	

- 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 \$22.26 to \$22.34. 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.
- 3. This transaction was executed in multiple trades at prices ranging from \$21.98 to \$22.67. 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.
- 4. Held by a trust for which the Reporting Person is a trust advisor who shares voting and investment control.
- 5. This transaction was executed in multiple trades at prices ranging from \$21.39 to \$22.12. 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.
- 6. All options are vested and exercisable.

Remarks:

/s/ Brian Keane, Attorney-in-11/14/2019 **fact**

** 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.