SEC Form 4

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							

Instruction 1(b)	y continue. See I.			pursuant to Section 16(a) of the Securities Exchange Act of 1934									
1. Name and Addr Madaus Mat	1 0	Person [*]	or Section 30(h) of the Investment Company Act of 2. Issuer Name and Ticker or Trading Symbol <u>Quanterix Corp</u> [QTRX]	5. Relationsh (Check all ap	oplicable)	Reporting Person(s) to Issuer ole)							
					ector	10% Owner							
(Last) (First) (Middle) C/O QUANTERIX CORPORATION,			3. Date of Earliest Transaction (Month/Day/Year) 07/01/2020	belo	cer (give title ow)	Other (specify below)							
900 MIDDLES	SEX TURNPIF	KΈ	4. If Amendment, Date of Original Filed (Month/Day,	/Year) 6. Individual Line)	or Joint/Group Fil	ling (Check Applicable							
(Street)				X For	m filed by One Re	eporting Person							
BILLERICA	MA	01821			m filed by More th son	nan One Reporting							
(City)	(State)	(Zip)											

Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned

1. Title of Security (Instr. 3)	2. Transaction Date (Month/Day/Year)	2A. Deemed Execution Date, if any (Month/Day/Year)	Transaction Code (Instr.		4. Securities Disposed Of 5)			5. Amount of Securities Beneficially Owned Following Reported	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
			Code	v	Amount	(A) or (D)	Price	Transaction(s) (Instr. 3 and 4)		(1150. 4)
Common Stock	07/01/2020		A		651 ⁽¹⁾	Α	\$27.8 ⁽²⁾	11,196 ⁽³⁾	D	
Common Stock								164,966	Ι	See Footnote ⁽⁴⁾

Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities)

1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	3A. Deemed Execution Date, if any (Month/Day/Year)	4. Transa Code (8)		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		6. Date Exerc Expiration Da (Month/Day/Y	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. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)	
				Code	v	(A)	(D)	Date Exercisable	Expiration Date	Title	Amount or Number of Shares				

Explanation of Responses:

1. Consists of stock granted in lieu of cash fees for service on the Company's Board of Directors and committees thereof for the second quarter of 2020, pursuant to an election made in December 2019. 2. Closing price of the Company's common stock on the Nasdaq Global Market on July 1, 2020.

2. Closing price of the Company's common stock

3. Includes 2,270 restricted stock units.

4. Held by a trust for which the Reporting Person is a trust advisor who shares voting and investment control.

Remarks:

<u>/s/ Brian Keane, Attorney-in-</u> Fact 07/06/2020

** 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.

Check this box if no longer subject to Section 16. Form 4 or Form 5 obligations may continue. See