SEC Form 4	
------------	--

П

FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

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

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

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

OMB APPROVAL OMB Number: 3235-0287 Estimated average burden

Estimated average burden	
hours per response:	0.5

1. Name and Address of Reporting Person*			2. Issuer Name and Ticker or Trading Symbol <u>Quanterix Corp</u> [QTRX]	5. Relationship of Reporting Person(s) to Issuer (Check all applicable)					
<u>Madaus Martin D</u>				X	Director	10% Owner			
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 113 HARTWELL AVENUE		()	3. Date of Earliest Transaction (Month/Day/Year) 08/07/2018		Officer (give title below)	Other (specify below)			
			4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Indiv Line)	Individual or Joint/Group Filing (Check Applicabl				
(Street)				X	Form filed by One Re	porting Person			
LEXINGTON	MA	02421			Form filed by More the Person	an One Reporting			
(City)	(State)	(Zip)							

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

Table 1 Hon Denvarie Geournies Adquireu, Disposed oi, or Benenolary 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.					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	08/07/2018		Α		2,270 ⁽¹⁾	Α	\$ <mark>0</mark>	2,270 ⁽²⁾	D		
Common Stock								224,966	Ι	See footnote ⁽³⁾	

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)		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. 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				
Stock Option (Right to Buy)	\$14.8	08/07/2018		A		7,900		12/31/2018	08/07/2028	Common Stock	7,900	\$0	7,900	D	

Explanation of Responses:

1. Represents restricted stock units that vest as to 100% of the shares on December 31, 2018. Each restricted stock unit represents the right to receive one share of the Issuer's common stock.

2. Represents 2,270 restricted stock units.

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

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

** Signature of Reporting Person

Date

08/09/2018

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.