FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

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

OMB APPRO	VAL
OMB Number:	3235-0287
Estimated average burde	en
hours per response:	0.5
	OMB Number: Estimated average burde

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* Madaus Martin D (Last) (First) (Middle)							2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX] 3. Date of Earliest Transaction (Month/Day/Year)								k all app Dired	olicable) ctor er (give title		erson(s) to Is 10% C Other below	Owner (specify	
C/O QUANTERIX CORPORATION,						07/01/2019														
113 HARTWELL AVENUE							4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Individual or Joint/Group Filing (Check Applicable					
(Street) LEXINGTON MA 02421														Line) X Form filed by One Reporting Person Form filed by More than One Reporting Person						
(City)	(S	ate) (Zip)																	
	Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																			
1. Title of Security (Instr. 3) 2. Transacti Date (Month/Day						Execution (Year) if any		Deemed ecution Date, ny onth/Day/Year)		3. 4. Securitie Transaction Code (Instr. 8)				and 5) Securi Benefi		cially Following	Forn (D) c	n: Direct or Indirect nstr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
									Code	v	Amount	(A) or (D)	Price	Trai		ction(s) 3 and 4)			(111311. 4)	
Common Stock 07/0				07/01/	2019				A		476(1)	A	\$35.	38 ⁽²⁾	5,737(3)			D		
Common Stock															22	4,966		T I	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)	tive Conversion Date Execution Date, or Exercise (Month/Day/Year) if any					Transaction Code (Instr.		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		6. Date Exercisable and Expiration Date (Month/Day/Year)			and t of ies ying ive y (Instr. 3	Dei Sec (Ins	Price of rivative curity str. 5)		Owners Form: Direct (or Indir (I) (Inst	Ownership	11. Nature of Indirect Beneficial Ownership (Instr. 4)	
					Code	v	(A)	(D)	Date Exercis	sable	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 2019, pursuant to an election made in December 2018.
- 2. Closing price of the Company's common stock on the Nasdaq Global Market on July 1, 2019.
- 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:

/s/ Brian Keane, Attorney-in-07/03/2019 **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.