Instruction 1(b).

FORM 4

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

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	e: 0.5									

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				2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]										ck all app Direc	licable) tor	r 10% Own		Owner	
1		CORPORATION	∕iiddle) √,		3. Date of Earliest Transaction (Month/Day/Year) 01/04/2021								Officer (give title below) Other (specify below)						
900 MIDDLESEX TURNPIKE (Street) BILLERICA MA 01821				4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Individual or Joint/Group Filing (Check Applicable Line) X Form filed by One Reporting Person Form filed by More than One Reporting						
(City)	(St	ate) (Z	Zip)		Person														
	Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																		
1. Title of Security (Instr. 3) 2. Transactio Date (Month/Day/Y			/ear) I	2A. Deemed Execution Date, if any (Month/Day/Year)		е,	3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 a			nd 5) Securitie Beneficia		es Form ally (D) of Following (I) (II		Direct Indirect str. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)		
								Code V		Amount	(A) or (D)	Price		Transaction(s) (Instr. 3 and 4)				(301 4)	
Common Stock 04/01/202			!1			A		294(1)	A	\$61.45	15 ⁽²⁾ 14,6		690 ⁽³⁾		D				
Common Stock												164,966		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)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	3A. Deemed Execution Date, if any (Month/Day/Year)	4. Trans Code 8)	action (Instr.	5. Nu of Deriv Secu Acqu (A) of Dispo of (D) (Instr	rities ired r osed) r. 3, 4	Exp	iration	a Date Amo sy/Year) Sect Und Deri Sect		7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		Price of erivative ecurity istr. 5)	9. Numbo derivativ Securitie Beneficia Owned Followin Reported Transact (Instr. 4)	e s ally g	10. Ownershi Form: Direct (D) or Indirec (I) (Instr.	Beneficial Ownership t (Instr. 4)	
				Code	v	(A)	Date Exercise			Expiration e 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 first quarter of 2021.
- 2. Closing price of the Company's common stock on the Nasdaq Global Market on April 1, 2021.
- 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-

04/02/2021

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.