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							
haiina man naamanaa	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

Name and Address of Reporting Person* Madaus Martin D						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]								elationship o eck all applic X Director	able) r	g Pers	10% Ov	vner
`	ANTERIX	irst) CORPORATIOI FURNPIKE	(Middle)		01	3. Date of Earliest Transaction (Month/Day/Year) 01/03/2023								below)	(give title		Other (s below)	
(Street) BILLER (City)		IA State)	01821 (Zip)		_ 4. _	4. If Amendment, Date of Original Filed (Month/Day/Year)								dividual or Joint/Group Filing (Check Applicable Form filed by One Reporting Person Form filed by More than One Reporting Person				
		Ta	ble I - Noi	n-Der	ivativ	ve Se	ecurities	Acq	uired,	Dis	oosed of	, or Ber	eficiall	y Owned				
1. Title of Security (Instr. 3) 2. Trans. Date (Month/I				2A. Deemed Execution Date, if any (Month/Day/Year)		Transaction Dis			es Acquire Of (D) (Inst		5. Amount of		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	Transacti	Transaction(s) (Instr. 3 and 4)			(111501.4)
Common	ommon Stock 01/03			03/20	/2023		А		8,083 ⁽¹⁾ A		\$0.00	26,5	26,526 ⁽²⁾		D			
Common	Common Stock												254	254,633			Held by trust.	
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)	ate Execution Da	ransaction Code (Instr.		Derivative		6. Date Exercisable and Expiration Date (Month/Day/Year)			7. Title and Amor of Securities Underlying Derivative Secur (Instr. 3 and 4)		8. Price of Derivative Security (Instr. 5)	9. Number derivative Securities Beneficial Owned Following Reported	e s ally g	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership ct (Instr. 4)	
					Code	v	(A)	(D)	Date Exercisa		Expiration Date	Title	Amount or Number of Shares		Transacti (Instr. 4)	on(s)		
Stock Option (Right to Buy)	\$12.86	01/03/2023			A		12,124 ⁽³⁾		12/31/20)23	01/03/2033	Common Stock	12,124	\$0.00	12,12	24	D	

Explanation of Responses:

- 1. Represents restricted stock units that vest as to 100% of the shares on December 31, 2023 granted pursuant to the Quanterix Corporation Amended and Restated 2018 Non-Employee Director Compensation Policy. Each restricted stock unit represents the right to receive one share of the Issuer's common stock.
- 2. Includes 8,083 restricted stock units.
- 3. Represents non-qualified stock option to purchase 12,124 shares of common stock granted pursuant to the Quanterix Corporation Amended and Restated 2018 Non-Employee Director Compensation Policy.

Remarks:

/s/ Brian Keane, Attorney-in-01/04/2023

** 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.