FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPI	OMB APPROVAL									
OMB Number: 3235-02										
Estimated average burden										
hours per response:	0.5									

Section 16. Form 4 or Form 5 obligations may continue. See Instruction 1(b)

Check this box if no longer subject to

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]									k all applic	10		. ,	Ssuer Owner	
1	ANTERIX	CORPORATION	(Middle)			3. Date of Earliest Transaction (Month/Day/Year) 01/03/2022								Officer (give title Other (spec below) below)						
900 MIDDLESEX TURNPIKE (Street)					- 4.1	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						
BILLER:	ICA M	IA .	01821		_										Form filed by More than One Reporting Person					
(City)	(S	tate)	(Zip)																	
		Tal	ole I - N	on-Deri	ivativ	e Se	curities	Ac	quire	d, Di	sposed o	f, or Be	nefici	ally	Owned					
1. Title of Security (Instr. 3) 2. Transacti Date (Month/Day					Execution Date,			3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 a			nd 5) Securities Beneficial Owned Fo		ly	Form:	nership Direct Indirect tr. 4)	7. Nature of Indirect Beneficial Ownership		
						Code	v	Amount	(A) or (D)	Price		Reported Transactio (Instr. 3 an				(Instr. 4)				
Common Stock 01/03/20						022			A		2,668(1)	A	\$0.0	\$0.00)16 ⁽²⁾		D		
Common Stock 01/03/20					/2022)22			A		427(3)	A	\$42.4	42.41 ⁽⁴⁾ 18,4		43(2)		D		
Common Stock															164,966		I		See Footnote ⁽⁵⁾	
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. Deen Executio if any (Month/D	n Date,	4. Transa Code (i 8)				6. Date Exerc Expiration Da (Month/Day/Y		ate Amoun Year) Securit Underly		nt of ties ying tive Security		8. Price of Derivative Security (Instr. 5)	9. Numb derivativ Securiti Benefici Owned Followir Reporte Transac	ve es ially ng	10. Ownersh Form: Direct (D or Indire (I) (Instr.	Beneficial Ownership ct (Instr. 4)	
					Code	v	(A)	(D)	Date Exerci	sable	Expiration Date	Title	Amou or Numb of Share	er		(Instr. 4)				
Stock Option (Right to Buy)	\$42.41	01/03/2022			A		4,002 ⁽⁶⁾		12/31/	/2022	01/03/2032	Common Stock	4,00	02 \$0.00 4,		4,00	02	D		

Explanation of Responses:

- 1. Represents restricted stock units that vest as to 100% of the shares on December 31, 2022 granted pursuant to the Quanterix Corporation 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 2,668 restricted stock units.
- 3. Consists of stock granted in lieu of cash fees for service on the Company's Board of Directors and committees thereof for the fourth quarter of 2021.
- 4. Closing price of the Company's common stock on the Nasdaq Global Market on January 3, 2022.
- $5. \ Held \ by \ a \ trust \ for \ which \ the \ Reporting \ Person \ is \ a \ trust \ advisor \ who \ shares \ voting \ and \ investment \ control.$
- 6. Represents non-qualified stock option to purchase 4,002 shares of common stock granted pursuant to the Quanterix Corporation 2018 Non-Employee Director Compensation Policy.

Remarks:

/s/ Brian Keane, Attorney-in-

01/05/2022

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.