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											
hours per response	. 05										

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

1. Name and Address of Reporting Person* MEISTER PAUL M					2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]									Relationship of Reporting Person(s) to Issuer (Check all applicable) N Director 10% Owner					
		CORPORATION	Middle)		3. Date of Earliest Transaction (Month/Day/Year) 04/01/2022								Λ		cer (give title		Other (specify below)		
900 MIDDLESEX TURNPIKE						4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Individual or Joint/Group Filing (Check Applicable Line)					
(Street) BILLER	ICA M	A 0	1821											X	Form filed by One Reporting Person Form filed by More than One Reporting Person				
(City)	(St	ate) (Z	Zip)																
		Table	I - No	on-Deriva	tive S	Secui	rities	Acc	quirec	d, Dis	posed of	, or B	enefic	ially	Own	ed			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/N					·	Execution Date,		3. Transaction Code (Instr. 8) 4. Securities Acq Disposed Of (D) (5. Amount of Securities Beneficially Owned Follo Reported		ties cially I Following	Form (D) o	vnership n: Direct r Indirect nstr. 4)	7. Nature of Indirect Beneficial Ownership	
								Code	v	Amount	(A) or (D)	Price		Transa	nsaction(s) str. 3 and 4)			(Instr. 4)	
Common	Stock			04/01/20	022		Α		492(1)	A	\$30.4	30.47(2)		17,414 ⁽³⁾		D			
	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			insaction of Deri Sec Acq (A) Dispersion of (Instr. Sec Acq (A) Dispersion of (I (Instruction and		osed) r. 3, 4	Expiration I (Month/Day		ate	7. Title and Amount of Securities Underlying Derivative Security (Ins 3 and 4)		t				10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)		

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 2022.
- 2. Closing price of the Company's common stock on the Nasdaq Global Market on April 1, 2022.
- 3. Includes 2,668 restricted stock units.

Remarks:

/s/ Brian Keane, Attorney-in-

fact

04/05/2022

** 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.