FORM 4

to Section 16. Form 4 or Form 5 obligations may continue. See

Instruction 1(b).

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington,	D.C.	20549
-------------	------	-------

Check this box if no longer subject	STATEMENT OF CHANGES IN BENEFICIAL	OWNERSHIP
to Section 16. Form 4 or Form 5		
obligations may continue. Con		

OMB APPROVAL 3235-0287 Estimated average burden hours per response: 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* MEISTER PAUL M					2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]									5. Relationship of Reporting Person(s) to Issuer (Check all applicable)						
MEISTER PAUL W					-\-					-				X	Direc	tor		10% O	wner	
		CORPORATION	Middle)		3. Date of Earliest Transaction (Month/Day/Year) 01/04/2021										Office below	cer (give title ow)		Other (below)	specify	
900 MIDDLESEX TURNPIKE					4. If Amendment, Date of Original Filed (Month/Day/Year)									6. Individual or Joint/Group Filing (Check Applicable						
(Street)	ICA 14		1001												X Form filed by One Reporting Person					
BILLER 	ICA M	A 0	1821											Form filed by More than One Reporting Person						
(City)	(St	ate) (Z	Zip)																	
		Table	I - No	on-Deriva	tive S	Secui	rities	Acc	quired	d, Dis	sposed of	, or E	Benefic	ially	Own	ed				
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/					Execution Date,		ate,	3. Transaction Code (Instr. 8) 4. Securities Acquire Disposed Of (D) (Inst						5. Amount of Securities Beneficially Owned Following Reported		wnership n: Direct or Indirect nstr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)			
								Code	v	Amount	(A) or (D)	Price		Transa	nnsaction(s) str. 3 and 4)			(111501.4)		
Common Stock 04/01/20					021			A		223(1)	A	\$61.4	61.45(2)		13,430(3)		D			
		Tal	ble II						,		osed of, convertib			•	Owne	d				
1. Title of Derivative Security (Instr. 3)	rivative conversion or Exercise Price of Derivative Security Date (Month/Day/Year) Execution Date, if any (Month/Day/Year)			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) Date Expiration Exercisable Date			7. Title and Amount of Securities Underlying Derivative Security (Instr 3 and 4) Amoun or Numbe of Title Shares		_		9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s) (Instr. 4)		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 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.

Remarks:

/s/ Brian Keane, Attorney-in-

<u>fact</u>

04/02/2021

** 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.