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, I	D.C.	20549
---------------	------	-------

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL							
OMB Number: 3235-0287							
Estimated average burden							
hours per response:	0.5						

Instruc	tion 1(b).			Filed	pursua or Se	ant to S ection 3	ection 80(h) o	16(a) f the I	of the s	Securi ent Co	ties Exchang Impany Act o	e Act of f 1940	1934			, , , , , , , , , , , , , , , , , , ,	эропас.	
1. Name and Address of Reporting Person* MEISTER PAUL M				2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]								Relationshi Check all app X Direc	,	ng Per	rson(s) to Is			
1	QUANTERIX CORPORATION,					3. Date of Earliest Transaction (Month/Day/Year) 04/01/2020								Officer (give title Other (s below) below)				specify
900 MIDDLESEX TURNPIKE					4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Individual or Joint/Group Filing (Check Applicable Line)					
(Street) BILLER (City)			1821 Zip)											X Form	Form filed by One Reporting Person Form filed by More than One Reporting Person			
		Table	I - No	n-Deriva	tive \$	Secui	rities	Acc	uired	l, Dis	sposed of	, or B	enefici	ally Own	ed			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day)				Execution Dat				3. Transaction Code (Instr. 8) 4. Securities Acq Disposed Of (D)					d 5) Securi Benefi	icially d Following	Form (D) o	n: Direct or Indirect nstr. 4)	7. Nature of Indirect Beneficial Ownership	
									Code	v	Amount	(A) or (D)	Price	Transa	action(s) 3 and 4)			(Instr. 4)
Common Stock 04/01/2				04/01/20	020			Α		715(1)	A	\$19.2	1 ⁽²⁾ 9	,736 ⁽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)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	if any	emed tion Date, n/Day/Year)	4. Transaction Code (Instr. 8) Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)			Expira (Monti	ation D h/Day/	or		nt of ties ying tive ty (Instr. 1) Amount or Number of	8. Price of Derivative Security (Instr. 5)	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 2020, pursuant to an election made in December 2019.
- 2. Closing price of the Company's common stock on the Nasdaq Global Market on April 1, 2020.
- 3. Includes 2,270 restricted stock units.

Remarks:

/s/ Brian Keane, Attorney-in-

<u>fact</u>

04/03/2020

** 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.