FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

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

Check this box if no longer subject to Section 16. Form 4 or Form 5 obligations may continue. See Instruction 1(b).

	tion 1(b).	. de .		Filed	pursua or Se	nt to Section 3	ection 80(h) o	16(a) f the Ir	of the S ovestme	ecurit nt Co	ies Exchang mpany Act o	e Act o f 1940	f 1934			nours	perre	sponse:	0.5
1. Name and Address of Reporting Person* MEISTER PAUL M				2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]									ationship k all app Direc	,	ng Per	rson(s) to Is			
(Last) (First) (Middle) C/O QUANTERIX CORPORATION,					3. Date of Earliest Transaction (Month/Day/Year) 10/01/2020									Office below	er (give title v)		Other (s 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	ICA M	A 0	1821											X		filed by One filed by Mo		J	
(City)	(St	ate) (Z	<u>Z</u> ip)																
		Table	I - No	n-Deriva	tive S	Secui	rities	Acq	uired,	Dis	posed of	, or B	enefic	cially	Own	ed			
1. Title of Security (Instr. 3) 2. Transact Date (Month/Day					Exec		Deemed cution Date, y nth/Day/Year)				es Acquired (A Of (D) (Instr. 3,			5. Amount of Securities Beneficially Owned Following		6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)		7. Nature of Indirect Beneficial Ownership	
									Code	v	Amount	(A) (D)	Pric	е	Report Transa (Instr.	ed ction(s) 3 and 4)			(Instr. 4)
Common Stock 10/01/2					2020		A		396(1)	A	\$34	.7 ⁽²⁾	10	0,626(3)		D			
		Tal									osed of, convertib				Owne	d			
1. Title of Derivative Security (Instr. 3)	2. Onversion or Exercise Price of Derivative Security 3. Transaction Date (Month/Day/Year) (Month/Day/Year) 3A. Deemed Execution Date, if any (Month/Day/Year)				ransaction of ode (Instr. Derivative			6. Date Exercisable and Expiration Date (Month/Day/Year)			7. Title and Amount of Securities Underlying Derivative Security (Inst 3 and 4)		nt		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)	
				Code	v	(A)	(D)	Date Exercis	able	Expiration Date	Title	of Shares							

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 third 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 October 1, 2020.
- 3. Includes 2,270 restricted stock units.

Remarks:

/s/ Brian Keane, Attorney-in-

fact

10/05/2020

** 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.