FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

	OMB APPRO	VAL							
	OMB Number:	3235-0287							
	Estimated average burden								
l	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).

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]									ck all app	olicable) ctor		Owner	
(Last) (First) (Middle) C/O QUANTERIX CORPORATION,						3. Date of Earliest Transaction (Month/Day/Year) 10/01/2019									Offic belo	er (give title w)	Oth belo	er (specify w)	
113 HARTWELL AVENUE							4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Individual or Joint/Group Filing (Check Applicable Line)				
(Street) LEXINGTON MA 02421															X Form filed by One Reporting Person Form filed by More than One Reporting Person				
(City)	(St	ate) (Zip)																
Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned															_				
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day)						Execution Date,						es Acquired (A) o of (D) (Instr. 3, 4 a			Secur Benef	icially d Following	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	of Indirect	
						Code	v	Amount	(A) ((D)	Pr	ice	Trans	action(s) 3 and 4)		(,				
Common	Stock	2019	019		A		642 ⁽¹⁾ A		\$	21.41(2	6	,157 ⁽³⁾	D						
		Та	ble II -								osed of, convertib				Owned				
1. Title of Derivative Security (Instr. 3)	ve Conversion Date Execution Date, or Exercise (Month/Day/Year) if any		on Date, Day/Year)			5. Nu of Deriv Secu Acqu (A) of Dispo of (D) (Instrand 5	rative rities ired r osed)	6. Date Exerc Expiration Da (Month/Day/\) Date Exercisable		te	7. Title and Amount of Securities Underlying Derivative Security (Instr. and 4) Amount or Number of Title Shares		De Se (In	Price of rivative curity str. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	Ownershi Form: Direct (D) or Indirec (I) (Instr. 4	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 third quarter of 2019, pursuant to an election made in December 2018.
- 2. Closing price of the Company's common stock on the Nasdaq Global Market on October 1, 2019.
- 3. Includes 2,270 restricted stock units.

Remarks:

/s/ Brian Keane, Attorney-in-

fact

** Signature of Reporting Person Date

10/03/2019

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.