SEC Form 4	
FORM	4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

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

Check this box to indicate that a transaction was made pursuant to a

contract, instruction or written plan for the purchase or sale of equity securities of the issuer that is

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934 or Section 30(h) of the Investment Company Act of 1940

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

	tisfy the affirmative tions of Rule 10b5- uction 10.					
1. Name and Add <u>MEISTER</u>	ress of Reporting F PAUL M	² erson [*]	2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]		tionship of Reporting Po all applicable) Director	10% Owner
	(First) ERIX CORPOR		3. Date of Earliest Transaction (Month/Day/Year) 10/01/2024	Officer (give title Other (speci below) below)	Other (specify below)	
900 MIDDLE	SEX TURNPIK	E	4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Indiv Line)	vidual or Joint/Group Fili	ing (Check Applicable
(Street) BILLERICA	MA	01821		v	Form filed by One Re Form filed by More th Person	
(City)	(State)	(Zip)				

Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned

1. Title of Security (Instr. 3)	2. Transaction Date (Month/Day/Year)	2A. Deemed Execution Date, if any (Month/Day/Year)	Code (4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 and 5)		Securities Beneficially		7. Nature of Indirect Beneficial Ownership (Instr. 4)	
			Code	v	Amount	(A) or (D)	Price	Transaction(s) (Instr. 3 and 4)		(1130. 4)
Common Stock	10/01/2024		Α		1,200(1)	Α	\$12.5 ⁽²⁾	295,729 ⁽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)	3A. Deemed Execution Date, if any (Month/Day/Year)	4. Transa Code (8)		of Deriv Secu Acqu (A) o Dispe of (D	r osed) r. 3, 4	6. Date Exerc Expiration Da (Month/Day/Y	ate	Deriv	unt of rities rlying ative rity (Instr.	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)
				Code	v	(A)	(D)	Date Exercisable	Expiration Date	Title	Amount or Number 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 2024.

2. Closing price of the Company's common stock on the Nasdaq Global Market on October 1, 2024.

3. Includes 2,947 restricted stock units.

Remarks:

/s/ Brian Keane, Attorney-in-

10/02/2024

** Signature of Reporting Person Date

fact

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.