FORM 4

Check this box if no longer subject

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	

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

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]										all app Direc	tor	ng Pers	10% O	wner
		CORPORATION	Middle N,	e)		3. Date of Earliest Transaction (Month/Day/Year) 08/10/2022									Office below	er (give title		Other (below)	specify
900 MIDDLESEX TURNPIKE				4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Individual or Joint/Group Filing (Check Applicable Line)					pplicable	
(Street) BILLER	ICA M.	A 0	1821										X	Form filed by One Reporting Person Form filed by More than One Reporting Person					
(City)	(St	ate) (Ž	Zip)																
	Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																		
Date			2. Transaction Date (Month/Day/Y	Execution (ear) if any		med on Date, Day/Year)		3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 a			and 5) Sec Ber Ow		curities neficially		Direct	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
									Code	v	Amount	(A) or (D)	Price		Transa	action(s) 3 and 4)			(111511. 4)
Common Stock 08/10/20				08/10/202	22				P		186,347	A	\$8.50	85(1) 204)4,654 ⁽²⁾		D	
Common Stock 08/10/202			22				P		47,957	A	\$8.60	\$8.6073(3)		252,611(2)		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)	erivative Conversion Date Execution Date, or Exercise (Month/Day/Year)			ransaction Code (Instr.)		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		Expiration Date (Month/Day/Year)		7. Title and Amount of Securitles Underlying Derivative Security (Inst 3 and 4)		Der Sec (Ins	vative (urity (str. 5) I	9. Number of derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	y D (I	0. ownership orm: birect (D) r Indirect) (Instr. 4)	Beneficial Ownership t (Instr. 4)		
Fundamentia.					Code	v	(A)	(D)	Date Exer	e rcisable	Expiration Date	Title	Amount or Number of Shares						

- 1. This transaction was executed in multiple trades at prices ranging from \$7.58 to \$8.57. The price reported reflects the weighted average sale price. The reporting person undertakes to provide upon request to the SEC staff, the issuer or any security holder of the issuer full information regarding the number of shares and prices at which the transaction was effected.
- 3. This transaction was executed in multiple trades at prices ranging from \$8.58 to \$8.75. The price reported reflects the weighted average sale price. The reporting person undertakes to provide upon request to the SEC staff, the issuer or any security holder of the issuer full information regarding the number of shares and prices at which the transaction was effected.

Remarks:

/s/ Brian Keane, Attorney-in-

08/10/2022

fact

** 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.