EC Form 4						
FOI	RM 4	UNITED ST	TATES SECURITIES AND EXCHANGE CO Washington, D.C. 20549	OMMISSION	OME	3 APPROVAL
Section 16. Fo	y continue. See		Filed pursuant to Section 16(a) of the Securities Exchange Act of 19 or Section 30(h) of the Investment Company Act of 1940	OMB Number Estimated at hours per re	verage burden	
1. Name and Address of Reporting Person [*] MEISTER PAUL M			2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]	(Check all appli X Directo Officer	or (give title	10% Owner Other (specify
	(First) RIX CORPORA EX TURNPIKI	1. A	3. Date of Earliest Transaction (Month/Day/Year) 01/04/2021	below)		below)
(Street)			4. If Amendment, Date of Original Filed (Month/Day/Year)	Line)	Joint/Group Filing filed by One Repo	g (Check Applicable orting Person
BILLERICA (City)	MA (State)	01821 (Zip)	_	Form Perso	filed by More thar n	n One Reporting

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 (Instr.		4. Securities Acquir Disposed Of (D) (Ins			5. Amount of Securities Beneficially Owned Following Reported	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)		
			Code	v	Amount	(A) or (D)	Price	Transaction(s) (Instr. 3 and 4)		(1150.4)		
Common Stock	01/04/2021		A		2,270 ⁽¹⁾	A	\$0.00	12,896(2)	D			
Common Stock	01/04/2021		A		311 ⁽³⁾	A	\$44.11(4)	13,207 ⁽²⁾	D			

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)	ction Derivative		6. Date Exerc Expiration Da (Month/Day/Y	ate	7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		8. Price of Derivative Security (Instr. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s)	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		(Instr. 4)		
Stock Option (Right to Buy)	\$44.1	01/04/2021		A		7,900 ⁽⁵⁾		12/31/2021	01/04/2031	Common Stock	7,900	\$0.00	7,900	D	

Explanation of Responses:

1. Represents restricted stock units that vest as to 100% of the shares on December 31, 2021 granted pursuant to the Quanterix Corporation 2018 Non-Employee Director Compensation Policy. Each restricted stock unit represents the right to receive one share of the Issuer's common stock.

2. Includes 2,270 restricted stock units.

3. Consists of stock granted in lieu of cash fees for service on the Company's Board of Directors and committees thereof for the fourth quarter of 2020.

4. Closing price of the Company's common stock on the Nasdaq Global Market on January 4, 2021.

5. Represents non-qualified stock option to purchase 7,900 shares of common stock granted pursuant to the Quanterix Corporation 2018 Non-Employee Director Compensation Policy.

Remarks:

/s/ Brian Keane, Attorney-in-01/05/2021

** Signature of Reporting Person

Reminder: Report on a separate line for each class of securities beneficially owned directly or indirectly.

 \ast 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.

				Date (Month/Day/Year)			Execution Date, if any (Month/Day/Year)		Transaction Code (Instr. 8)		Disposed O	f (D) (Instr.	3, 4 and 5)	Securities Beneficially Owned Following Reported		Form: Direct (D) or Indirect (I) (Instr. 4)		Indir Bene Own (Inst	
									Code	v	Amount	(A) or (D)	Price	Transact (Instr. 3 a	ion(s)			(inst	
nmon Stock					4/2021	-			A		2,270 ⁽¹⁾	A	\$0.00	12,896 ⁽²⁾			D		
nmon Stock				01/04/2021					A		3 11 ⁽³⁾	A	\$44.11(4)) 13,207 ⁽²⁾			D		
	Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities)																		
tle of vative urity r. 3) 2. Conversion or Exercise Price of Derivative Security		3. Transaction Date (Month/Day/Year)	Date Execution Da			ction Instr.	5. Number Derivativ Securitie Acquired or Dispos of (D) (In 3, 4 and 9	6. Date Exercisa Expiration Date (Month/Day/Yea		ite	7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		3. Price of Derivative Security (Instr. 5)	9. Number derivative Securities Beneficiall Owned Following Reported	y	10. Ownershij Form: Direct (D) or Indirect (I) (Instr. 4	Be Ov (In		
					Code	v	(A)	(D)	Date Exercis	able	Expiration Date	Title	Amount or Number of Shares		Transactio (Instr. 4)	n(s)			
k on ht to	\$44.1	01/04/2021			Α		7,900 ⁽⁵⁾		12/31/2	.021	01/04/2031	Common Stock	7,900	\$0.00	7,900		D		

fact

Date