FORM 4

Check this box if no longer subject

UNITED STATES SECURITIES AND EXCHANGE COMMISSION Washington, D.C. 20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL

L								
OMB Number:	3235-0287							
Estimated average burden								
hours per response:	0.5							

to Section 16. F	Form 4 or Form					Estimate	d average bur	rden
Obligations may Instruction 1(b)		F	Filed pursuant to Section 16(a) of the Securities Exchange Act of 19	934		hours pe	er response:	0.5
(_)		1	or Section 30(h) of the Investment Company Act of 1940					
1. Name and Addr <u>Fry John J</u> (Last) C/O QUANTE 900 MIDDLES	(First) RIX CORPO	(Middle) DRATION,	2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX] 3. Date of Earliest Transaction (Month/Day/Year) 10/01/2020		ck all applicabl Director Officer (giv below)	le) ve title	10% (Other below	Owner (specify /)
			4. If Amendment, Date of Original Filed (Month/Day/Year)		lividual or Join	t/Group F	iling (Check	Applicable
(Street)				- I '	Form filed	by One F	Reporting Pe	rson
BILLERICA	ME	01821	_		 5. Relationship of Reporting Person(s. (Check all applicable) Director Officer (give title Officer (give title Officer (give title General Counsel & Sect 6. Individual or Joint/Group Filing (Checking) X Form filed by One Reporting Form filed by More than One Person 	1 0		
(City)	(State)	(Zip)						
		Table I - Non-Der	rivative Securities Acquired, Disposed of, or Ber	neficiall	y Owned			
	()							

1. Title of Security (Instr. 3)	2. Transaction Date (Month/Day/Year)	2A. Deemed Execution Date, if any (Month/Day/Year)	Code (ransaction Disposed Of (I Code (Instr.				5. Amount of Securities Beneficially Owned Following Reported	Form: Direct (D) or Indirect (I)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
			Code	v	Amount	(A) or (D)	Price	Transaction(s) (Instr. 3 and 4)	(Instr. 4)	(1150. 4)
Common Stock	10/01/2020		S		143(1)	D	\$34.1224(2)	31,227 ⁽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)			vative nities nired r osed) r. 3, 4	Expiration Date (Month/Day/Year)		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) (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. The sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 468 RSUs.

2. This transaction was executed in multiple trades at prices ranging from \$33.881 to \$34.51. 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. Includes 26,148 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorney-10/02/2020

in-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.