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

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

to Section 16.	Form 4 or Form y continue. See).	5	Filed pursuant to Section 16(a) of the Securities Exchange Act of 193 or Section 30(h) of the Investment Company Act of 1940	34		Estimated av hours per res	erage burden sponse:	0.5
1. Name and Address of Reporting Person [*] Fry John J (Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE		(Middle) DRATION,	2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX] 3. Date of Earliest Transaction (Month/Day/Year) 01/05/2021		ck all applicable Director Officer (giv below)	Officer (give title Other (sp		
(Street) BILLERICA (City)	MA (State)	01821 (Zip)	4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Indi Line) X	Form filed	by One Repo) (Check Applic orting Person n One Reportin	
		Table I - Non-Dei	rivative Securities Acquired, Disposed of, or Bend	eficially	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 (D Code (Instr.				Securities Beneficially Owned Following	Form: Direct (D) or Indirect (I)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
			Code	v	Amount	(A) or (D)	Price	Reported Transaction(s) (Instr. 3 and 4)	(Instr. 4)	(instr. 4)
Common Stock	01/05/2021		S		153(1)	D	\$44.1317 ⁽²⁾	31,885 ⁽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)	Code (4. 5. Number Transaction Code (Instr. Derivative		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 \$43.91 to \$44.62. 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 24,744 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorney-01/06/2021

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.