FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington D.C. 20549

vvasiiiigtoii,	D.C. 20349

STATEMENT	OF CHANGES	IN BENEFICIAL	OWNERSHIP

OMB APPROVAL										
OMB Number:	3235-0287									
Estimated average burden										
hours por rosponso:	0.5									

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

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

Name and Address of Reporting Person* Fry John J						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]								Check	all app	o of Reportir licable) tor er (give title	ng Pei	rson(s) to Is 10% O Other (wner
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						3. Date of Earliest Transaction (Month/Day/Year) 06/11/2020								X	belov	below) General Counsel &		below)	
(Street) BILLERICA ME 01821 (City) (State) (Zip)					4. If <i>i</i>	4. If Amendment, Date of Original Filed (Month/Day/Year)								i. Indiv ine) X					
		Table	I - I	Non-Deriva	tive	Secui	rities	Ac	quir	ed, D	isposed o	f, or I	3enefic	ially	Own	ed			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Yea			ar) E	2A. Deemed Execution Date, if any (Month/Day/Year)		Ī	3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 an			nd 5) Secu Bend		urities I eficially (ned Following I		wnership m: Direct or rect (I)	7. Nature of Indirect Beneficial Ownership (Instr. 4)		
								c	Code	v	Amount	(A) or (D)	Price		Transaction(s) (Instr. 3 and 4)		(Instr. 4)		(msu. 4)
Common Stock 06/11/202			06/11/2020	0			S		1,842(1)	D	\$24.7653(2)		31,803 ⁽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)	crivative conversion or Exercise (Month/Day/Year) Frice of Derivative Security Execution Date, if any (Month/Day/Year)		4. Transa Code 8)		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		Expiration Date (Month/Day/Year)			Amo Secu Unde Deriv Secu 3 and	Amount or Number of	nt er				10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)		

Explanation of Responses:

- 1. The sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 5,625 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$24.7502 to \$24.81. 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 28,020 restricted stock units.

Remarks:

<u>/s/ Brian Keane, as Attorney-in-Fact</u>

06/12/2020

** 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.