FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington,	D.C.	20549
-------------	------	-------

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

## 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*  Fry John J						2. Issuer Name <b>and</b> Ticker or Trading Symbol Quanterix Corp [ QTRX ]										all app Direc		ng Pe	rson(s) to Is 10% Ov Other (s	wner
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						3. Date of Earliest Transaction (Month/Day/Year) 09/01/2021									Λ	Gen	,	sel 8	below) el & Secretary	
(Street) BILLERICA MA 01821 (City) (State) (Zip)						4. If Amendment, Date of Original Filed (Month/Day/Year)									6. Indiv Line) X					
Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																				
1. Title of Security (Instr. 3)  2. Transact Date (Month/Date					Executio ear) if any				3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 an				nd 5) Sed Ber Ow		curities neficially ned Following		m: Direct or rect (I)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
							Code V		Am	ount	(A) or (D)	Price		Transa	Transaction(s) (Instr. 3 and 4)			(3 4)		
Common	Common Stock 09/01/20							T	S	П	3	300(1)	D	\$49.78	B2 <sup>(2)</sup>	33,399(3)			D	
Common Stock 09/				09/01/202	21				S			50 <sup>(1)</sup>	D	\$50.74	85 <sup>(4)</sup>	<sup>4)</sup> 33,349 <sup>(3)</sup>			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)	Exe if an	Deemed cution Date, ry nth/Day/Year)		action (Instr.	5. Nu of Deriv Secu Acqu (A) o Dispo of (D (Instrand 5	vative irities iired r osed ) r. 3, 4	Exp (Mo	piration onth/Da	n Date		Amo Secu Unde Deriv Secu 3 and	le and unt of rities rrlying rative rity (Instr. 1 4) Amount or Number	Der Sec (Ins	Price of rivative curity str. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	y	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership t (Instr. 4)
					Code	ode V (A) (D						Date	Title	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 700 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$49.32 to \$50.22. 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 20,270 restricted stock units.
- 4. This transaction was executed in multiple trades at prices ranging from \$50.39 to \$51.21. 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, as Attorney-

09/03/2021

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.