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(h)

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

	OMB APPROVAL								
	OMB Number: 3235-0287								
	Estimated average burden								
ı	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

Name and Address of Reporting Person* Fry John J						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]									(Check	ionship of Reportir all applicable) Director Officer (give title		g Person(s) to Issi 10% Owr Other (sp		wner
(Last)		3. Date of Earliest Transaction (Month/Day/Year) 03/01/2022									X	below			below)					
900 MIDDLESEX TURNPIKE (Street) BILLERICA MA 01821					4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)									6. Indiv Line) X	-/				
(City)	(City) (State) (Zip) Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Ye				ear) i	2A. Deemed Execution Date,		3. Ti C	3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 ar			d (A) or	5. Am Secui Bener		nount of 6 rities F ficially (I		n: Direct r ect (I)	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
								c	ode	v	Amount		(A) or (D)	Price	Trans		action(s) 3 and 4)		nstr. 4)	(111341. 4)
Common Stock				03/01/202	22				S		272 ⁽¹)	D	\$31.20)66 ⁽²⁾	30,852 ⁽³⁾			D	
Common Stock 03/01/20				03/01/202	!2				S		99(1		D	\$32.15		30,753(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)	Exe if ar	Deemed cution Date, ny nth/Day/Year)		saction e (Instr.	5. Num of Derivat Securit Acquir (A) or Dispos of (D) (Instr. 3 and 5)	ive ies ed ed	Expiration Date (Month/Day/Year)			Amo Secu Unde Deriv	Amount	Deri Sec (Ins		9. Number derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	Owners Form: Direct (I or Indire (I) (Instr	Ownership	11. Nature of Indirect Beneficial Ownership (Instr. 4)	
					Code	e V	(A) (D)		Date Exercisal			ation	Title	or Number of Shares	r					

Explanation of Responses:

- 1. The sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 776 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$30.84 to \$31.83. 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 15,075 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorney-

in-Fact

03/03/2022

** 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.