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

Name and Address of Reporting Person* Fry John J						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]										all app Direc	ionship of Reporting all applicable) Director Officer (give title		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) 04/01/2021										Gen	,	sel &	below) el & Secretary	
(Street)	(Street) BILLERICA MA 01821					4. If Amendment, Date of Original Filed (Month/Day/Year)									6. Indiv Line) X	vidual or Joint/Group Filing (Check Applicable Form filed by One Reporting Person Form filed by More than One Reporting Person				on
(City)	(3)				<u> </u>															
1. Title of Security (Instr. 3) 2. Trai				2. Transaction	2A. Deen		med on Date,		3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 an			d (A) or	5. Amo Securit Benefic		ount of ties cially d Following	Forr (D) o Indii	n: Direct or rect (I)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
								Code	v	Amount		(A) or (D)	Price		Transaction(s) (Instr. 3 and 4)		(Instr. 4)		(111341.4)	
Common	Stock		04/01/202	21				S			60(1)	D	\$61.9	61.9196 ⁽²⁾		34,973(3)		D		
Common Stock 04/02					1	1					1	126 ⁽¹⁾	D	\$62.69	62.6984(4)		34,847(3)		D	
Common Stock 04/					1				S			15(1)	D	\$63.36		34,832(3)			D	
		Tal	ble I	II - Derivati (e.g., pu												Owne	d			
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, ıy nth/Day/Year)		ransaction of Code (Instr. Derivativ				piration	kercisable and n Date ay/Year)		Amo Secu Unde Deriv Secu	7. Title and Amount of Securities Underlying Derivative Security (Inst 3 and 4)		rice of ivative urity tr. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction(: (Instr. 4)	y	Ownership Form: Direct (D) or Indirect (I) (Instr. 4	Beneficial Ownership (Instr. 4)
					Code	Code V (A) (C			Date) Exercisab			Expiratior Date	n Title	Amoun or Numbe 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 700 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$61.30 to \$62.27. 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 23,770 restricted stock units.
- 4. This transaction was executed in multiple trades at prices ranging from \$62.30 to \$63.21. 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-

04/05/2021

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.