Ch

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 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]									heck	ionship of Reportin all applicable) Director Officer (give title		ng Person(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) 11/02/2021									X	below) General Counse		sel &	below)	`
(Street) BILLER (City)	ICA M.	A 0	1821 Zip)	L	4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)									Indiv ne) X	fividual or Joint/Group Filing (Check Applicable Form filed by One Reporting Person Form filed by More than One Reporting Person				
	Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																			
1. Title of Security (Instr. 3) 2. Transact Date				2. Transaction	2A. Deem		ned n Date,	3. Ti C	3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 an			(A) or	5. An Secu Bene Own		ount of ties cially d Following	Forn (D) o Indir	n: Direct or rect (I)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
								С	Code V		Amount	(A) or (D)	P	Price		Reported (In Transaction(s) (Instr. 3 and 4)		(IIISt	Instr. 4)	(111301.4)
Common Stock 11				11/02/202	21				S		239(1)	D	\$	\$52.3974	974 ⁽²⁾		32,770 ⁽³⁾		D	
Common Stock				11/02/202	21				S		92(1)	D	\$	\$53.1449	9(4)	32,678			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, ry nth/Day/Year)		saction e (Instr.	5. Num of Derivat Securit Acquin (A) or Dispos of (D) (Instr. 3 and 5)	ive ies ed ed	Exp	Date Exercisable and piration Date on the piration Date on the piration of the			7. Title and Amount of Securities Underlying Derivative Security (Ins 3 and 4)		Deri Sec		9. Number of derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	Owners Form: Direct (or Indir (I) (Insti	Ownership	11. Nature of Indirect Beneficial Ownership (Instr. 4)
					Code	de V (A) (D)				e rcisabl	Expiratio e Date	n Title		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 700 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$51.86 to \$52.85. 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 18,870 restricted stock units.
- 4. This transaction was executed in multiple trades at prices ranging from \$52.96 to \$53.78. 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-

11/04/2021

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.