FORM 4

Instruction 1(b).

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

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

Check this box if no longer subject	STATE
to Section 16. Form 4 or Form 5	
obligations may continue. See	

MENT 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]									Check	all app	o of Reportir licable) tor er (give title	ng Pe	erson(s) to Is 10% Ov Other (s	wner
· `		st) (M CORPORATION TURNPIKE	Middle <mark>N</mark> ,	e)		3. Date of Earliest Transaction (Month/Day/Year) 02/02/2022								X	belov	elow) General Couns		below)	·	
(Street) BILLER (City)	ICA M	A 0	182: Zip)			4. If Amendment, Date of Original Filed (Month/Day/Year)						Í	ine) X	Form Form Perso	fual or Joint/Group Filing (Check Applicable Form filed by One Reporting Person Form filed by More than One Reporting Person					
		Table	I - I	Non-Deriva	tive	Secui	rities	Ac	qui	red, D)isp	posed o	f, or I	Benefic	ially	Own	ed			
Date				2. Transaction Date (Month/Day/Ye	Execution (ear) if any				3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 an				nd 5) Secu Bene Own		icially d Following	Fori (D) (Indi	m: Direct or irect (I)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
									Code	v	Am	nount	(A) or (D)	or Price		Reported Transaction(s) (Instr. 3 and 4)		(Instr. 4)		(111501.4)
Common	Stock			02/02/202	2				S		2	272(1)	D	\$31.200	31.2066 ⁽²⁾ 31,704 ⁽³⁾ D					
Common	Stock			02/02/202	2				S			99(1)	D	\$32.1	15	31,605 ⁽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)	Derivative Conversion Date Security or Exercise (Month/Day/Year) Execution Date,				4. Transa Code 8)	action (Instr.	5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		Expiration Date (Month/Day/Yea			Date //Year) Amount of Securities Underlying Derivative Security (Inst 3 and 4)		unt of rities erlying rative rity (Instr. 1 4)	Deri Sec	rice of ivative urity tr. 5)	9. Number derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	,	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)
					Code	v	(A)	(D)	Da Ex	ate kercisab	le	Expiration Date	Title	Amount or 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 \$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 16,770 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorney-

in-Fact

02/04/2022

** 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.