FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

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

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

3235-0287 ES IN BENEFICIAL OWNERSHIP Estimated average burden hours per response: Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934 or Section 30(h) of the Investment Company Act of 1940

OMB APPROVAL

0.5

Name and Address of Reporting Person* Fry John J					2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]											all app	ionship of Reportir all applicable) Director Officer (give title		10% O	
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE				3. Date of Earliest Transaction (Month/Day/Year) 02/01/2021										Λ	below) below) General Counsel & Secretary				y	
(Street) BILLERICA MA 01821 (City) (State) (Zip)					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				
		Table	I - I	Non-Deriva	tive	Secur	ities	Ac	quir	ed, D	ispos	ed o	f, or E	Benefic	cially	Own	ed			
Date				2. Transaction Date (Month/Day/Ye	Execution (ear)				3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 ar				nd 5) Sec Ben Owi		urities neficially ned Following		wnership n: Direct or rect (I) tr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
				Code					v	Amoun	:	(A) or (D)	Price		Transa	eported (In ransaction(s) nstr. 3 and 4)		u. 4)	(IIISU. 4)	
Common Stock				02/01/202	21				S		68(1)	D	\$67.92	298(2)	31,817(3)			D	
Common Stock				02/01/202	21				S		95 ⁽¹)	D	\$68.49	925 ⁽⁴⁾	31,722(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	Exercise (Month/Day/Year) if any (Month/Day/Year) ivative				action (Instr.	5. Num of Derive Securi Acqui (A) or Dispo of (D) (Instr. and 5	ative rities ired sed	Expiration Date (Month/Day/Year)			7. Title and Amount of Securities Underlying Derivative Security (Inst 3 and 4)		Deri Sec (Ins	rice of ivative urity tr. 5)			10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)	

Explanation of Responses:

- 1. The sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 468 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$67.25 to \$68.23. 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.
- 4. This transaction was executed in multiple trades at prices ranging from \$68.27 to \$68.63. 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-

02/02/2021

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.