FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION Washington, D.C. 20549

OMB APPROVAL

OMB Number:	3235-0287								
Estimated average	Estimated average burden								
hours per response:	. 0.5								

		Iable I - Non-L	Derivative Securities Acquired, Disposed of, or Ber	(Check all applicable) Director 10% Owner Director 10% Owner X Officer (give title below) Other (specify below) General Counsel & Secretary 6. Individual or Joint/Group Filing (Check Applicable Line) X Form filed by One Reporting Person Form filed by More than One Reporting Person				
		Table I Non F	Derivative Securities Acquired Dispessed of an Per	acticially Owned				
(City)	(State)	(Zip)		Person				
(Street) BILLERICA	MA	01821		X Form filed by One Reporting Person				
			4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Individual or Joint/Group Filing (Check Applicable				
C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE								
(Last)	(Last) (First) (Middle)		3. Date of Earliest Transaction (Month/Day/Year) 08/01/2022	, , ,				
1. Name and Addre Fry John J	ess of Reportir	ng Person [*]	2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]	X Officer (give title Other (specify				
Instruction 1(b).			Filed pursuant to Section 16(a) of the Securities Exchange Act of 19 or Section 30(h) of the Investment Company Act of 1940	934 0.5				

1. Title of Security (Instr. 3)	2. Transaction Date (Month/Day/Year)	2A. Deemed Execution Date, if any (Month/Day/Year)	3. Transa Code (8)		4. Securities Disposed Of			Securities	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership
			Code	v	Amount	(A) or (D)	Price	Transaction(s) (Instr. 3 and 4)	(msu. 4)	(Instr. 4)
Common Stock	08/01/2022		S		401(1)	D	\$15.6147 ⁽²⁾	37,671 ⁽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)	3A. Deemed Execution Date, if any (Month/Day/Year)		saction e (Instr. (A) or Disposed of (D) (Instr. 3, 4 and 5)		6. Date Exercisable and Expiration Date (Month/Day/Year)		7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		8. Price of Derivative Security (Instr. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction(S) (Instr. 4)	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)	
				Code	v	(A)	(D)	Date Exercisable	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 776 RSUs.

2. This transaction was executed in multiple trades at prices ranging from \$15.58 to \$15.6412. 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 20,232 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorney-08/03/2022

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.

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

in-Fact ** Signature of Reporting Person

Check this box if no longer subject to Section 16. Form 4 or Form 5