Instruction 1(b).

FORM 4

Check this box if no longer subject

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

wasnington,	D.C. 20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

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

to Section 16. Form 4 or Form 5 obligations may continue. See 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	,	ng Pe	10% O	
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						3. Date of Earliest Transaction (Month/Day/Year) 08/03/2021								X	below		sel 8	below)	
(Street) BILLERICA MA 01821 (City) (State) (Zip)					4. If A	4. If Amendment, Date of Original Filed (Month/Day/Year)								. Indiv ine) X	·				
		Table	I - N	on-Deriva	tive S	Secui	rities	Ac	quire	d, Di	sposed of	f, or E	Benefic	ially	Own	ed			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Yell)					Year)	Execution I		on Date,		3. Transaction Code (Instr. 8) 4. Securities Disposed Of				and 5) Secu Bene		cially I Following	Forn (D) (wnership m: Direct or Indirect nstr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
										v	Amount	(A) or (D)	Price	Tra		action(s) . 3 and 4)			
Common Stock 08/03/202					21				S		344(1)	D	\$53.78	.788(2)		33,699(3)		D	
		Tal	ble II	- Derivati (e.g., pu							osed of, convertib				Owne	d			
1. Title of Derivative Security (Instr. 3)	ve Conversion Date Execution Date, or Exercise (Month/Day/Year) if any			Transaction Code (Instr.		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		6. Date Exercisable and Expiration Date (Month/Day/Year)			e and unt of rities rlying ative rity (Instr. 4)	Der Sec (Ins	rice of ivative urity tr. 5)	9. Number derivative Securities Beneficiall Owned Following Reported Transaction (Instr. 4)	у	Ownership Form: Direct (D) or Indirect (I) (Instr. 4	Beneficial Ownership (Instr. 4)		
					Code	v	(A)	(D)	Date Exer	cisable	Expiration Date	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 \$53.43 to \$53.99. 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 20,970 restricted stock units

Remarks:

/s/ Brian Keane, as Attorneyin-Fact

08/03/2021

** 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.