Instruction 1(b).

FORM 4

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

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

STATEMENT	OF	CHANGES	IN BEN	NEFICIAL	OWNERS	HIP

OMB APP	OMB APPROVAL										
OMB Number:	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	tionship of Repor all applicable) Director Officer (give titl		10%		Issuer Owner (specify	
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE				3. Date of Earliest Transaction (Month/Day/Year) 04/01/2022										X				below)		
(Street) BILLERICA MA 01821 (City) (State) (Zip)			4. If Amendment, Date of Original Filed (Month/Day/Year)									6. Indiv Line) X	, i							
		Table	1-1	Non-Deriva	tive	Secui	rities	Ac	quir	ed, D	ispose	d of	f, or E	Benefic	cially	Own	ed			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Yes				2A. Deemed Execution Date, if any (Month/Day/Year)		,	3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 ar				nd 5) Se Be Ov		5. Amount of Securities Beneficially Dwned Following Reported		wnership n: Direct or rect (I) tr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)		
						-	Code	v	Amount		(A) or (D)	Price		Transaction(s) (Instr. 3 and 4)		(11151	u. 4)	(111341. 4)		
Common Stock 04/01/20			04/01/202	22				S		524(1)		D	D \$29.7929 ⁽²⁾		30,229(3)			D		
		Tal	ble	II - Derivati (e.g., pu)wne	d			
1. Title of Derivative Security (Instr. 3)	Derivative Conversion Date Execution Date, Security or Exercise (Month/Day/Year) if any					action (Instr.			Expiration Date (Month/Day/Year)			7. Title and Amount of Securities Underlying Derivative Security (Insi 3 and 4)		Deri Sec (Ins	rice of vative derivative Securities (r. 5) Beneficiall Owned Following Reported Transactic (Instr. 4)		Ownershi Form: y Direct (D) or Indirec (I) (Instr. 4		Beneficial Ownership t (Instr. 4)	
					Code	v	(A)	(D)	Dat Exe	e ercisabl	Expira Date	ition	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 776 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$29.25 to \$30.00. 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 14,299 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorneyin-Fact

** Signature of Reporting Person Date

04/05/2022

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.