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

Washington, D.C. 20549

<b>STATEMENT</b>	OF CHANGES IN	N BENEFICIAL	<b>OWNERSHIP</b>
• .,	0. 0		•

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

1. Name and Address of Reporting Person*  Fry John J					_ <u>Q</u> ı	2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [ QTRX ]								eck all appli Direct Office	ationship of Reporting Per call applicable) Director Officer (give title below)		10% Ov Other (s below)	ner
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						3. Date of Earliest Transaction (Month/Day/Year) 02/18/2020									General Counsel & Secretary			
(Street) BILLER (City)			01821 (Zip)		_	f Ame	ndment, I	Date o	of Origina	l Filed	d (Month/Da	ay/Year)	6. I Lin	X Form	iled by One	e Repo	(Check Apporting Person  One Repor	1
		Tab	le I - Noi	n-Deri	vativ	e Se	curities	s Ac	quired	, Dis	posed o	of, or Be	neficial	ly Owned	l			
Date			saction /Day/Ye	action 2A. Deemed Execution Date, if any (Month/Day/Year)		Code	Transaction Disposed Of (I Code (Instr. 5)		ties Acquir I Of (D) (Ins	Acquired (A) or (D) (Instr. 3, 4 and		5. Amount of Securities Beneficially Owned Following Reported		: Direct r Indirect str. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)			
									v	Amount	(A) c (D)	Price	Transac	action(s) . 3 and 4)			(	
Common Stock 02/18			8/202	2020		A		11,143	11,143 <sup>(1)</sup> A \$		0 33,	33,645(2)		D				
		7	Гable II -								osed of, converti			Owned				
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)	Date,	4. Transaction Code (Instr. 8)		ı of E		Expiration	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)	e S Illy	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership (Instr. 4)
					Code	v	(A)	(D)	Date Exercisa		Expiration Date	Title	Amount or Number of Shares					
Stock Option (right to buy)	\$27.76	02/18/2020			A		26,000		(3)		02/18/2030	Common Stock	26,000	\$0.00	26,000	0	D	

## **Explanation of Responses:**

- 1. Represents shares of stock issuable on vesting of restricted stock units. The restricted stock units vest as to 25% of the shares on February 18, 2021 and vest as to an additional 2.083% of the shares per month for 36 months thereafter.
- 2. Includes 33,645 restricted stock units.
- 3. The option vests as to 25% of the shares on February 18, 2021 and vests as to an additional 2.083% of the shares per month for 36 months thereafter.

## Remarks:

/s/ John J Fry

02/20/2020

\*\* Signature of Reporting Person

Date

Reminder: Report on a separate line for each class of securities beneficially owned directly or indirectly.

- $^{\star}$  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.