П

FORM 4

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

OMB APPROVAL

Check this box if no longer subject to	0
Section 16. Form 4 or Form 5	
obligations may continue. See	
Instruction 1(b).	

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

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 Number: 3235-028									
Estimated average burden									
hours per response:	0.5								

		n*	2. Issuer Name and Ticker or Trading Symbol <u>Quanterix Corp</u> [QTRX]		tionship of Reporting Pers all applicable) Director Officer (give title	rson(s) to Issuer 10% Owner Other (specify	
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE		(<i>'</i> ,	3. Date of Earliest Transaction (Month/Day/Year) 02/04/2020		below) SVP, Commercial &	below)	
(Street) BILLERICA (City)	MA (State)	01821 (Zip)	4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Indiv Line) X	ridual or Joint/Group Filing Form filed by One Repo Form filed by More thar Person	orting Person	

Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned

1. Title of Security (Instr. 3)	2. Transaction Date (Month/Day/Year)	if any						Securities	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
			Code	v	Amount	(A) or (D)	Price	Transaction(s) (Instr. 3 and 4)		(mour 4)
Common Stock	02/04/2020		S		55 ⁽¹⁾	D	\$26.71	14,184 ⁽²⁾	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)	4. Transa Code (8)		of Exp		Expiration Da	6. Date Exercisable and Expiration Date (Month/Day/Year)		and ht of ties ying tive ty (Instr. 3	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 148 RSUs.

2. Includes 11,245 restricted stock units.

Remarks:

<u>/s/ Brian Keane, Attorney-in-</u>

fact

02/04/2020

Date

** Signature of Reporting Person

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.