FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington,	D.C. 20549
-------------	------------

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

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL 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 ar	2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]									(Check	all app Direc	licable)	ng Person(s) to Is 10% Ov Other (s							
(Last) C/O QUA		3. Date of Earliest Transaction (Month/Day/Year) 06/16/2020									X	below) SVP R&D at				below)				
(Street) BILLERICA MA 01821 (City) (State) (Zip)					4. If Amendment, Date of Original Filed (Month/Day/Year)									6. Indiv Line) X	·					
		Table	I - No	n-Deriva	tive S	Secu	rities	Acq	uired,	, Dis	posed of	, or E	Benef	icially	Own	ed				ı
1. Title of Security (Instr. 3) 2. Transacti Date (Month/Day				Execut //Year) if any		Deemed ution Date, / th/Day/Year)				4. Securitie Disposed 0 5)			4 and Second Ben Owr		amount of curities neficially ned Following ported		n: Direct or Indirect nstr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)		
									Code	v	Amount	(A) (D)	or Pr	ice	Transa	saction(s) . 3 and 4)			(111341. 4)	ı
Common Stock 06/16/20					2020				S		1,667(1)	D \$		26.35	49,525(2)			D		ı
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. Transaction Code (Instr. 8)		of	r osed (, 3, 4	6. Date Exerci Expiration Dat (Month/Day/Ye		ite	7. Title and Amount of Securities Underlying Derivative Security (Ins 3 and 4)		Der Sec (Ins	Price of rivative curity str. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	Owner Form Direct or Ind (I) (In	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)	
					Code	v	(A)	(D)	Date Exercis	sable	Expiration Date	Title	Numb of Share							

Explanation of Responses:

- 1. The sale of shares reported on this Form 4 was made pursuant to a Rule 10b5-1 trading plan adopted on December 3, 2019.
- 2. Includes 13,006 restricted stock units.

Remarks:

/s/ Brian Keane, Attorney-in-

fact

** Signature of Reporting Person

Date

06/18/2020

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.