SEC Form 4

FORM 4

Check this box if no longer subject to Section 16. Form 4 or Form 5

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549 STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL

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

Instruction 1(b).	Filed	pursuant to Section 16(a) of the Securities Exchange Act of 1934		nours per response:		0.5	
			or Section 30(h) of the Investment Company Act of 1940					
	ress of Reporting Per KY E KEVIN	son*	2. Issuer Name and Ticker or Trading Symbol <u>Quanterix Corp</u> [QTRX]		ationship of Re < all applicable Director		on(s) to Issuer 10% Owner	
	(First) CRIX CORPORAT SEX TURNPIKE	(Middle) TION,	3. Date of Earliest Transaction (Month/Day/Year) 07/16/2020	x	Officer (give below) Chairma	10		у
(Street) BILLERICA (City)	MA (State)	01821 (Zip)	4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Indiv Line) X	Form filed b	by One Repor	(Check Applica rting Person One Reporting	

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

1. Title of Security (Instr. 3)	2. Transaction Date (Month/Day/Year)	2A. Deemed Execution Date, if any (Month/Day/Year)	3. Transaction Code (Instr. 8)					Securities Beneficially		7. Nature of Indirect Beneficial Ownership (Instr. 4)
			Code	v	Amount	(A) or (D)	Price	Transaction(s) (Instr. 3 and 4)		(11150.4)
Common Stock	07/16/2020		S ⁽¹⁾		4,500	D	\$27 . 94	1,066,840 ⁽²⁾	D	

Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities) 5. Number 1. Title of Derivative 3A. Deemed Execution Date, 6. Date Exercisable and Expiration Date 7. Title and Amount of 8. Price of Derivative 9. Number of derivative 10. Ownership 11. Nature of Indirect 3. Transaction Date Conversion Transaction (Month/Dav/Year) Derivative or Exercise if any Co (Month/Day/Year) 8) Code (Instr. (Month/Day/Year) Securities Security (Instr. 5) Securities Form: Beneficial Price of Securities Underlying Beneficially Direct (D) Ownership

(Instr. 3)		Price of Derivative Security		(Month/Day/Year)	8)		Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		(Underlying Derivative Security (Instr. 3 and 4)		(Instr. 5)	Beneficially Owned Following Reported Transaction(s) (Instr. 4)	Direct (D) or Indirect (I) (Instr. 4)	Ownership (Instr. 4)	,
		Code			v	(A)	(D)	Date Exercisable	Expiration Date	Title	Amount or Number of Shares						

Explanation of Responses:

1. The sale of shares was effected pursuant to a Rule 10b5-1 trading plan adopted on December 5, 2019.

2. Includes 178,132 restricted stock units and 39,815 shares of common stock that are subject to performance based vesting.

Remarks:

Security

/s/ Brian Keane, as Attorney-07/17/2020 in-Fact

** 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.