SEC Form 4	
------------	--

П

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

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 APPROVAL OMB Number: 3235-0287

L	Estimated average burden	
	hours per response:	0.5

1. Name and Address of Reporting Person [*] Connolly John M			2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]		tionship of Reporting Pe all applicable) Director	erson(s) to Issuer 10% Owner	
(Last) (First) (Middle) C/O BAIN CAPITAL VENTURE INVESTORS, LLC		, , , , , , , , , , , , , , , , , , ,	3. Date of Earliest Transaction (Month/Day/Year) 08/07/2018		Officer (give title below)	Other (specify below)	
200 CLARENDON STREET			4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Individual or Joint/Group Filing (Check Applicat Line)			
(Street)				X	Form filed by One Re	porting Person	
BOSTON	MA	02116			Form filed by More th Person	an One Reporting	
(City)	(State)	(Zip)					

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)	Transaction Code (Instr.		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 and 5)			5. Amount of Securities Beneficially Owned Following Reported	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)		(1150.4)
Common Stock	08/07/2018		A		2,270 ⁽¹⁾	Α	\$0.00	2,270 ⁽²⁾	D	

Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities)

	(-3)														
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)		n of 🛛		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)	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				
Stock Option (Right to Buy)			12/31/2018	08/07/2028	Common Stock	7,900	\$0.00	7,900	D						

Explanation of Responses:

1. Represents restricted stock units that vest as to 100% of the shares on December 31, 2018. Each restricted stock unit represents the right to receive one share of the Issuer's common stock.

2. Represents 2,270 restricted stock units.

Remarks:

/s/ John M. Connolly

** Signature of Reporting Person

08/09/2018

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.

Date