FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

asnington, D.C. 20549

OMB APPROVAL										
OMB Number:	3235-028									
Estimated average b	ourden									

0.5

hours per response:

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

Name and Address of Reporting Person* Connolly John M					2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]							(Ch	elationship ceck all applic	able) r	g Perso	10% Ow	ner
(Last) (First) (Middle) C/O BAIN CAPITAL VENTURE INVESTORS, LLC					3. Date of Earliest Transaction (Month/Day/Year) 01/02/2020								Officer below)	Officer (give title below)		Other (s below)	pecify
200 CLARENDON STREET					4. If Amendment, Date of Original Filed (Month/Day/Year)							Individual or Joint/Group Filing (Check Applicable Line)					
(Street) BOSTON MA 02116													Form fi	Form filed by One Reporting Person Form filed by More than One Reporting Person			
(City)	(S	tate)	(Zip)														
		Tal	ole I - Non	-Deriva	tive S	ecurities	Acq	uired,	Dis	posed o	f, or Ber	neficiall	y Owned				
Date			2. Transac Date (Month/Da	Execution Date,		Execution Date, f any		Transaction Disposed Code (Instr. 5)		ties Acquired (A) o d Of (D) (Instr. 3, 4		5. Amour Securitie Beneficia Owned F Reported	es ally Following	Form:	Direct I Indirect E str. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
							Code	v	Amount	Amount (A) or (D)		Transact	ansaction(s) nstr. 3 and 4)		(
Common Stock 01/02				01/02/	2/2020		A		2,270	1) A	\$0.00	9,510(2)			D		
			Table II - C			curities <i>i</i> Ils, warra							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	ate, Transacti Code (Ins				6. Date Exercisable and Expiration Date (Month/Day/Year)		е	7. Title and Amount of Securities Underlying Derivative Secu (Instr. 3 and 4)		8. Price of Derivative Security (Instr. 5)	9. Number derivative Securities Beneficial Owned Following Reported	lly	Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership (Instr. 4)
				Cod	de V	(A)		Date Exercisal		Expiration Date	Title	Amount or Number of Shares		Transactio (Instr. 4)	אונס)		
Stock Option (Right to Buy)	\$23.12	01/02/2020		A		7,900 ⁽³⁾		12/31/202	20	01/02/2030	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, 2020 granted pursuant to the Quanterix Corporation 2018 Non-Employee Director Compensation Policy. Each restricted stock unit represents the right to receive one share of the Issuer's common stock.
- 2. Includes 2,270 restricted stock units.
- 3. Represents non-qualified stock option to purchase 7,900 shares of common stock granted pursuant to the Quanterix Corporation 2018 Non-Employee Director Compensation Policy.

Remarks:

<u>/s/ John M. Connolly</u> <u>01/06/2020</u>

** 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.