

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

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

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

Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934 or Section 30(h) of the Investment Company Act of 1940

<p>1. Name and Address of Reporting Person*</p> <p><u>ARCH VENTURE FUND VI LP</u></p> <hr/> <p>(Last) (First) (Middle)</p> <p><u>8755 WEST HIGGINS ROAD, SUITE 1025</u></p> <hr/> <p>(Street)</p> <p><u>CHICAGO IL 60631</u></p> <hr/> <p>(City) (State) (Zip)</p>	<p>2. Issuer Name and Ticker or Trading Symbol</p> <p><u>Quanterix Corp [QTRX]</u></p> <hr/> <p>3. Date of Earliest Transaction (Month/Day/Year)</p> <p><u>12/11/2017</u></p> <hr/> <p>4. If Amendment, Date of Original Filed (Month/Day/Year)</p>	<p>5. Relationship of Reporting Person(s) to Issuer (Check all applicable)</p> <p>Director <input checked="" type="checkbox"/> 10% Owner</p> <p>Officer (give title below) <input type="checkbox"/> Other (specify below)</p> <hr/> <p>6. Individual or Joint/Group Filing (Check Applicable Line)</p> <p><input checked="" type="checkbox"/> Form filed by One Reporting Person</p> <p><input type="checkbox"/> Form filed by More than One Reporting Person</p>
--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------	-------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------	-------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------

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)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 and 5)			5. Amount of Securities Beneficially Owned Following Reported Transaction(s) (Instr. 3 and 4)	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			
Common Stock	12/11/2017		C		398,257	A	(1)	398,257	D ⁽²⁾	
Common Stock	12/11/2017		C		1,045,426	A	(1)	1,443,683	D ⁽²⁾	
Common Stock	12/11/2017		C		477,042	A	(1)	1,920,725	D ⁽²⁾	
Common Stock	12/11/2017		C		109,769	A	(1)	2,030,494	D ⁽²⁾	
Common Stock	12/11/2017		C		381,507	A	(1)	2,412,001	D ⁽²⁾	
Common Stock	12/11/2017		C		1,271,684	A	(1)	1,271,684	D ⁽³⁾	
Common Stock	12/11/2017		P		266,000 ⁽⁴⁾	A	\$15	1,537,684	D ⁽³⁾	

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)		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		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				
Series A-1 Preferred Stock	(1)	12/11/2017		C			1,280,000	(5)	(6)	Common Stock	398,257	(1)	0	D ⁽²⁾	
Series A-2 Preferred Stock	(1)	12/11/2017		C			3,360,000	(5)	(6)	Common Stock	1,045,426	(1)	0	D ⁽²⁾	
Series B Preferred Stock	(1)	12/11/2017		C			1,533,214	(5)	(6)	Common Stock	477,042	(1)	0	D ⁽²⁾	
Series C Preferred Stock	(1)	12/11/2017		C			352,800	(5)	(6)	Common Stock	109,769	(1)	0	D ⁽²⁾	
Series D Preferred Stock	(1)	12/11/2017		C			1,226,158	(5)	(6)	Common Stock	381,507	(1)	0	D ⁽²⁾	
Series D Preferred Stock	(1)	12/11/2017		C			4,087,193	(5)	(6)	Common Stock	1,271,684	(1)	0	D ⁽³⁾	

<p>1. Name and Address of Reporting Person*</p> <p><u>ARCH VENTURE FUND VI LP</u></p> <hr/> <p>(Last) (First) (Middle)</p> <p><u>8755 WEST HIGGINS ROAD, SUITE 1025</u></p> <hr/> <p>(Street)</p> <p><u>CHICAGO IL 60631</u></p> <hr/> <p>(City) (State) (Zip)</p>	<p>1. Name and Address of Reporting Person*</p>
--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------	-------------------------------------------------

[ARCH Venture Fund VIII Overage, L.P.](#)

(Last) (First) (Middle)
8755 WEST HIGGINS ROAD, SUITE 1025

(Street)
CHICAGO IL 60631

(City) (State) (Zip)

1. Name and Address of Reporting Person*

[ARCH VENTURE PARTNERS VI LP](#)

(Last) (First) (Middle)
8755 WEST HIGGINS ROAD, SUITE 1025

(Street)
CHICAGO IL 60631

(City) (State) (Zip)

1. Name and Address of Reporting Person*

[ARCH VENTURE PARTNERS VI LLC](#)

(Last) (First) (Middle)
8755 WEST HIGGINS ROAD, SUITE 1025

(Street)
CHICAGO IL 60631

(City) (State) (Zip)

1. Name and Address of Reporting Person*

[ARCH Venture Partners VIII, LLC](#)

(Last) (First) (Middle)
8755 WEST HIGGINS ROAD, SUITE 1025

(Street)
CHICAGO IL 60631

(City) (State) (Zip)

1. Name and Address of Reporting Person*

[BYBEE CLINTON](#)

(Last) (First) (Middle)
8755 WEST HIGGINS ROAD, SUITE 1025

(Street)
CHICAGO IL 60631

(City) (State) (Zip)

1. Name and Address of Reporting Person*

[NELSEN ROBERT](#)

(Last) (First) (Middle)
8755 WEST HIGGINS ROAD, SUITE 1025

(Street)
CHICAGO IL 60631

(City) (State) (Zip)

Explanation of Responses:

1. The shares of preferred stock automatically converted into common stock immediately prior to the closing of the Issuer's initial public offering using a calculation defined in the Issuer's Restated Certificate of Incorporation, as amended, based on the initial purchase price and the conversion rate in effect at the time of conversion. The shares of preferred stock automatically converted on a 1-for-3.214 basis and such conversion rate is reflected in the amount of common stock underlying the security.
2. The shares are held of record by ARCH Venture Fund VI, L.P. ("ARCH Fund VI"). The sole general partner of ARCH Fund VI is ARCH Venture Partners VI, L.P. ("ARCH Partners VI"), which may be deemed to beneficially own the shares held by ARCH Fund VI. The sole general partner of ARCH Partners VI is ARCH Venture Partners VI, LLC ("ARCH VI LLC"), which may be deemed to beneficially own the shares held by ARCH Fund VI. ARCH Partners VI and ARCH VI LLC disclaim beneficial ownership of such shares, except to the extent of any pecuniary interest therein. As a managing director of ARCH VI LLC, the Reporting Person may be deemed to beneficially own the shares held by ARCH Fund VI. The Reporting Person disclaims beneficial ownership of such shares, except to the extent of any pecuniary interest therein.
3. The shares are held of record by ARCH Venture Fund VIII Overage, L.P. ("ARCH Fund Overage"). The sole general partner of ARCH Fund Overage is ARCH Venture Partners VIII, LLC ("ARCH VIII LLC"), which may be deemed to beneficially own the shares held by ARCH Fund Overage. ARCH VIII LLC disclaims beneficial ownership of such shares, except to the extent of any pecuniary interest therein. As a

managing director of ARCH VIII LLC, the Reporting Person may be deemed to beneficially own the shares held by ARCH Fund Overage. The Reporting Person disclaims beneficial ownership of such shares, except to the extent of any pecuniary interest therein.

4. Reflects shares purchased in the Issuer's initial public offering.

5. These shares were convertible into the Issuer's common stock in accordance with the Issuer's Restated Certificate of Incorporation, as amended, at any time after the issuance of such shares, at the holder's election.

6. Not applicable.

/s/ Mark McDonnell, as
Attorney-in-Fact for ARCH 12/11/2017
Venture Fund VI, L.P.

/s/ Mark McDonnell, as
Attorney-in-Fact for ARCH 12/11/2017
Venture Fund VIII Overage, L.P.

/s/ Mark McDonnell, as
Attorney-in-Fact for ARCH 12/11/2017
Venture Partners VI, L.P.

/s/ Mark McDonnell, as
Attorney-in-Fact for ARCH 12/11/2017
Venture Partners VI, LLC

/s/ Mark McDonnell, as
Attorney-in-Fact for ARCH 12/11/2017
Venture Partners VIII, LLC

/s/ Mark McDonnell, as
Attorney-in-Fact for Clinton 12/11/2017
Bybee

/s/ Mark McDonnell, as
Attorney-in-Fact for Robert 12/11/2017
Nelsen

** 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.