

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

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

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

1. Name and Address of Reporting Person* <u>BAIN CAPITAL VENTURE INVESTORS, LLC</u> (Last) (First) (Middle) 200 CLARENDON STREET (Street) BOSTON MA 02116 (City) (State) (Zip)	2. Issuer Name and Ticker or Trading Symbol <u>Quanterix Corp [QTRX]</u>	5. Relationship of Reporting Person(s) to Issuer (Check all applicable) Director <input checked="" type="checkbox"/> 10% Owner Officer (give title below) Other (specify below)
	3. Date of Earliest Transaction (Month/Day/Year) 12/11/2017	
	4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Individual or Joint/Group Filing (Check Applicable Line) Form filed by One Reporting Person <input checked="" type="checkbox"/> Form filed by More than One Reporting Person

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		394,273	A	(1)	394,273	I	See Footnotes ⁽²⁾ (3)(4)(5)
Common Stock	12/11/2017		C		1,034,971	A	(1)	1,429,244	I	See Footnotes ⁽²⁾ (3)(4)(5)
Common Stock	12/11/2017		C		472,271	A	(1)	1,901,515	I	See Footnotes ⁽²⁾ (3)(4)(5)
Common Stock	12/11/2017		C		108,670	A	(1)	2,010,185	I	See Footnotes ⁽²⁾ (3)(4)(5)
Common Stock	12/11/2017		C		125,900	A	(1)	2,136,085	I	See Footnotes ⁽²⁾ (3)(4)(5)

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			1,267,200	(6)	(7)	Common Stock	394,273	(1)	0	I	See Footnotes ⁽²⁾⁽³⁾ (4)(5)
Series A-2 Preferred Stock	(1)	12/11/2017		C			3,326,401	(6)	(7)	Common Stock	1,034,971	(1)	0	I	See Footnotes ⁽²⁾⁽³⁾ (4)(5)
Series B Preferred Stock	(1)	12/11/2017		C			1,517,880	(6)	(7)	Common Stock	472,271	(1)	0	I	See Footnotes ⁽²⁾⁽³⁾ (4)(5)
Series C Preferred Stock	(1)	12/11/2017		C			349,271	(6)	(7)	Common Stock	108,670	(1)	0	I	See Footnotes ⁽²⁾⁽³⁾ (4)(5)
Series D Preferred Stock	(1)	12/11/2017		C			404,632	(6)	(7)	Common Stock	125,900	(1)	0	I	See Footnotes ⁽²⁾⁽³⁾ (4)(5)

1. Name and Address of Reporting Person*
BAIN CAPITAL VENTURE INVESTORS, LLC
 (Last) (First) (Middle)
 200 CLARENDON STREET
 (Street)
 BOSTON MA 02116
 (City) (State) (Zip)

1. Name and Address of Reporting Person*
BAIN CAPITAL VENTURE PARTNERS 2005, L.P.
 (Last) (First) (Middle)
 200 CLARENDON STREET

(Street)
BOSTON MA 02116

(City) (State) (Zip)

1. Name and Address of Reporting Person*

[BAIN CAPITAL VENTURE FUND 2005, L.P.](#)

(Last) (First) (Middle)
200 CLARENDON STREET

(Street)
BOSTON MA 02116

(City) (State) (Zip)

1. Name and Address of Reporting Person*

[BCIP Associates III, LLC](#)

(Last) (First) (Middle)
200 CLARENDON STREET

(Street)
BOSTON MA 02116

(City) (State) (Zip)

1. Name and Address of Reporting Person*

[BCIP Associates III](#)

(Last) (First) (Middle)
200 CLARENDON STREET

(Street)
BOSTON MA 02116

(City) (State) (Zip)

1. Name and Address of Reporting Person*

[BCIP Associates III-B, LLC](#)

(Last) (First) (Middle)
200 CLARENDON STREET

(Street)
BOSTON MA 02116

(City) (State) (Zip)

1. Name and Address of Reporting Person*

[BCIP Associates III-B](#)

(Last) (First) (Middle)
200 CLARENDON STREET

(Street)
BOSTON MA 02116

(City) (State) (Zip)

1. Name and Address of Reporting Person*

[KRUPKA MICHAEL A](#)

(Last) (First) (Middle)
C/O BAIN CAPITAL VENTURE INVESTORS, LLC
200 CLARENDON STREET

(Street)
BOSTON MA 02116

(City) (State) (Zip)

1. Name and Address of Reporting Person*

[Agarwal Ajay](#)

(Last)	(First)	(Middle)
C/O BAIN CAPITAL VENTURE INVESTORS, LLC		
200 CLARENDON STREET		
(Street)		
BOSTON	MA	02116
(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. Bain Capital Venture Investors, LLC ("BCVI") is the general partner of Bain Capital Venture Partners 2005, L.P. ("BCVP"), which is the general partner of Bain Capital Venture Fund 2005, L.P. ("Fund 2005"). As a result, BCVP may be deemed to share voting and dispositive power with respect to the securities held by Fund 2005. BCVP disclaims beneficial ownership of such securities except to the extent of its pecuniary interest therein. Immediately prior to the closing of the Issuer's initial public offering on December 11, 2017, the 1,106,387 shares of the Issuer's Series A-1 Preferred Stock, 2,904,268 shares of the Issuer's Series A-2 Preferred Stock, 1,325,258 shares of the Issuer's Series B Preferred Stock, 304,948 shares of the Issuer's Series C Preferred Stock and 353,283 shares of the Issuer's Series D Preferred Stock held by Fund 2005 automatically converted into 344,239, 903,630, 412,339, 94,881 and 109,921 shares of the Issuer's common stock, respectively.
3. Immediately prior to the closing of the Issuer's initial public offering on December 11, 2017, the 157,639 shares of the Issuer's Series A-1 Preferred Stock, 413,802 shares of the Issuer's Series A-2 Preferred Stock, 188,823 shares of the Issuer's Series B Preferred Stock, 43,449 shares of the Issuer's Series C Preferred Stock and 50,336 shares of the Issuer's Series D Preferred Stock held by BCIP Associates III, LLC ("BCIP III"), whose manager is BCIP Associates III ("BCIPA III"), whose managing partner is Boylston Coinvestors, LLC ("Boylston"), automatically converted into 49,047, 128,749, 58,750, 13,518 and 15,663 shares of the Issuer's common stock, respectively. BCIPA III may be deemed to share voting and dispositive power with respect to the securities held by BCIP III. BCIPA III disclaims beneficial ownership of such securities except to the extent of its pecuniary interest therein.
4. Immediately prior to the closing of the Issuer's initial public offering on December 11, 2017, the 3,174 shares of the Issuer's Series A-1 Preferred Stock, 8,331 shares of the Issuer's Series A-2 Preferred Stock, 3,799 shares of the Issuer's Series B Preferred Stock, 874 shares of the Issuer's Series C Preferred Stock and 1,013 shares of the Issuer's Series D Preferred Stock held by BCIP Associates III-B, LLC ("BCIP III-B" and together with Fund 2005 and BCIP III, the "Bain Capital Entities"), whose manager is BCIP Associates III-B ("BCIPA III-B"), whose managing partner is Boylston, automatically converted into 987, 2,592, 1,182, 271 and 316 shares of the Issuer's common stock, respectively. BCIPA III-B may be deemed to share voting and dispositive power with respect to the securities held by BCIP III-B. BCIPA III-B disclaims beneficial ownership of such securities except to the extent of its pecuniary interest therein.
5. The governance, investment strategy and decision-making process with respect to the investments held by the Bain Capital Entities is directed by the Executive Committee of BCVI, which consists of Michael A. Krupka and Ajay Agarwal. As a result, BCVI and Messrs. Krupka and Agarwal may be deemed to share voting and dispositive power with respect to the securities held by the Bain Capital Entities. BCVI and Messrs. Krupka and Agarwal disclaim beneficial ownership of such securities except to the extent of their pecuniary interest therein.
6. 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.
7. Not applicable.

Remarks:

Bain Capital Venture Investors, LLC, By: /s/ Michael A. Krupka, Name: Michael A. Krupka, Title: Managing Director 12/11/2017

Bain Capital Venture Partners 2005, L.P., By: Bain Capital Venture Investors, LLC, its general partner, By: /s/ Michael A. Krupka, Name: Michael A. Krupka, Title: Managing Director 12/11/2017

Bain Capital Venture Fund 2005, L.P., By: Bain Capital Venture Partners 2005, L.P., its general partner, By: Bain Capital Venture Investors, LLC, its general partner, By: /s/ Michael A. Krupka, Name: Michael A. Krupka, Title: Managing Director 12/11/2017

BCIP Associates III, LLC, By: BCIP Associates III, its manager, By: Boylston Coinvestors, LLC, its managing partner, By: /s/ Michael A. Krupka, Name: Michael A. Krupka, Title: Authorized Signatory 12/11/2017

BCIP Associates III, By: Boylston Coinvestors, LLC, its managing partner, By: /s/ Michael A. Krupka, Name: Michael A. Krupka, Title: Authorized Signatory 12/11/2017

BCIP Associates III-B, LLC, By: BCIP Associates III-B, its manager, By: Boylston Coinvestors, LLC, its managing partner, By: /s/ Michael A. Krupka, Name: Michael A. Krupka, Title: Authorized Signatory 12/11/2017

BCIP Associates III-B, By: Boylston Coinvestors, LLC, its managing partner, By: /s/ Michael A. Krupka, Name: Michael A. Krupka, Title: Authorized Signatory 12/11/2017

/s/ Michael A. Krupka 12/11/2017

/s/ Ajay Agarwal 12/11/2017

** Signature of Reporting Person Date

* 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.