# **UNITED STATES** SECURITIES AND EXCHANGE COMMISSION WASHINGTON, D.C. 20549

## **SCHEDULE 13G**

**Under the Securities Exchange Act of 1934** 

(Amendment No. 1)\*

Quanterix Corporation	
(Name of Issuer)	
Common Stock	
(Title of Class of Securities)	
T47000404	
74766Q101	
(CUSIP Number)	
D. J. 24 2040	
December 31, 2018	
(Date of Event Which Requires Filing of this Statement)	

Check the appropriate box to designate the rule pursuant to which this Schedule is filed:

☐ Rule 13d-1(b)☐ Rule 13d-1(c)

**⊠** Rule 13d-1(d)

\*The remainder of this cover page shall be filled out for a reporting person's initial filing on this form with respect to the subject class of securities, and for any subsequent amendment containing information which would alter disclosures provided in a prior cover page.

The information required in the remainder of this cover page shall not be deemed to be "filed" for the purpose of Section 18 of the Securities Exchange Act of 1934 ("Act") or otherwise subject to the liabilities of that section of the Act, but shall be subject to all other provisions of the Act (however, see the Notes).

CUSIP No. 74	4766Q101			13G	Page 2 of 15 Pages				
		•							
		NAMES OF REPORTING PERSONS							
1	I.R.S.	I.R.S. IDENTIFICATION NOS. OF ABOVE PERSONS (ENTITIES ONLY)  ARCH Venture Fund VI, L.P.							
	ARCH								
	CHEC	CHECK THE APPROPRIATE BOX IF A MEMBER OF A GROUP*							
2					(a)□				
					(b)□				
3	SEC U	SE ONI	L <b>Y</b>						
3									
	CITIZ	ENSHII	P OR PLACE (	OF ORGANIZATION					
4	Delawa	are							
			COLEVOT	ING POWER					
		5	SOLE VOI	ING POWER					
		J	0						
			SHARED V	OTING POWER					
NUMBER OF SI BENEFICIAL	LLY	6	3,949,684						
OWNED BY E REPORTING PI			SOLE DISP	OSITIVE POWER					
WITH	ZKSON	7	0						
	-		CHAREDD	ICDOCITIVE DOWED					
		8	SHARED D	ISPOSITIVE POWER					
		U	3,949,684						
	AGGR	REGATE	E AMOUNT BI	ENEFICIALLY OWNED BY EACH REPORTING	G PERSON				
9	3,949,6	584							
	CHECK IF THE AGGREGATE AMOUNT IN ROW (9) EXCLUDES CERTAIN SHARES*								
10									
11	PERCENT OF CLASS REPRESENTED BY AMOUNT IN ROW (9)								
11	17.8%								
	ТҮРЕ	OF REI	PORTING PE	RSON*					
12	PN								

CUSIP No. 74	4766Q101			13G	Page 3 of 15 Pages				
		NAMES OF REPORTING PERSONS							
1	I.R.S. IDENTIFICATION NOS. OF ABOVE PERSONS (ENTITIES ONLY)  ARCH Venture Fund VIII Overage, L.P.								
	ARCH Venture Fund VIII Overage, L.P.								
	CHEC	CHECK THE APPROPRIATE BOX IF A MEMBER OF A GROUP*							
2					(a)□				
	OF C. I	ICE ON	T 77		(b)□				
3	SEC U	JSE ON	LY						
J									
	CITIZ	ENSHI	P OR PLACE	OF ORGANIZATION					
4	Delaw	are							
			SOLE VOT	ING POWER					
		5		INGTOWER					
		J	0						
				OTING POWER					
NUMBER OF SI BENEFICIAL OWNED BY E	LLY	6	3,949,684						
REPORTING PI	ERSON		ERSON		SOLE DISE	POSITIVE POWER			
WITH	7	7	0						
	-		SHARED D	ISPOSITIVE POWER					
		8	3,949,684						
				ENEFICIALLY OWNED BY EACH REPORTING	C DEDSON				
9	AGGREGATE AMOUNT B			ENEFICIALLI OWNED DI EACH REPURIIN	GFERSUN				
	3,949,684								
1.0	CHECK IF THE AGGREGATE AMOUNT IN ROW (9) EXCLUDES CERTAIN SHARES*								
10									
	PERC	ENT O	F CLASS REP	RESENTED BY AMOUNT IN ROW (9)					
11				• •					
		17.8%							
12	TYPE	OF RE	PORTING PE	RSON*					
14	PN	PN							

CUSIP No. 74	4766Q101			13G	Page 4 of 15 Pages			
		NAMES OF REPORTING PERSONS						
1	I.R.S.	IDENTI	IFICATION NO	OS. OF ABOVE PERSONS (ENTITIES ONLY)				
	ARCH	ARCH Venture Partners VI, L.P.						
	CHEC	CHECK THE APPROPRIATE BOX IF A MEMBER OF A GROUP*						
2					(a)□			
	SEC I	ISE ONI	IV		(b)□			
3	SECU	SE UNI	LI					
	CITIZ	ENSHI	P OR PLACE	OF ORGANIZATION				
4	Delaw	are						
			SOLE VOT	ING POWER				
		5	0					
			-					
NUMBER OF SI	IADEC			OTING POWER				
BENEFICIAL	LLY	6	3,949,684					
OWNED BY E			SOLE DISF	POSITIVE POWER				
WITH		7	0					
				ACDOCATE NO LIED				
		8	SHARED D	DISPOSITIVE POWER				
		U	3,949,684					
	AGGF	REGATI	E AMOUNT BI	ENEFICIALLY OWNED BY EACH REPORTING	G PERSON			
9	3,949,684							
	CHECK IF THE AGGREGATE AMOUNT IN ROW (9) EXCLUDES CERTAIN SHARES*							
10								
11	PERCENT OF CLASS REPRESENTED BY AMOUNT IN ROW (9)							
11	17.8%							
	TYPE	OF RE	PORTING PEI	RSON*				
12	PN							
	111							

CUSIP No. 74	766Q101	-		13G	Page 5 of 15 Pages				
		NAMES OF REPORTING PERSONS I.R.S. IDENTIFICATION NOS. OF ABOVE PERSONS (ENTITIES ONLY)							
1				·					
	ARCH	ARCH Venture Partners VI, LLC							
_	CHEC	CHECK THE APPROPRIATE BOX IF A MEMBER OF A GROUP*							
2					(a)□				
	SEC I	JSE ONI	I V		(b)□				
3	SEC C	SE ON	LI						
	CITIZ	ENSHI	P OR PLACE	OF ORGANIZATION					
4	Delaw	are							
			SOLE VOT	ING POWER					
		5		1.010.121					
			0						
NUMBER OF SE	TADEC	•	SHARED V	OTING POWER					
NUMBER OF SI BENEFICIAL	LLY	v	3,949,684						
OWNED BY E			SOLE DISF	POSITIVE POWER					
WITH	LKSON	7	0						
		8	SHARED D	ISPOSITIVE POWER					
		O	3,949,684						
_	AGGF	REGATE	E AMOUNT BI	ENEFICIALLY OWNED BY EACH REPORTING	G PERSON				
9	3,949,684								
10	CHECK IF THE AGGREGATE AMOUNT IN ROW (9) EXCLUDES CERTAIN SHARES*								
44	PERCENT OF CLASS REPRESENTED BY AMOUNT IN ROW (9)								
11	17.8%								
	TYPE	OF RE	PORTING PEI	RSON*					
12	00								
	00								

CUSIP No. 74	766Q101			13G	Page 6 of 15 Pages			
		NAMES OF REPORTING PERSONS I.R.S. IDENTIFICATION NOS. OF ABOVE PERSONS (ENTITIES ONLY)						
1	I.R.S.	IDENTI	IFICATION N					
	ARCH	ARCH Venture Partners VIII, LLC						
	CHEC	CHECK THE APPROPRIATE BOX IF A MEMBER OF A GROUP*						
2					(a)□			
	SEC I	ISE ONI	IV		(b)□			
3	SECU	SE UNI	LI					
	CITIZ	ENSHI	P OR PLACE	OF ORGANIZATION				
4	Delaw	are						
			SOLE VOT	TING POWER				
		5	0					
			-					
NUMBER OF SI	IADEC			OTING POWER				
BENEFICIAL	LLY	6	3,949,684					
OWNED BY E			SOLE DISE	POSITIVE POWER				
WITH		7	0					
				ACDOCATE NO LIED				
		8	SHARED D	DISPOSITIVE POWER				
		U	3,949,684					
	AGGF	REGATI	E AMOUNT B	ENEFICIALLY OWNED BY EACH REPORTING	G PERSON			
9	3,949,684							
	CHECK IF THE AGGREGATE AMOUNT IN ROW (9) EXCLUDES CERTAIN SHARES*							
10								
11	PERCENT OF CLASS REPRESENTED BY AMOUNT IN ROW (9)							
11	17.8%							
	TYPE	OF RE	PORTING PE	RSON*				
12	00							

CUSIP No. 74	4766Q101			13G	Page 7 of 15 Pages									
		NAMES OF REPORTING PERSONS  LD S. IDENTIFICATION NOS. OF ABOVE BEDSONS (ENTITIES ONLY)												
1	I.R.S. IDENTIFICATION NOS. OF ABOVE PERSONS (ENTITIES ONLY)  Keith Crandell													
	Keith (													
	CHEC	CHECK THE APPROPRIATE BOX IF A MEMBER OF A GROUP*												
2					(a)□ 									
	SEC I	JSE ON	IV		(b) □									
3	SEC C	JSE UNI	Lĭ											
_	CITIZ	ZENSHI	P OR PLACE	OF ORGANIZATION										
4	United	l States o	of America											
			SOLE VOT	ING POWER										
		5		INGTOWER										
			10,170											
		•	SHARED V	OTING POWER										
NUMBER OF SE BENEFICIAL	LLY	6	3,949,684											
OWNED BY E REPORTING PI		PERSON		PERSON	PERSON							SOLE DISE	POSITIVE POWER	
WITH								7	10,170					
				SHARED D	ISPOSITIVE POWER									
		8	3,949,684											
9	AGGREGATE AMOUNT B			ENEFICIALLY OWNED BY EACH REPORTING	G PERSON									
3	3,959,	854												
	CHECK IF THE AGGREGATE AMOUNT IN ROW (9) EXCLUDES CERTAIN SHARES*													
10														
11	PERCENT OF CLASS REPRESENTED BY AMOUNT IN ROW (9)													
	17.8%													
	TYPE	OF RE	PORTING PE	RSON*										
12	IN	IN												

CUSIP No. 74	766Q101			13G	Page 8 of 15 Pages				
		•							
		NAMES OF REPORTING PERSONS							
1	I.R.S. IDENTIFICATION NOS. OF ABOVE PERSONS (ENTITIES ONLY)  Clinton Bybee								
	Clinton								
	CHEC	CHECK THE APPROPRIATE BOX IF A MEMBER OF A GROUP*							
2					(a)□				
	SEC I	JSE ON	I V		(b)□				
3	SEC C	SE UNI	LI						
	CITIZ	ENSHI	P OR PLACE (	OF ORGANIZATION					
4	United	l States o	of America						
			SOLE VOT	ING POWER					
		5	0						
	_		-						
NUMBER OF SI	TADES	C	SHARED V	OTING POWER					
BENEFICIAL	LLY	Y G	3,949,684						
OWNED BY E			SOLE DISP	POSITIVE POWER					
WITH		7		7		0			
				MCDOCITIVE DOWED					
		8		DISPOSITIVE POWER					
		U	3,949,684						
	AGGF	REGATI	E AMOUNT BI	ENEFICIALLY OWNED BY EACH REPORTING	G PERSON				
9	3,949,0	684							
	CHECK IF THE AGGREGATE AMOUNT IN ROW (9) EXCLUDES CERTAIN SHARES*								
10									
11	PERC	ENT OI	F CLASS REPI	RESENTED BY AMOUNT IN ROW (9)					
11	17.8%								
	TYPE	OF RE	PORTING PEI	RSON*					
12	IN								

CUSIP No. 74	766Q101			13G	Page 9 of 15 Pages			
		NAMES OF REPORTING PERSONS						
1	I.R.S. IDENTIFICATION NOS. OF ABOVE PERSONS (ENTITIES ONLY)  Robert Nelsen							
	Robert							
	CHEC	CHECK THE APPROPRIATE BOX IF A MEMBER OF A GROUP*						
2					(a)□			
	SEC I	ISE ONI	I V		(b)□			
3	SEC C	SE UNI	LI					
	CITIZ	ENSHI	P OR PLACE (	OF ORGANIZATION				
4	United	States o	of America					
			SOLE VOT	TING POWER				
		5	0					
NUMBER OF SE	IADEC	C	SHARED V	OTING POWER				
BENEFICIAI	LLY	LY	3,949,684					
OWNED BY E REPORTING PI			SOLE DISP	POSITIVE POWER				
WITH	220011	7	0					
				MCDOCITIVE DOWED				
		8		DISPOSITIVE POWER				
		U	3,949,684					
	AGGF	REGATE	E AMOUNT BI	ENEFICIALLY OWNED BY EACH REPORTING	G PERSON			
9	3,949,0	684						
	CHECK IF THE AGGREGATE AMOUNT IN ROW (9) EXCLUDES CERTAIN SHARES*							
10								
11	PERCENT OF CLASS REPRESENTED BY AMOUNT IN ROW (9)							
11	17.8%							
	ТҮРЕ	OF RE	PORTING PEI	RSON*				
12	IN							

CUSIP No. 74766Q101	13G	Page 10 of 15 Pages

## Item 1(a). Name of Issuer

Quanterix Corporation (the "Issuer").

### Item 1(b). Address of Issuer's Principal Executive Offices

113 Hartwell Avenue, Lexington, MA 02421

### Item 2(a). Name of Person Filing

ARCH Venture Fund VI, L.P. ("ARCH Venture Fund VI"); ARCH Venture Fund VIII Overage, L.P. ("AVF VIII Overage"); ARCH Venture Partners VI, L.P. ("AVP VI LP"); ARCH Venture Partners VI, LLC ("AVP VI LLC"); ARCH Venture Partners VIII, LLC ("AVP VIII LLC") (collectively, the "Reporting Entities" and individually, each a "Reporting Entity"); and Keith Crandell ("Crandell"), Robert Nelsen ("Nelsen") and Clinton Bybee ("Bybee") (collectively, the "Managing Directors" and individually, each a "Managing Director"). The Reporting Entities and the Managing Directors collectively are referred to as the "Reporting Persons".

### Item 2(b). Address of Principal Business Office or, if none, Residence

8755 W. Higgins Avenue, Suite 1025, Chicago, IL 60631.

#### Item 2(c). Citizenship

ARCH Venture Fund VI, AVF VIII Overage and AVP VI LP are limited partnerships organized under the laws of the State of Delaware. AVP VI LLC and AVP VIII LLC are limited liability companies organized under the laws of the State of Delaware. Each Managing Director is a US citizen.

#### Item 2(d). Title of Class of Securities

Common stock, par value \$0.001 per share.

#### Item 2(e). CUSIP Number

74766Q101.

#### Item 3. If this statement is filed pursuant to Rule 13d-1(b), or 13d-2(b) or (c), Check Whether the Person Filing is a:

Not Applicable.

## Item 4. Ownership

(a) Amount beneficially owned:

ARCH Venture Fund VI is the record owner of 2,412,000 shares of Common Stock (the "ARCH VI Shares") and AVF VIII Overage is the record holder of 1,537,684 share of Common Stock (the "ARCH Overage Shares"; combined with ARCH VI Shares, the "Record Shares") as of December 31, 2018. AVP VI LP, as the sole general partner of ARCH Venture Fund VI, may be deemed to beneficially own the ARCH VI Shares. AVP VI LLC, as the sole general partner of AVP VI LP, may be deemed to beneficially own the ARCH VI Shares. AVP VIII LLC, as the sole general partner of AVF VIII Overage, may be deemed beneficially own the ARCH Overage Shares. As managing directors of AVP VI LLC and AVP VIII LLC, each Managing Director may also be deemed to share the power to direct the disposition and vote of the Record Shares. In addition, Crandell holds options to 15,800 shares of Common Stock, of which 10,170 are vested as of December 31, 2018, and 0 will vest within 60 days of December 31, 2018 (the "Vested Option Shares").

(b) Percent of class:

In the aggregate, the Reporting Persons beneficially own approximately 17.8% of the outstanding shares of common stock of the Issuer, based upon 22,197,779 shares of common stock outstanding as of November 1, 2018, as reported on the Issuer's Form 10-Q as filed with the Securities and Exchange Commission on November 7, 2018. For Crandell, the Vested Option Shares were included in the number of shares of common stock outstanding.

CUSIP	No. 747	766Q101	13G	Page 11 of 15 Pages			
	(c)	Number of shares as to which	such person has:				
		(i) Sole power to vote or to o	lirect the vote:				
		See line 5 of the cover she	eets.				
		(ii) Shared power to vote or t	to direct the vote:				
		See line 6 of the cover she	eets.				
		(iii) Sole power to dispose or	to direct the disposition:				
		See line 7 of the cover sh	eets.				
		(iv) Shared power to dispose	or to direct the disposition:				
		See line 8 of the cover sh					
Each Repor	rting Pers	on disclaims beneficial ownersh	ip of such shares of Common Stock except for the sh	ares, if any, such Reporting Person holds of record.			
Item 5.	Ownership of Five Percent or Less of a Class						
	Not Ap	pplicable.					
Item 6.	Owner	rship of More Than Five Perce	nt on Behalf of Another Person				
	Not Applicable.						
Item 7.	Identií	fication and Classification of th	ne Subsidiary Which Acquired the Security Being	Reported on By the Parent Holding Company			
	Not Applicable.						
Item 8.	Identif	fication and Classification of M	lembers of the Group				
	Not Ap	oplicable.					
Item 9.	Notice	of Dissolution of Group					
	Not Ap	oplicable.					
Item 10.	Certifi	ication					
	Not Ap	oplicable.					

## **SIGNATURE**

After reasonable inquiry and to the best of my knowledge and belief, I certify that the information set forth in this statement is true, complete and correct.

Dated: February 14, 2019

ARCH VENTURE FUND VI, L.P.
By: ARCH Venture Partners VI, L.P. its General Partner
By: ARCH Venture Partners VI, LLC its General Partner
By: *
Keith Crandell Managing Director
ARCH VENTURE PARTNERS VI, L.P.
By: ARCH Venture Partners VI, LLC its General Partner
By: *
Keith Crandell Managing Director
ARCH VENTURE PARTNERS VI, LLC
By:*
Keith Crandell Managing Director
*
Keith Crandell
*
Robert Nelsen
Trobert Treasen
*
Clinton Bybee
ARCH VENTURE FUND VIII OVERAGE, L.F
By: ARCH Venture Partners VIII, LLC its General Partner
By:*
Keith Crandell Managing Director
ARCH VENTURE PARTNERS VIII, LLC
D *

Keith Crandell Managing Director \* By: /s/ Mark McDonnell

Mark McDonnell as

Attorney-in-Fact

This Amendment No. 1 to Schedule 13G was executed by Mark McDonnell pursuant to Powers of Attorney attached hereto as Exhibit 2 and Exhibit 2.1 and incorporated herein by reference.

#### Exhibit 1

## **AGREEMENT**

Pursuant to Rule 13d-1-(k)(1) under the Securities Exchange Act of 1934, the undersigned hereby agree that only one statement containing the information required by Schedule 13G need be filed with respect to the ownership by each of the undersigned of shares of stock of Quanterix Corporation.

an original.

	This Agreement may be executed in any num	ber of counterparts, each of which shall be deemed
Dated:	February 14, 2019	ARCH VENTURE FUND VI, L.P.
		By: ARCH Venture Partners VI, L.P. its General Partner
		By: ARCH Venture Partners VI, LLC its General Partner
		By:*
		Keith Crandell Managing Director
		ARCH VENTURE PARTNERS VI, L.P.
		By: ARCH Venture Partners VI, LLC its General Partner
		By: *  Keith Crandell  Managing Director
		ARCH VENTURE PARTNERS VI, LLC
		By:*
		Keith Crandell Managing Director
		*
		Keith Crandell
		*
		Robert Nelsen
		*
		Clinton Bybee
		ARCH VENTURE FUND VIII OVERAGE, L.P.
		By: ARCH Venture Partners VIII, LLC its General Partner
		By:*
		Keith Crandell Managing Director
		ARCH VENTURE PARTNERS VIII, LLC

Keith Crandell

Managing Director

* By:	/s/ Mark McDonnell
Ü	Mark McDonnell as
	Attorney-in-Fact

This Agreement was executed by Mark McDonnell pursuant to Powers of Attorney attached hereto as <u>Exhibit 2 and Exhibit 2.1</u> and incorporated herein by reference.

Exhibit 2

#### **POWERS OF ATTORNEY**

KNOW ALL MEN BY THESE PRESENTS, that each person whose signature appears below hereby constitutes and appoints Mark McDonnell his true and lawful attorney-in-fact, with full power of substitution, to sign any and all instruments, certificates and documents that may be necessary, desirable or appropriate to be executed on behalf of himself as an individual or in his capacity as a direct or indirect general partner, director, officer or manager of any partnership, corporation or limited liability company, pursuant to section 13 or 16 of the Securities Exchange Act of 1934, as amended (the "Exchange Act"), and any and all regulations promulgated thereunder, and to file the same, with all exhibits thereto, and any other documents in connection therewith, with the Securities and Exchange Commission, and with any other entity when and if such is mandated by the Exchange Act or by the Financial Industry Regulatory Authority, granting unto said attorney-in-fact full power and authority to do and perform each and every act and thing necessary, desirable or appropriate, fully to all intents and purposes as he might or could do in person, thereby ratifying and confirming all that said attorney-in-fact, or his substitutes, may lawfully do or cause to be done by virtue hereof.

ARCH VENTURE FUND VI, L.P.

IN WITNESS WHEREOF, this Power of Attorney has been signed as of the 10th day of November, 2010.

/s/ Keith Crandell Keith Crandell

<u>/s/ Robert Nelsen</u> Robert Nelsen

/s/ Clinton Bybee Clinton Bybee

Exhibit 2.1

#### **POWERS OF ATTORNEY**

KNOW ALL MEN BY THESE PRESENTS, that each person whose signature appears below hereby constitutes and appoints Mark McDonnell his true and lawful attorney-in-fact, with full power of substitution, to sign any and all instruments, certificates and documents that may be necessary, desirable or appropriate to be executed on behalf of himself as an individual or in his capacity as a direct or indirect general partner, director, officer or manager of any partnership, corporation or limited liability company, pursuant to section 13 or 16 of the Securities Exchange Act of 1934, as amended (the "Exchange Act"), and any and all regulations promulgated thereunder, and to file the same, with all exhibits thereto, and any other documents in connection therewith, with the Securities and Exchange Commission, and with any other entity when and if such is mandated by the Exchange Act or by the Financial Industry Regulatory Authority, granting unto said attorney-in-fact full power and authority to do and perform each and every act and thing necessary, desirable or appropriate, fully to all intents and purposes as he might or could do in person, thereby ratifying and confirming all that said attorney-in-fact, or his substitutes, may lawfully do or cause to be done by virtue hereof.

IN WITNESS WHEREOF, this Power of Attorney has been signed as of the 26th day of July, 2017.

ARCH VENTURE FUND VIII OVERAGE, L.P.

By: ARCH Venture Partners VIII, LLC its General Partner

By: <u>/s/ Keith Crandell</u>

Managing Director

ARCH VENTURE PARTNERS VIII, L.P.

By: ARCH Venture Partners VIII, LLC its General Partner

By: /s/ Keith Crandell

Managing Director