FORM 4

## **UNITED STATES SECURITIES AND EXCHANGE COMMISSION**

Washington, D.C. 20549

## STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

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

Check this box if no longer subject to Section 16. Form 4 or Form 5 obligations may continue. See

Instruc	ion 1(b).			Filed							ities Exchangompany Act o		f 1934							
1. Name and Address of Reporting Person*  ARCH VENTURE FUND VI LP				2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [ QTRX ]							Relationship of Reporting Person(s) to Issuer (Check all applicable)     Director X 10% Owner									
(Last) (First) (Middle) 8755 WEST HIGGINS ROAD, SUITE 1025					3. Date of Earliest Transaction (Month/Day/Year) 11/20/2020							Officer (give title Other (specify below) below)								
(Street) CHICAGO IL 60631					4. If Amendment, Date of Original Filed (Month/Day/Year)						Individual or Joint/Group Filing (Check Applicable Line)     Form filed by One Reporting Person     Form filed by More than One Reporting Person									
(City)	(St	ate) (2	Zip)												1 0100					
		Table	I - No	on-Deriva	tive	Secu	rities	Acc	quired	l, Dis	sposed of	, or E	Benef	icia	lly Own	ed				
1. Title of Security (Instr. 3)				2. Transaction Date (Month/Day/Year)		Execution Date,		3. Transaction Code (Instr. 8)		4. Securities Acquired (A) Disposed Of (D) (Instr. 3, 4 5)				s illy ollowing	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)		7. Nature of Indirect Beneficial Ownership (Instr. 4)			
								Code	v	Amount	(A) o (D)	) or Price Transaction(s		on(s)			(IIISU	. 4)		
Common Stock			11/20/2020				<b>J</b> <sup>(1)</sup>		229,006	D	4	<b>60</b>	2,182	182,995 I		I	See Footnote <sup>(2)</sup>			
Common Stock				11/20/2020				<b>J</b> <sup>(1)</sup>		145,994	D	\$	<b>60</b>	1,391,690 I			See Footnote <sup>(3)</sup>			
Common Stock				11/20/2020				J <sup>(1)(4)</sup>		854	A	\$	<b>50</b>	85	54	D <sup>(5)</sup>				
Common Stock				11/20/2020				J <sup>(1)(4)</sup>		853	A	1	\$ <del>0</del>	85	53	I	)(6)			
Common Stock				11/24/2020				<b>J</b> <sup>(1)</sup>		277,514	D	\$	<b>50</b>	1,905	5,481		I	See Footnote <sup>(2)</sup>		
Common Stock				11/24/2020				<b>J</b> <sup>(1)</sup>		176,920	D	\$	<b>60</b>	1,214	1,770		I	See Footnote <sup>(3)</sup>		
Common Stock 11/24/2				11/24/2	.020						1,034	A	1	<b>60</b>	1,8	88	I	<b>)</b> <sup>(5)</sup>		
Common Stock 11/			11/24/2	11/24/2020				J <sup>(1)(4)</sup>		1,036	A	1	\$ <del>0</del>	1,8	89	D <sup>(6)</sup>				
		Та	ble II								osed of, convertib				y Owned	t				
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	Execu	eemed ition Date,	emed 4.		5. Number of		6. Date Exer		cisable and	7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		tr.	8. Price of Derivative Security (Instr. 5)	9. Numbo derivativ Securitie Beneficia Owned Followin Reported Transact (Instr. 4)	e es ally g d ion(s)	10. Ownersh Form: Direct (D or Indirec (I) (Instr.	nip of Ind Benef Owne ct (Instr.	11. Nature of Indirect Beneficial Ownership Instr. 4)
					Code	v	(A)	(D)	Date Exerci	sable	Expiration Date	Title	Amou or Numb of Share	er						
		Reporting Person																		
(Last) 8755 WI		(First) NS ROAD, SUI	`	liddle)																
(Street)						- 1														

60631

(Zip)

(Middle)

**CHICAGO** 

(Last)

IL

1. Name and Address of Reporting  $\mathsf{Person}^{^\star}$ 

(State)

ARCH Venture Fund VIII Overage, L.P.

(First)

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) 8755 WEST HIGO	(First) GINS ROAD, SUITE	(Middle)						
(Street) CHICAGO	IL	606031						
(City)	(State)	(Zip)						
1. Name and Address of Reporting Person*  ARCH VENTURE PARTNERS VI LLC								
(Last) 8755 WEST HIGO	(First) GINS ROAD, SUITE	(Middle) 1025						
(Street) CHICAGO	IL	606031						
(City)	(State)	(Zip)						
1. Name and Address of Reporting Person*  ARCH Venture Partners VIII, LLC								
(Last)	(First)	(Middle)						
8755 WEST HIGO	GINS ROAD, SUITE	1025						
(Street) CHICAGO	IL	606031						
(City)	(State)	(Zip)						
Name and Address of Reporting Person*     BYBEE CLINTON								
(Last)	(First)	(Middle)						
C/O ARCH VENTURE PARTNERS IX, LLC								
8755 W. HIGGINS ROAD, SUITE 1025								
(Street) CHICAGO	IL	60631						
(City)	(State)	(Zip)						
1. Name and Address of Reporting Person*  NELSEN ROBERT								
(Last) 8755 W. HIGGINS	(First) S ROAD, SUITE 102	(Middle)						
(Street) CHICAGO	IL	60631						
(City)	(State)	(Zip)						

## Explanation of Responses:

- 1. Distribution of Common Stock held by a limited partnership to its partners for no consideration.
- 2. These shares are directly held by ARCH Venture Fund VI, L.P. ("ARCH VI"). ARCH Venture Partners VI, L.P. ("GPLP"), the sole general partner of ARCH VI, may be deemed to beneficially own the shares held by ARCH VI. ARCH Venture Partners VI, LLC ("GPLLC"), as the sole general partner of GPLP, may be deemed to beneficially own the shares held by ARCH VI. As managing directors of GPLLC, each of Robert Nelsen and Clinton Bybee (the "ARCH Managing Directors"), may be deemed to beneficially own the shares held by ARCH VI. Each of GPLP, GPLLC and the ARCH Managing Directors disclaims beneficial ownership of the securities reported herein, except to the extent of their pecuniary interest therein, if any.
- 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 managing directors of ARCH VIII LLC, each of the ARCH Managing Directors may be deemed to beneficially own the shares held by ARCH Overage. Each of ARCH VIII LLC and the ARCH Managing Directors disclaims beneficial ownership of the securities reported herein, except to the extent of their pecuniary interest therein, if any.
- 4. Change from indirect to direct ownership of shares previously reported as beneficially owned by the Reporting Person.
- 5. Shares owned by Nelsen.
- 6. Shares owned by Bybee.

/s/ Mark McDonnell, as Attorney-in-Fact for ARCH Venture Fund VI, L.P.	11/24/2020
/s/ Mark McDonnell, as Attorney-in-Fact for ARCH Venture Fund VIII Overage, L.P.	11/24/2020
/s/ Mark McDonnell, as Attorney-in-Fact for ARCH Venture Partners VI, L.P.	11/24/2020
/s/ Mark McDonnell, as Attorney-in-Fact for ARCH Venture Partners VI, LLC	11/24/2020
/s/ Mark McDonnell, as Attorney-in-Fact for ARCH Venture Partners VIII, LLC	11/24/2020
/s/ Mark McDonnell, as Attorney-in-Fact for Clinton Bybee	11/24/2020
/s/ Mark McDonnell, as Attorney-in-Fact for Robert Nelsen	11/24/2020
** Signature of Reporting Person	Date

Reminder: Report on a separate line for each class of securities beneficially owned directly or indirectly.

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.

 $<sup>^{\</sup>star}$  If the form is filed by more than one reporting person, see Instruction 4 (b)(v).

<sup>\*\*</sup> Intentional misstatements or omissions of facts constitute Federal Criminal Violations See 18 U.S.C. 1001 and 15 U.S.C. 78ff(a).

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  $10 \, \mathrm{th}$  day of November, 2010.

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: /s/ Keith Crandell Managing Director

ARCH VENTURE PARTNERS VI, L.P.

By: ARCH Venture Partners VI, LLC its General Partner

By: /s/ Keith Crandell Managing Director

ARCH VENTURE PARTNERS VI, LLC

By: /s/ Keith Crandell Managing Director

/s/ Keith Crandell Keith Crandell

/s/ Robert Neslon Robert Nelsen

/s/ Clinton Bybee Clinton Bybee

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: /s/ Keith Crandell Managing Director

ARCH VENTURE PARTNERS VIII, L.P.

By: ARCH Venture Partners VIII, LLC

its General Partner

By: /s/ Keith Crandell Managing Director