SEC Form 4	
------------	--

FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

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).

## 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

1. Name and Address of Reporting Person*				suer Name <b>and</b> Tic <u>anterix Corp</u>		Symbol	5. Relationship of Reporting Person(s) to Issuer (Check all applicable)					
CRANDELL KEITH				r_			X	Director	Х	10%	Owner	
(Last)	(First)	(Middle)		ate of Earliest Trans 03/2022	saction (Montl	n/Day/Year)		Officer (give title below)		Othe belov	r (specify v)	
		ID, 5011L 1025		Amenderent Dete			C. In die	idual an laint/Cras		2h e el c d	Ameliaahla	
				Amenument, Date	oi Originai File	ed (Month/Day/Year)	6. Individual or Joint/Group Filing (Check Applicable Line)					
(Street) CHICAGO	IL	60631					X	Form filed by O	ne Reporti	ng Per	son	
	IL	00031						Form filed by Me Person	ore than C	ne Re	porting	
(City)	(State)	(Zip)										
	Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned											
Date			2. Transaction Date (Month/Day/Year)	2A. Deemed Execution Date, if any	3. Transaction Code (Instr.	4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 a	nd 5)	5. Amount of Securities Beneficially	6. Owner Form: Di (D) or Inc	rect	7. Nature of Indirect Beneficial	

			Code (Instr. 8)					Securities Beneficially Owned Following	(D) or Indirect (I) (Instr. 4)	Indirect Beneficial Ownership	
			Code	v	Amount	(A) or (D)	Price	Reported Transaction(s) (Instr. 3 and 4)		(Instr. 4)	
Common Stock	11/16/2021		J <sup>(1)</sup>		189,311	D	\$0.00	177,247	Ι	See footnote <sup>(2)</sup>	
Common Stock	11/16/2021		J <sup>(1)</sup>		120,689	D	\$0.00	113,003	Ι	See Footnote <sup>(3)</sup>	
Common Stock	11/16/2021		<b>J</b> (1)(4)		9,952	A	\$0.00	105,964	D		
Common Stock	11/18/2021		J <sup>(1)</sup>		177,247	D	\$0.00	0.00	I	See footnote <sup>(2)</sup>	
Common Stock	11/18/2021		J <sup>(1)</sup>		113,003	D	\$0.00	0.00	I	See Footnote <sup>(3)</sup>	
Common Stock	11/18/2021		J <sup>(1)(4)</sup>		9,318	A	\$0.00	115,282	D		
Common Stock	01/03/2022		A		2,668 <sup>(5)</sup>	A	\$0.00	117 <b>,9</b> 50 <sup>(6)</sup>	D		
Common Stock	01/03/2022		A		<b>307</b> <sup>(7)</sup>	A	\$42.41(8)	118,257(6)	D		

## 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. Transa Code ( 8)		Derivative		6. Date Exercisable and Expiration Date (Month/Day/Year)		I 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	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		Transaction(s) (Instr. 4)		
Stock Option (Right to Buy)	\$42.41	01/03/2022		A		4,002 <sup>(9)</sup>		12/31/2022	01/03/2032	Common Stock	4,002	\$0.00	4,002	D	

## 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 VITUF 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. The Reporting Person is a managing director of the GPLLC and shares voting and investment control with respect to these shares. The Reporting Person may be deemed to beneficially own the shares held by ARCH VI. The Reporting Person disclaims beneficial ownership of the securities reported herein, except to the extent of his pecuniary interest therein, if any.

3. These shares are directly held by ARCH Venture Fund VIII Overage, L.P. ("ARCH Overage"). ARCH Venture Partners VIII, LLC ("ARCH VIII LLC") the sole general partner of ARCH Overage, may be deemed to beneficially own the shares held by ARCH Overage. The Reporting Person is a managing director of the ARCH VIII LLC and shares voting and investment control with respect to these shares. The Reporting Person may be deemed to beneficially own the shares held by ARCH Overage. The Reporting Person disclaims beneficial ownership of the securities reported herein, except to the extent of his pecuniary interest therein, if any.

4. Change from indirect to direct ownership of shares previously reported as beneficially owned by the Reporting Person.

5. Represents restricted stock units that vest as to 100% of the shares on December 31, 2022 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.

6. Includes 2,668 restricted stock units.

7. Consists of stock granted in lieu of cash fees for service on the Company's Board of Directors and committees thereof for the fourth quarter of 2021.

8. Closing price of the Company's common stock on the Nasdaq Global Market on January 3, 2021.

9. Represents non-qualified stock option to purchase 4,002 shares of common stock granted pursuant to the Quanterix Corporation 2018 Non-Employee Director Compensation Policy.

Remarks:

/s/ Brian Keane, as Attorney-in-01/05/2022

Fact for Keith L. Crandell

\*\* 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.