Instruction 1(b).

FORM 4

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

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington,	D.C.	20549
-------------	------	-------

STATEMENT	OF CHANGES II	N BENEFICIAL	OWNERSHIP

	OMB APPROVAL									
- 1										
	OMB Number: 3235-028									
	Estimated average bu	ırden								
	hours per response:	0.5								

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*  OLSON LAURIE J					2. Issuer Name <b>and</b> Ticker or Trading Symbol Quanterix Corp [ QTRX ]									k all applic Directo	ctor		10% Ov	mer		
· ·	ANTERIX	CORPORATION	(Middle) <mark>V</mark> ,			3. Date of Earliest Transaction (Month/Day/Year) 04/27/2021									Officer below)	(give title		Other (s below)	specify	
900 MID	DLESEX	TURNPIKE			4.1	4. If Amendment, Date of Original Filed (Month/Day/Year)								Individual or Joint/Group Filing (Check Applicable Line)						
(Street) BILLER	ICA M	A	01821											X		•		orting Perso one Repo		
(City)	(S	tate)	(Zip)		-									Person						
	Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																			
Date				2. Trans Date (Month/		action 2A. Deemed Execution Date, if any (Month/Day/Year)		Transaction Dispose Code (Instr. 5)		Disposed	ities Acquired (A) or d Of (D) (Instr. 3, 4 and			5. Amou Securitie Beneficia Owned F Reported	s ally following	Forn ly (D) o		7. Nature of Indirect Beneficial Ownership (Instr. 4)		
								Code	v	Amount	(A) or (D) Pr		се	Transact (Instr. 3	action(s)					
Common Stock 04/27					7/202	7/2021		A		4,540	$A^{(1)}$ A		0.00	4,540(2)			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 Date Date (Month/Day/Year) Price of Derivative Security  3. Transaction Execution Date Execution Date, if any (Month/Day/Year) (Month/Day/Year)  4. Transaction Execution Date, if any (Month/Day/Year) (Month/Day/Year)			5. Numb of Derivati Securiti Acquire (A) or Dispose of (D) (II 3, 4 and	ve les ed ed nstr.	Expiration	Expiration Date of: (Month/Day/Year) Un De			7. Title and Amour of Securities Underlying Derivative Security (Instr. 3 and 4)		s. Price of Derivative Security Instr. 5)	9. Number derivative Securities Beneficial Owned Following Reported Transactic (Instr. 4)	Own For Illy Dire or Ii	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)				
					Code	v	(A)	(D)	Date Exercisal		Expiration Date	Title	Amou or Numl of Share	per						
Stock Option (right to buy)	\$77.79	04/27/2021			A		15,800		(3)	(	)4/27/2031	Common Stock	15,8	00	\$0.00	15,800	0	D		

## **Explanation of Responses:**

- 1. Represents shares of stock issuable on vesting of restricted stock units. The restricted stock units vest as to one-third of shares on each of April 27, 2022, April 27, 2023 and April 27, 2024.
- 2. Includes 4,540 restricted stock units
- 3. The option vests as to one-third of the shares on April 27, 2022, and the remaining two-thirds of the shares vest monthly thereafter over two years in 24 successive equal installments.

## Remarks:

/s/ Brian Keane, as Attorney-in-04/28/2021 **Fact** 

\*\* Signature of Reporting Person

Date

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

- $^{\star}$  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.