FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

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

## TATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

**OMB APPROVAL** OMB Number: Estimated average burden 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

Name and Address of Reporting Person*     Orticerio Ernest						2. Issuer Name <b>and</b> Ticker or Trading Symbol Quanterix Corp [ QTRX ]								eck all appli Direct Office	ationship of Reporting call applicable) Director Officer (give title		10% Ow Other (s	ner
•	ANTERIX	CORPORATION	(Middle) <mark>N</mark> ,				ate of Earliest Transaction (Month/Day/Year) 0/2018							below	SVP Finance and Co		below) orporate Dev	
113 HARTWELL AVENUE				4.1	If Amendment, Date of Original Filed (Month/Day/Year)								6. Individual or Joint/Group Filing (Check Applicable					
(Street)	TON M	ÍΑ	02421										Line	X Form	•		orting Person	
(City)	(S	tate)	(Zip)		-									Perso		re mai	гопе керог	urig
		Tab	le I - Nor	n-Deri	vativ	e Se	curities	s Ac	quired,	Dis	osed o	f, or Be	neficial	ly Owned	I			
Dat			Date	e nth/Day/Year) i		2A. Deemed Execution Date, if any (Month/Day/Year)		Transaction Disp Code (Instr. 5)		Disposed	ırities Acquired (A) or ed Of (D) (Instr. 3, 4 an		Benefici	es ally Following	Form (D) o	: Direct r Indirect str. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
									Code	v	Amount	(A) o (D)	Price	Transac	nnsaction(s) str. 3 and 4)			(msu. 4)
Common Stock 05			05/1	10/201	/2018		A		7,131 <sup>(1)</sup> A		\$0	53,802(2)			D			
		7	Гable II -									or Ben ble secu		Owned				
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date Month/Day/Year)  (Month/Day  (Month/Da	Date,	4. Transa Code ( 8)				6. Date Ex Expiration (Month/Da	Date		of Securit Underlyin	g Security	8. Price of Derivative Security (Instr. 5)	9. Number derivative Securities Beneficia Owned Following Reported Transacti (Instr. 4)	e s illy	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership t (Instr. 4)	
					Code	v	(A)	(D)	Date Exercisab		expiration Date	Title	Amount or Number of Shares					
Stock Option (right to buy)	\$19.12	05/10/2018			A		24,311		(3)	O	5/10/2028	Common Stock	24,311	\$0	24,31	.1	D	

## **Explanation of Responses:**

- 1. Represents shares of common stock issuable upon vesting of restricted stock units. The restricted stock units vest as to 25% of the shares on May 10, 2019 and vest as to an additional 2.083% of the shares per
- 2. Represents 46,671 shares of common stock and 7,131 restricted stock units.
- 3. The option vests as to 25% of the shares on May 10, 2019 and vests as to an additional 2.083% of the shares per month for 36 months thereafter.

/s/ Brian Keane, as Attorney-in-05/14/2018 **Fact** 

\*\* 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.