FORM 4

## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

## OMB APPROVAL

- 1		
	OMB Number:	3235-0287
	Estimated average burd	en
	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*						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [ QTRX ]								5. Relationship of Reporting Person(s) to Issuer (Check all applicable)					
<u>DEKKERS MARIJN E</u>						Zamiram gorb [ drint ]								X Directo	or		10% Ov	ner	
(Last)	(F	First)	(Middle)			3. Date of Earliest Transaction (Month/Day/Year)								Officer below)	(give title		Other (s below)	pecify	
C/O QUANTERIX CORPORATION,						12/11/2017													
113 HARTWELL AVENUE						4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Individual or Joint/Group Filing (Check Applicable					
(Street)					_								Lin	,	iled by One	Reno	rting Persor	,	
LEXINGTON MA 02421													X Form filed by One Reporting Person  Form filed by More than One Reporting  Person						
(City)	(\$	State)	(Zip)																
		Tal	ble I - Nor	ı-Deriv	vativ	e Se	curi	ties Acc	quired,	Dis	posed o	f, or Be	neficial	ly Owned					
1. Title of Security (Instr. 3)  2. Transa Date (Month/E					/Day/Year)		2A. Deemed Execution Date, if any (Month/Day/Year)		Transaction Dis		Disposed	urities Acquired (A) sed Of (D) (Instr. 3,		Benefici	es ally Following	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)		7. Nature of Indirect Beneficial Ownership	
									Code	v	Amount	(A) o (D)	r Price	Transact (Instr. 3	ion(s)			Instr. 4)	
Common Stock 12/11.					1/201	/2017		С		77,37	78 A		77,378			D			
			Table II -								osed of, onvertib			Owned					
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	Conversion Date or Exercise (Month/Day/Year) Price of Derivative				ction Instr.	5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		6. Date Exercisable Expiration Date (Month/Day/Year)		е	7. Title and An of Securities Underlying Derivative Sec (Instr. 3 and 4)		8. Price of Derivative Security (Instr. 5)		e s ally g	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership (Instr. 4)	
					Code	v	(A)	(D)	Date Exercisa		Expiration Date		Amount or Number of Shares		(Instr. 4)	5.1(3)			
Series D-1 Preferred Stock	(1)	12/11/2017			С			248,694	(2)		(3)	Common Stock	77,378	(1)	0		D		

## **Explanation of Responses:**

- 1. The shares of preferred stock automatically converted into common stock immediately prior to the closing of the Issuer's initial public offering using a calculation defined in the Issuer's Restated Certificate of Incorporation, as amended, based on the initial purchase price and the conversion rate in effect at the time of conversion. The shares of preferred stock automatically converted on a 1-for-3.214 basis and such conversion rate is reflected in the amount of common stock underlying the security.
- 2. These shares were convertible into the Issuer's common stock in accordance with the Issuer's Restated Certificate of Incorporation, as amended, at any time after the issuance of such shares, at the holder's election.
- 3. Not applicable.

/s/ Brian Keane, as Attorney-in-Fact 12/11/2017

\*\* 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.