(City)

(State)

(Zip)

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

OMB APPROVAL	

OMB Number:	3235-0287
Estimated average burder	n
hours per response:	0.5

⊫

5. Relationship of Reporting Person(s) to Issuer (Check all applicable)

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

1. Name and Address of Reporting Person*

FORM 4

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

2. Issuer Name and Ticker or Trading Symbol

<u>Quanterix Corp</u> [QTRX]

<u>Flagsh</u>	<u>ip Ventur</u>	<u>es Fund 2004</u>	<u>L P</u>		Qua	antei	<u>1x Corp</u>	[QTRX]					Director		Х	10% O	wner
				. Date of Earliest Transaction (Month/Day/Year) 2/11/2017							Officer (give title Other (specify below) below)						
(Street) CAMBRIDGE MA 02142 (City) (State) (Zip)					If Amendment, Date of Original Filed (Month/Day/Year)							 6. Individual or Joint/Group Filing (Check Applicable Line) Form filed by One Reporting Person X Form filed by More than One Reporting Person 					
(0.13)		,		n Dori	(otivo	500	urition A	oquirod I		nocod	of or P	noficially	, Ownod				
Table I - Non-Derivative S 1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Year)			2A Ex (r) if a	3. Transac	3. Transaction Code (Instr.			red (A) or str. 3, 4 and 5	5. Amount of Securities Beneficially Owned Following		6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)		7. Nature of Indirect Beneficial Ownership				
								Code	v	Amount	(A) (D)	or Price	Reported Transactio (Instr. 3 a	on(s)			(Instr. 4)
Commor	Stock			12/1	L/2017		C			398,257 A		A (1)	398,257		D ⁽²⁾		
Commor	Stock			12/1	L/2017	T		С		1,045,426 A		A (1)	1,443	1,443,683		D ⁽²⁾	
Commor	Stock			12/1	l/2017			С		477,0	42	A (1)	1,920),725	D ⁽²⁾		
Commor	Stock			12/1	l/2017			С		109,7	70	A (1)	2,030	2,030,495		D ⁽²⁾	
			Table II -	Deriva	tive S	Secu	rities Aco	quired, Di	isp	osed of	, or Be	neficially	Owned				
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	3A. Deemed Execution Da if any (Month/Day/Y	te, 4. Tra	nsactio de (Insti	5. De Se Ac or (D	Number of rivative curities quired (A) Disposed of ((Instr. 3, 4 d 5)	6. Date Exe Expiration (Month/Day	ercis Date	e Securities Under		nd Amount of s Underlying e Security	8. Price of Derivative Security (Instr. 5)	derivati Securiti	curities Form neficially Dire whed or In llowing (I) (II ported		Beneficial Ownershi t (Instr. 4)
				Co	de V	(A)) (D)	Date Exercisable		Expiration Date	Title	Amount or Number of Shares		(Instr. 4)			
Series A-1 Preferred Stock	(1)	12/11/2017		0	:		1,280,000	(3)		(4)	Common Stock	398,257	(1)	0		D ⁽²⁾	
Series A-2 Preferred Stock	(1)	12/11/2017		0	;		3,360,000	(3)		(4)	Common Stock	1,045,42	6 (1)	0	1	D ⁽²⁾	
Series B Preferred Stock	(1)	12/11/2017		0	;		1,533,214	(3)	T	(4)	Common Stock	477,042	(1)	0		D ⁽²⁾	
Series C Preferred Stock	(1)	12/11/2017		(:		352,800	(3)	T	(4)	Common Stock	109,770	(1)	0		D ⁽²⁾	
Flagsh (Last) 55 CAM (Street) CAMBE (City) 1. Name a Flagsh (Last)	ip Ventur IBRIDGE F RIDGE nd Address o ip Ventur	f Reporting Person [*] es Fund 2004 (First) PARKWAY, SUI MA (State) f Reporting Person [*] es General Pa (First) PARKWAY, SUI	LP (Middl FE 800E 0214 (Zip) rttner LLC (Middl	2													
(Street) CAMBI	RIDGE	MA	0214	2													

1. Name and Address of Reporting Person* AFEYAN NOUBAR								
(Last)	(First)	(Middle)						
55 CAMBRIDGE PARKWAY, SUITE 800E								
(Street) CAMBRIDGE	MA	02142						
(City)	(State)	(Zip)						
1. Name and Address of Reporting Person [*] KANIA EDWIN M JR								
(Last)	(First)	(Middle)						
55 CAMBRIDGE PARKWAY, SUITE 800E								
(Street) CAMBRIDGE	MA	02142						
(City)	(State)	(Zip)						

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. Held by Flagship Ventures Fund 2004, L.P. Flagship Ventures General Partner LLC (the "GP") is the General Partner of Flagship Ventures Fund 2004, L.P. Noubar B. Afeyan, Ph.D. and Edwin M. Kania, Jr. are the Managers of the GP. The GP, Dr. Afeyan and Mr. Kania may be deemed to possess voting and investment control over all shares held by Flagship Ventures Fund 2004, L.P. Each of the reporting persons disclaims beneficial ownership of the shares except to the extent of his or its pecuniary interest therein.

3. 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. 4. Not applicable.

> FLAGSHIP VENTURES FUND 2004, L.P. By: Flagship Ventures General Partner LLC By: /s/ 12/11/2017 Noubar B. Afeyan, Ph.D., <u>Manager</u> **FLAGSHIP VENTURES** GENERAL PARTNER LLC By: 12/11/2017 /s/ Noubar B. Afeyan, Ph.D., Manager /s/ Noubar B. Afeyan, Ph.D. 12/11/2017 /s/ Edwin M. Kania, Jr. 12/11/2017 ** 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.