FORM 3

### UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

# INITIAL STATEMENT OF BENEFICIAL OWNERSHIP OF SECURITIES

OMB APPROVAL							
OMB Number: 3235-							
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*     bioMerieux/ADR			2. Date of Ever Requiring State (Month/Day/Ye 12/06/2017	ement	3. Issuer Name and Ticker or Trading Symbol Quanterix Corp [ QTRX ]							
(Last) CHEMIN DE	(First) L'ORME	(Middle)			4. Relationship of Reporting Pers (Check all applicable)  Director X	. ,	(Mon	Amendment, Da th/Day/Year)	te of Original Filed			
(Street) MARCY- L'ETOILE	IO (State)	69280			Officer (give title below)	Other (spec below)		cable Line) Form filed by	Group Filing (Check  One Reporting Person  More than One erson			
(City) (State) (Zip)												
Table I - Non-Derivative Securities Beneficially Owned												
				Amount of Securities     Beneficially Owned (Instr. 4)			Nature of Indirect Beneficial Ownership nstr. 5)					
Table II - Derivative Securities Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities)												
1. Title of Derivative Security (Instr. 4)  2. Date Exercisable and Expiration Date (Month/Day/Year)			ate	Underlying Derivative Security (Instr. 4)		4. Conversion or Exercise Price of	5. Ownership Form: Direct (D)	6. Nature of Indirect Beneficial Ownership (Instr. 5)				
			Date Exercisable	Expiration Date	Title	Amount or Number of Shares	Derivative Security	or Indirect (I) (Instr. 5)				
Series C Preferred Stock		(1)	(2)	Common Stock	2,055,632	(3)	D					

#### **Explanation of Responses:**

- 1. These shares are 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.
- 2. Not applicable.
- 3. The shares of preferred stock will convert 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 will convert on a 1-for-3.214 basis and such conversion rate is reflected in the amount of common stock underlying the security.

## Remarks:

Exhibit 24.1 - Power of Attorney

/s/ Brian Keane, Attorney-in-Fact for bioMerieux S.A. 12/06/2017

\*\* 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 5 (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.

#### **POWER OF ATTORNEY**

Know all by these presents, that the undersigned hereby constitutes and appoints Joseph Driscoll, Chief Financial Officer of Quanterix Corporation (the "Company"), Ernest Orticerio, Senior Vice President, Finance and Corporate Development of the Company, Brian Keane, General Counsel of the Company, and Megan Gates, John Condon, Brian Shea and Allyson Wilkinson, each of Mintz, Levin, Cohn, Ferris, Glovsky and Popeo, P.C., signing singly, with full power of substitution, the undersigned's true and lawful attorney-in-fact to:

- (1) execute for and on behalf of the undersigned, forms and authentication documents for EDGAR Filing Access;
- (2) do and perform any and all acts for and on behalf of the undersigned which may be necessary or desirable to complete and execute any such forms and authentication documents;
- (3) execute for and on behalf of the undersigned, in the undersigned's capacity as an officer, director and/or 10% shareholder of the Company, Forms 3, 4 and 5 in accordance with Section 16(a) of the Securities Exchange Act of 1934 and the rules thereunder;
- (4) do and perform any and all acts for and on behalf of the undersigned which may be necessary or desirable to complete and execute any such Form 3, 4 or 5 and timely file such form with the United States Securities and Exchange Commission and any stock exchange or similar authority; and
- take any other action of any type whatsoever in connection with the foregoing which, in the opinion of such attorney-in-fact, may be of benefit to, in the best interest of, or legally required by, the undersigned, it being understood that the documents executed by such attorney-in-fact, on behalf of the undersigned pursuant to this Power of Attorney, shall be in such form and shall contain such terms and conditions as such attorney-in-fact may approve in such attorney-in-fact's discretion.

The undersigned hereby grants to each such attorney-in-fact full power and authority to do and perform any and every act and thing whatsoever requisite, necessary, or proper to be done in the exercise of any of the rights and powers herein granted, as fully to all intents and purposes as the undersigned might or could do if personally present, with full power of substitution or revocation, hereby ratifying and confirming all that such attorney-in-fact, or such attorney-in-fact's substitute or substitutes, shall lawfully do or cause to be done by virtue of this power of attorney and the rights and powers herein granted. The undersigned acknowledges that the foregoing attorneys-in-fact, in serving in such capacity at the request of the undersigned, is not assuming, nor is the Company assuming, any of the undersigned's responsibilities to comply with Section 16 of the Securities Exchange Act of 1934.

This Power of Attorney shall remain in full force and effect until the undersigned is no longer required to file Forms 3, 4 and 5 with respect to the undersigned's holdings of and transactions in securities issued by the Company, unless earlier revoked by the undersigned in a signed writing delivered to the foregoing attorneys-in-fact.

IN WITNESS WHEREOF, the undersigned has caused this Power of Attorney to be executed this 1st day of December 2017.

## BIOMÉRIEUX S.A.

By: /s/ Alexandre Mérieux
Name: Alexandre Mérieux