
**UNITED STATES
SECURITIES AND EXCHANGE COMMISSION
WASHINGTON, D.C. 20549**

SCHEDULE 13G

**Under the Securities Exchange Act of 1934
(Amendment No.)***

Quanterix Corporation
(Name of Issuer)

Common Stock
(Title of Class of Securities)

74766Q101
(CUSIP Number)

December 31, 2017
(Date of Event Which Requires Filing of this Statement)

Check the appropriate box to designate the rule pursuant to which this Schedule is filed:

- Rule 13d-1(b)
- Rule 13d-1(c)
- Rule 13d-1(d)

* The remainder of this cover page shall be filled out for a reporting person's initial filing on this form with respect to the subject class of securities, and for any subsequent amendment containing information which would alter disclosures provided in a prior cover page.

The information required in the remainder of this cover page shall not be deemed to be "filed" for the purpose of Section 18 of the Securities Exchange Act of 1934 ("Act") or otherwise subject to the liabilities of that section of the Act, but shall be subject to all other provisions of the Act (however, see the Notes).

1	NAMES OF REPORTING PERSONS	
	bioMérieux SA	
2	CHECK THE APPROPRIATE BOX IF A MEMBER OF A GROUP* (a) <input type="checkbox"/> (b) <input type="checkbox"/>	
3	SEC USE ONLY	
4	CITIZENSHIP OR PLACE OF ORGANIZATION	
	France	
NUMBER OF SHARES BENEFICIALLY OWNED BY EACH REPORTING PERSON WITH	5	SOLE VOTING POWER
		2,055,632
	6	SHARED VOTING POWER
		0
	7	SOLE DISPOSITIVE POWER
		2,055,632
	8	SHARED DISPOSITIVE POWER
		0
9	AGGREGATE AMOUNT BENEFICIALLY OWNED BY EACH REPORTING PERSON	
	2,055,632	
10	CHECK IF THE AGGREGATE AMOUNT IN ROW (9) EXCLUDES CERTAIN SHARES* <input type="checkbox"/>	
11	PERCENT OF CLASS REPRESENTED BY AMOUNT IN ROW (9)	
	9.4%	
12	TYPE OF REPORTING PERSON*	
	CO	

-
- Item 1(a). Name of Issuer**
Quanterix Corporation (the “Issuer”).
- Item 1(b). Address of Issuer’s Principal Executive Offices**
113 Hartwell Avenue, Lexington, MA 02421
- Item 2(a). Name of Person Filing**
bioMérieux SA
- Item 2(b). Address of Principal Business Office or, if none, Residence**
376 Chemin de l’Orme, 69280 Marcy l’Étoile, France
- Item 2(c). Citizenship**
France
- Item 2(d). Title of Class of Securities**
Common stock, par value \$0.001 per share.
- Item 2(e). CUSIP Number**
74766Q101.
- Item 3. If This Statement Is Filed Pursuant to Sections 240.13d-1(b) or 240.13d-2(b) or (c), Check Whether the Person Filing Is a(n):**
Not Applicable.
- Item 4. Ownership**
- (a) Amount beneficially owned:
2,055,632 shares of Common Stock.
- (b) Percent of class:
bioMérieux SA beneficially owns approximately 9.4% of the outstanding shares of common stock of the Issuer, based upon 21,820,840 shares of common stock outstanding as of December 11, 2017, which combines 21,179,560 Common Stock outstanding as reported on the Issuer’s 424B4

as filed with the Securities and Exchange Commission on December 7, 2017, plus 641,280 shares issued (underwriter's allocation) as reported on the Issuer's 8-K as filed with the Securities and Exchange Commission on December 15, 2017.

- (c) Number of shares as to which such person has:
 - (i) Sole power to vote or to direct the vote:
2,055,632
 - (ii) Shared power to vote or to direct the vote:
0
 - (iii) Sole power to dispose or to direct the disposition:
2,055,632
 - (iv) Shared power to dispose or to direct the disposition:
0

Item 5. Ownership of Five Percent or Less of a Class

Not Applicable.

Item 6. Ownership of More Than Five Percent on Behalf of Another Person

Not Applicable.

Item 7. Identification and Classification of the Subsidiary Which Acquired the Security Being Reported on By the Parent Holding Company

Not Applicable.

Item 8. Identification and Classification of Members of the Group

Not Applicable.

Item 9. Notice of Dissolution of Group

Not Applicable.

Item 10. Certifications

Not Applicable.

SIGNATURE

After reasonable inquiry and to the best of my knowledge and belief, I certify that the information set forth in this statement is true, complete and correct.

Dated: February 14, 2018

BIOMÉRIEUX SA

By /s/ Andy Parker

Name: Andy Parker

Title: Vice President - Human Resources & General Counsel
- Americas