## UNITED STATES SECURITIES AND EXCHANGE COMMISSION

WASHINGTON, D.C. 20549

S	CI	$\Pi$	${f F.D}$	T	JT.	E.	13	$\mathbf{G}$	Ά
$\mathbf{\mathcal{O}}$	$\sim$		-	$\sim$	_	_		$\mathbf{u}$	

Under the Securities Exchange Act of 1934 (Amendment No. 1 – Exit Filing)\*

## **Quanterix Corporation**

(Name of Issuer)

Common Stock (Title of Class of Securities)

> 74766Q101 (CUSIP Number)

November 7, 2019 (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).

CUSI	IP No. 74766Q1	101				
1	1 NAMES OF REPORTING PERSONS					
	bioMérieux SA					
2	2 CHECK THE APPROPRIATE BOX IF A MEMBER OF A GROUP*					
	(a) $\Box$ (b) $\Box$					
3	SEC USE ON	LY				
4	CITIZENSHI	P (	OR PLACE OF ORGANIZATION			
	France					
		5	SOLE VOTING POWER			
N	UMBER OF		0			
	SHARES NEFICIALLY	6	SHARED VOTING POWER			
	OWNED BY		0			
R	EACH EPORTING	7	SOLE DISPOSITIVE POWER			
	PERSON WITH		0			
	WIIH	8	SHARED DISPOSITIVE POWER			
			0			
9	AGGREGAT	E A	MOUNT BENEFICIALLY OWNED BY EACH REPORTING PERSON			
	0					
10	10 CHECK IF THE AGGREGATE AMOUNT IN ROW (9) EXCLUDES CERTAIN SHARES*					
11	11 PERCENT OF CLASS REPRESENTED BY AMOUNT IN ROW (9)					
	0%					
12	12 TYPE OF REPORTING PERSON*					
	СО					

Item 1(a).	Name of Issuer						
	Quanterix Corporation (the "Issuer").						
Item 1(b).	Address of Issuer's Principal Executive Offices						
	900 Middlesex Turnpike Billerica, MA 01821						
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:						
	0 shares of Common Stock.						
	(b) Percent of class:						

	(c) Number of shares as to which such person has:						
		(i)	Sole power to vote or to direct the vote:				
			0				
		(ii)	Shared power to vote or to direct the vote:				
			0				
		(iii)	Sole power to dispose or to direct the disposition:				
			0				
		(iv)	Shared power to dispose or to direct the disposition:				
			0				
Item 5.	Ownership of Five Percent or Less of a Class						
			ent is being filed to report the fact that as of the date hereof the reporting person has ceased to be the beneficial owner of more at of the class of securities, check the following:				
Item 6.	Ownership of More Than Five Percent on Behalf of Another Person						
	Not A	Applicat	ole.				
Item 7.	Identification and Classification of the Subsidiary Which Acquired the Security Being Reported on By the Parent Holding Company						
	Not A	Applicat	lle.				
Item 8.	Identification and Classification of Members of the Group						
		Applicat					
		••					
Item 9.	Notic	e of Dis	ssolution of Group				
	Not A	Applicat	ole.				

**Item 10.** 

**Certifications**Not Applicable.

## $\underline{SIGNATURE}$

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: November 18, 2019

## BIOMÉRIEUX SA

By /s/ Alexandre Mérieux

Name: Alexandre Mérieux Title: Chief Executive Officer