UNITED STATES SECURITIES AND EXCHANGE COMMISSION Washington, D.C. 20549

SCHEDULE 13G

INFORMATION TO BE INCLUDED IN STATEMENTS FILED PURSUANT TO RULES 13d -1(b), (c), AND (d) AND AMENDMENTS THERETO FILED PURSUANT TO RULE 13d-2. (Amendment No.)*

	Quanterix Corporation
	(Name of Issuer)
	Common Stock
	(Title of Class of Securities)
	74766Q101 (CUSIP Number)
	June 17, 2024
	(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)
	inder 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 quent amendment containing information which would alter the disclosures provided in a prior cover page.
	nation 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 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

CU	USIP N	No. 74766	6Q10	1	
	-				
	1.	. Names of Reporting Persons Portolan Capital Management, LLC			
	•				
	2.	Check tl	he Ap	propriate Box if a Member of a Group (See Instructions)	
	(a) Control in the properties Both in a strong (see instance of a strong of a				
(b)					
	3.	SEC Use	e Onl	у	
	4	Citizens Delawar		r Place of Organization	
	l.				
			5.	Sole Voting Power	
				1,918,030	
Νι	umber	of			
	nares		6.	Shared Voting Power	
	enefici wned b				
Еа	wheu t ach eportin		7.	Sole Dispositive Power	
	erson V			1,918,030	
			8.	Shared Dispositive Power	
	9.	Aggregate Amount Beneficially Owned by Each Reporting Person			
		1,918,03	30		
	10.	Check if	the A	Aggregate Amount in Row (9) Excludes Certain Shares (See Instructions) □	
	11.	Percent of 5.01%	of Cla	ass Represented by Amount in Row (9)	
	•				
	12.	Type of IA	Repo	rting Person (See Instructions)	

С	USIP N	No. 7476	6Q10	01	
	1.	Names of Reporting Persons George McCabe			
	•	•			
	2.	Check tl	ne Ap	propriate Box if a Member of a Group (See Instructions)	
		(a) 🗆			
		(b)			
	3. SEC Use Only			у	
	4 Citizenship or Place of Organization USA				
		5.	Sole Voting Power		
				1,918,030	
N	umber	of			
S	hares		6.	Shared Voting Power	
	enefici wned b				
E	ach		7.	Sole Dispositive Power	
Po	Reporting Person With			1,918,030	
Γ					
			8.	Shared Dispositive Power	
			1	•	
	9.	Aggregate Amount Beneficially Owned by Each Reporting Person 1,918,030			
	10.	Check if	the A	Aggregate Amount in Row (9) Excludes Certain Shares (See Instructions) □	
	11.	Percent 5.01%	of Cla	ass Represented by Amount in Row (9)	
	•				
	12.	Type of IN	Repo	rting Person (See Instructions)	

Item 1.					
	(a)	Name of Issuer Quanterix Corporation			
	(b)	900 Middlesex Turnpike Billerica, Massachusetts 01821			
Item 2.					
	(a)	Name of Person Filing This statement is being filed with respect to the shares of class A common stock ("Common Stock") of the Issuer beneficially owned (1) directly by Portolan Capital Management, LLC, a registered investment adviser, in its capacity as investment manager for various clients, and (2) indirectly by George McCabe, the Manager of Portolan Capital Management, LLC. Portolan Capital Management, LL and Mr. McCabe are sometimes individually referred to herein as a "Reporting Person" and collectively as the "Reporting Persons."			
	(b)	Address of Principal Business Office or, if none, Residence Portolan Capital Management, LLC and George McCabe 2 International Place, FL 26, Boston, MA 02110			
(c) Citizenship Portolan Capital Management, LLC – DE Mr. McCabe – USA					
	(d)	of Class of Securities amon Stock			
	(e)	CUSIP Number 74766Q101			
Item 3.	If th	is stat	ement is filed pursuant to §§240.13d-1(b) or 240.13d-2(b) or (c), check whether the person filing is a:		
	(a)		Broker or dealer registered under section 15 of the Act (15 U.S.C. 78o).		
	(b)		Bank as defined in section 3(a)(6) of the Act (15 U.S.C. 78c).		
	(c)		Insurance company as defined in section 3(a)(19) of the Act (15 U.S.C. 78c).		
	(d)		Investment company registered under section 8 of the Investment Company Act of 1940 (15 U.S.C. 80a-8).		
	(e)		An investment adviser in accordance with §240.13d-1(b)(1)(ii)(E);		
	(f)		An employee benefit plan or endowment fund in accordance with §240.13d-1(b)(1)(ii)(F);		
	(g)		A parent holding company or control person in accordance with §240.13d-1(b)(1)(ii)(G);		
	(h)		A savings association as defined in Section 3(b) of the Federal Deposit Insurance Act (12 U.S.C. 1813);		
	(i)		A church plan that is excluded from the definition of an investment company under section 3(c)(14) of the Investment Company Act of 1940 (15 U.S.C. 80a-3);		
	(j)		A non-U.S. institution in accordance with § 240.13d–1(b)(1)(ii)(J);		
	(k)		Group, in accordance with §240.13d-1(b)(1)(ii)(J).		
			3		

T. 4	
Item 4.	Ownership

Provide the following information regarding the aggregate number and percentage of the class of securities of the issuer identified in Item 1.

Reference is hereby made to Items 5-9 and 11 of pages 1 - 2 of this Schedule, which Items are incorporated by reference herein.

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

If this statement 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 than five percent of the class of securities, check the following \Box .

Not applicable.

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

Various persons have the right to receive or the power to direct the receipt of dividends from, or the proceeds from the sale of, the Common Stock of Quanterix Corporation. No one person's interest in the Common Stock of Quanterix Corporation is more than five percent of the total outstanding Common Stock.

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

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. Certification

By signing below I certify that, to the best of my knowledge and belief, the securities referred to above were not acquired and are not held for the purpose of or with the effect of changing or influencing the control of the issuer of the securities and were not acquired and are not held in connection with or as a participant in any transaction having that purpose or effect.

Signature

	June 27, 2024
	Date
	Portolan Capital Management, LLC
By:	/s/ George McCabe
-	George McCabe, Manager
	/s/ George McCabe
-	George McCabe

Intentional misstatements or omissions of fact constitute Federal Criminal Violations (See 18 U.S.C. 1001).