

SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

SCHEDULE 13G

UNDER THE SECURITIES EXCHANGE ACT OF 1934

(Amendment No. 5)*

Quanterix Corp

(Name of Issuer)

Common Stock, \$0.001 par value per share

(Title of Class of Securities)

(CUSIP Number)

06/30/2026

(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)

SCHEDULE 13G

CUSIP No.

- Names of Reporting Persons
- 1Portolan Capital Management, LLC
- Check the appropriate box if a member of a Group (see instructions)
- 2☐ (a)
- ☐ (b)
- 3Sec Use Only
- Citizenship or Place of Organization
- 4DELAWARE

Number of Shares Beneficially Owned by Each Reporting Person With:	5	Sole Voting Power
		1,899,308.00
		Shared Voting Power
	6	0.00
		Sole Dispositive Power
	7	1,899,308.00
		Shared Dispositive Power
	8	0.00
9	Aggregate Amount Beneficially Owned by Each Reporting Person	
		1,899,308.00
10	Check box if the aggregate amount in row (9) excludes certain shares (See Instructions)	
		☐
11	Percent of class represented by amount in row (9)	
		4.02 %
12	Type of Reporting Person (See Instructions)	
		IA

SCHEDULE 13G

CUSIP No.

1	Names of Reporting Persons	
	George McCabe	
	Check the appropriate box if a member of a Group (see instructions)	
2	☐	(a)
	☐	(b)
3	Sec Use Only	
4	Citizenship or Place of Organization	
	UNITED STATES	
		Sole Voting Power
	5	1,899,308.00
		Shared Voting Power
	6	0.00
		Sole Dispositive Power
	7	1,899,308.00
		Shared Dispositive Power
	8	0.00
9	Aggregate Amount Beneficially Owned by Each Reporting Person	
		1,899,308.00
10	Check box if the aggregate amount in row (9) excludes certain shares (See Instructions)	
		☐

11	Percent of class represented by amount in row (9)
	4.02 %
12	Type of Reporting Person (See Instructions)
	IN

SCHEDULE 13G

Item 1.

Name of issuer:

(a)

Quanterix Corp

Address of issuer's principal executive offices:

(b)

900 Middlesex Turnpike, Billerica, MA 01821

Item 2.

Name of person filing:

(a)

This statement is being filed with respect to the shares 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, LLC and Mr. McCabe are sometimes individually referred to herein as a "Reporting Person" and collectively as the "Reporting Persons."

Address or principal business office or, if none, residence:

(b)

The address of the principal business office of each of the Reporting Persons is 2 International Place, FL 26, Boston, MA 02110

Citizenship:

(c)

Portolan Capital Management, LLC - DE Mr. McCabe - USA

Title of class of securities:

(d)

Common Stock, \$0.001 par value per share

(e)

CUSIP No.:

Item 3. If this statement 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 associations 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). If filing as a non-U.S. institution in accordance with § 240.13d-1(b)(1)(ii)(J), please specify the type of institution:

(k)

☐ Group, in accordance with Rule 240.13d-1(b)(1)(ii)(K).

Item 4. Ownership

Amount beneficially owned:

(a)

The information required by Item 4(a) is set forth in Row 9 of the cover page for each of the Reporting Persons and is incorporated by reference herein.

(b)

Percent of class:

The information required by Item 4(b) is set forth in Row 11 of the cover page for each of the Reporting Persons and is incorporated by reference herein. %

(c) Number of shares as to which the person has:

(i) Sole power to vote or to direct the vote:

The information required by Item 4(c)(i) is set forth in Row 5 of the cover page for each of the Reporting Persons and is incorporated by reference herein.

(ii) Shared power to vote or to direct the vote:

0

(iii) Sole power to dispose or to direct the disposition of:

The information required by Item 4(c)(iii) is set forth in Row 7 of the cover page for each of the Reporting Persons and is incorporated by reference herein.

(iv) Shared power to dispose or to direct the disposition of:

0

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

☒ Ownership of 5 percent or less of a class

Item 6. Ownership of more than 5 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 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. Certifications:

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, other than activities solely in connection with a nomination under § 240.14a-11.

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.

Portolan Capital Management, LLC

Signature: George McCabe

Name/Title: George McCabe/Manager

Date: 08/14/2026

George McCabe

Signature: George McCabe

Name/Title: George McCabe

Date: 08/14/2026