

SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

SCHEDULE 13G

UNDER THE SECURITIES EXCHANGE ACT OF 1934

Quanterix Corp

(Name of Issuer)

Common

(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	
1	First Eagle Investment Management, LLC	
	Check the appropriate box if a member of a Group (see instructions)	
2	☐ (a)	
	☐ (b)	
3	Sec Use Only	
	Citizenship or Place of Organization	
4	DELAWARE	
Number of	Sole Voting Power	
Shares	5	
Beneficially	2,753,165.00	

Owned by Each Reporting Person With:	6	Shared Voting Power
		0.00
		Sole Dispositive Power
	7	2,753,165.00
		Shared Dispositive Power
	8	0.00
9	Aggregate Amount Beneficially Owned by Each Reporting Person	
		2,753,165.00
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)	
		5.8 %
12	Type of Reporting Person (See Instructions)	
		IA

SCHEDULE 13G

CUSIP No.

1	Names of Reporting Persons	
	FIRST EAGLE SMALL CAP OPPORTUNITY FUND	
	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	
	DELAWARE	
	Sole Voting Power	
	5	2,421,533.00
Number of Shares Beneficially Owned by Each Reporting Person With:	6	Shared Voting Power
		0.00
		Sole Dispositive Power
	7	2,421,533.00
		Shared Dispositive Power
	8	0.00
9	Aggregate Amount Beneficially Owned by Each Reporting Person	
		2,421,533.00
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)	
		5.1 %

IC

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) First Eagle Investment Management, LLC, FIRST EAGLE SMALL CAP OPPORTUNITY FUND
Address or principal business office or, if none, residence:
- (b) 1345 Avenue of the Americas, New York, NY 10105, United States
Citizenship:
- (c) DELAWARE
Title of class of securities:
- (d) Common
CUSIP No.:
- (e)

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) 2,753,165
Percent of class:
- (b) 5.8 %
- (c) Number of shares as to which the person has:
- (i) Sole power to vote or to direct the vote:
- 2,753,165
- (ii) Shared power to vote or to direct the vote:

0

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

2,753,165

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

0

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

Item 6. Ownership of more than 5 Percent on Behalf of Another Person.

If any other person is known to have the right to receive or the power to direct the receipt of dividends from, or the proceeds from the sale of, such securities, a statement to that effect should be included in response to this item and, if such interest relates to more than 5 percent of the class, such person should be identified. A listing of the shareholders of an investment company registered under the Investment Company Act of 1940 or the beneficiaries of employee benefit plan, pension fund or endowment fund is not required.

First Eagle Investment Management, LLC (FEIM), an investment adviser registered under Section 203 of the Investment Advisers Act of 1940, is deemed to be the beneficial owner of 2,753,165 shares, or 5.84% of the common stock believed to be outstanding as a result of acting as investment adviser to various clients. Clients of FEIM have the right to receive and the ultimate power to direct the receipt of dividends from, or the proceeds of the sale of, such securities. The First Eagle Small Cap Opportunity Fund, a registered investment company for which FEIM acts as investment adviser, may be deemed to beneficially own 2,421,533 of these 2,753,165 shares, or 5.14% of the Company's 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. Certifications:

By signing below I certify that, to the best of my knowledge and belief, the securities referred to above were acquired and are held in the ordinary course of business and 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.

First Eagle Investment Management, LLC

Signature: David O'Connor

Name/Title: General Counsel

Date: 08/06/2026

FIRST EAGLE SMALL CAP OPPORTUNITY FUND

Signature: David O'Connor

Name/Title: General Counsel

Date: 08/06/2026