UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

SCHEDULE 13G

Under the Securities Exchange Act of 1934 (Amendment No. 1)*

Quanterix Corp

(Name of Issuer)

Common Shares (Title of Class of Securities)

74766Q101 (CUSIP Number)

December 30, 2022 (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 the 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).

CUSIP No. 74766Q101

1.	. Names of Reporting Persons				
	Macquarie Group Limited				
2.	Check	the A	Appropriate Box if a Member of a Group (See Instructions)		
	$(a) \boxtimes (b) \square$				
3.	SEC U	se O	nly		
4.	Citizer	shin	or Place of Organization		
т.	4. Cluzenship of Place of Organization				
	Sydney	, Ne	w South Wales Australia		
		5.	Sole Voting Power		
Nur	nber of		0		
	hares	6.	Shared Voting Power		
Beneficially					
Owned by Each		7.	0 Sole Dispositive Power		
Reporting		,.	Sole Dispositive Former		
	erson		0		
With		8.	Shared Dispositive Power		
			0		
9.	Aggreg	gate 1	Amount Beneficially Owned by Each Reporting Person		
	0 deer	ad h	eneficially owned due to reporting person's ownership of Macquarie Investment Management Holdings Inc., Macquarie		
			Management Business Trust, whose individual holdings are shown on the following forms.		
10.	Check	if th	e Aggregate Amount in Row (9) Excludes Certain Shares (See Instructions)		
	П				
11.	_	t of (Class Represented by Amount in Row (9)		
10	0.00%	CD			
12.	Type o	i Kej	porting Person (See Instructions)		
	НС				

CUSIP No. 74766Q101

1	M	a C D	an arting Danson		
1.	1. Names of Reporting Persons				
	Macquarie Management Holdings Inc				
2.	2. Check the Appropriate Box if a Member of a Group (See Instructions)				
	(a) \boxtimes (b) \square				
3.	SEC U	se O	nly		
4.	4. Citizenship or Place of Organization				
١.	4. Citizenship of Trace of Organization				
State of Delaware					
Number of Shares Beneficially Owned by Each		5.	Sole Voting Power		
			0		
		6.	Shared Voting Power		
			0		
		7.	Sole Dispositive Power		
	oorting erson		0		
	With	8.	Shared Dispositive Power		
		0.	Shared Dispositive Fower		
			0		
9.	Aggreg	gate 1	Amount Beneficially Owned by Each Reporting Person		
	0 deem	ned b	eneficially owned due to reporting person's ownership of Macquarie Investment Management Business Trust		
10.	1 1 1 1 1 1 1				
	\boxtimes				
11.		t of (Class Represented by Amount in Row (9)		
	0.00%				
12.	Type of Reporting Person (See Instructions)				
	НС				

CUSIP No. 74766Q101

1.	1. Names of Reporting Persons				
	Macquarie Investment Management Business Trust				
2.	2. Check the Appropriate Box if a Member of a Group (See Instructions)				
	(a) ⊠ (b) □				
3.	3. SEC Use Only				
4.	4. Citizenship or Place of Organization				
	State of Delaware				
Number of Shares Beneficially Owned by Each Reporting Person		5.	Sole Voting Power		
			0		
		6.	Shared Voting Power		
			0		
		7.	Sole Dispositive Power		
			0		
,	With	8.	Shared Dispositive Power		
			0		
9.	Aggreg	gate 1	Amount Beneficially Owned by Each Reporting Person		
	0				
10.	0. Check if the Aggregate Amount in Row (9) Excludes Certain Shares (See Instructions)				
11.	Percen	t of (Class Represented by Amount in Row (9)		
	0.00%				
12.					
	IA				

Item 1.		
	(a)	Name of Issuer Quanterix Corp
	(b)	Address of Issuer's Principal Executive Offices 900 Middlesex Turnpike, Building 1 Billerica, MA 01821
Item 2.		
	(a)	Name of Person Filing This Schedule 13G is jointly filed by Macquarie Group Limited, Macquarie Management Holdings Inc, Macquarie Investment Management Business Trust
	(b)	Address of Principal Business Office or, if none, Residence The principal business address of Macquarie Group Limited is 50 Martin Place Sydney, New South Wales, Australia. The principal business address of Macquarie Management Holdings Inc. and Macquarie Investment Management Business Trust is 2005 Market Street, Philadelphia, PA 19103.
	(c)	Citizenship Macquarie Group Limited, Macquarie Investment Management Global Limited - Sydney, New South Wales, Australia Corporation Macquarie Management Holdings Inc., Macquarie Investment Management Business Trust– incorporated or formed under the laws of the State of Delaware.
	(d)	Title of Class of Securities Common Stock
	(e)	CUSIP Number 74766Q101
Item 3.	If this s	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. 780);
	(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 $\S 240.13d-1(b)(1)(ii)(K)$. If filing as a non-U.S. institution in accordance with $\S 240.13d-1(b)(1)(ii)(J)$, please specify the type of institution:
		Page 5 of 12

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.

- (a) Amount beneficially owned: See responses on the cover page hereto.
- (b) Percent of class:

See responses on the cover page hereto.

- (c) Number of shares as to which the person has:
 - (i) Sole power to vote or to direct the vote See responses on the cover page hereto.
 - (ii) Shared power to vote or to direct the vote
 - (iii) Sole power to dispose or to direct the disposition of See responses on the cover page hereto.
 - (iv) Shared power to dispose or to direct the disposition of

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 \boxtimes

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 or Control Person

See Exhibit A.

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 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.

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.

Macquarie Group Limited	February 14, 2023
•	Date
/s/ Michael Wainer	/s/ Charles Glorioso
Signature	Signature
Michael Wainer	Charles Glorioso
Associate Director	Division Director
After reasonable inquiry and to the best of my knowledge and belief, I certificorrect.	y that the information set forth in this statement is true, complete and
Macquarie Management Holdings, Inc.	February 14, 2023
	Date
/s/ Brian L. Murray	
Signature	_
Brian L. Murray Chief Compliance Officer	_
Macquarie Investment Management Business Trust	February 14, 2023 Date
/s/ Brian L. Murray	
Signature	_
Brian L. Murray Chief Compliance Officer	_

Page 7 of 12

EXHIBIT A

AGREEMENT TO FILE JOINT ACQUISITION STATEMENTS

AGREEMENT made this 2nd day of FEBRUARY, 2021 by and between Delaware Funds® by Macquarie listed on Annex A hereto, Macquarie Investment Management Business Trust, Macquarie Management Holdings, Inc, and the Macquarie Parties listed on Annex B hereto (collectively referred to as the "parties").

WHEREAS, the parties hereto may be deemed to be the direct or indirect beneficial owners of the same equity securities for the purpose of the reporting requirements of Section 13(d) of the Securities Exchange Act of 1934, as amended (the "Exchange Act"), and

WHEREAS, the regulations promulgated under Section 13(d) of the Exchange Act permit the joining of such beneficial owners in the filing of a single Joint Acquisition Statement reporting such ownership to the Securities and Exchange Commission.

NOW, THEREFORE, in consideration of the mutual covenants herein contained, and each of the parties hereto intending to be legally bound, it is agreed as follows:

- 1. In the event that any two or more parties shall be deemed to be the direct or indirect beneficial owners of the same equity security required to be reported to the Securities and Exchange Commission such parties may join together in the filing of a Joint Acquisition Statement with respect to that security. Additional persons who may after the date hereof be deemed to be the direct or indirect beneficial owners of the same equity security as a party hereto and required to be reported to the Securities and Exchange Commission (a "New Party") may be added as a party this agreement by signing a counterpart hereof. An amendment to this agreement is deemed effective upon the signature of such new party and the amendment of the applicable Annex which may be affixed to this agreement as amended. Each party hereto agrees that this agreement, as it may be amended from time to time as provided herein, is a valid and binding agreement of each such party.
- 2. With respect to each Joint Acquisition Statement in which a party joins, each party acknowledges that (a) it will be eligible under applicable regulations of the Securities and Exchange Commission to join in the filing and (b) it will be responsible for the timely filing of such statement and any amendments thereto and the completeness and accuracy of the information concerning such party; but each such party shall not be responsible for the completeness and accuracy of the information concerning the other parties making the filing, unless such party knows or has reason to believe that such information with respect to such other parties is inaccurate.
- 3. The parties consent to the inclusion of a copy of this agreement as an exhibit to any Joint Acquisition Statement filed on behalf of any of them.

IN WITNESS WITEREOF, the parties hereto have executed this agreement by their duty	authorized officers as of the date set forth above.	
DELAWARE FUNDS® BY MACQUARIE (listed on Annex A hereto)		
ATTEST BY:		
/s/ Brian L. Murray	/s/ David Connor	
Signature	Signature	
Brian L. Murray Chief Compliance Officer	David Connor General Counsel	
MACQUARIE INVESTMENT MANAGEMENT BUSINESS TRUST	Solicial Counsel	
/s/ Brian L. Murray	/s/ David Connor	
Signature	Signature	
Brian L. Murray Chief Compliance Officer	David Connor General Counsel	
MACQUARIE MANAGEMENT HOLDINGS, INC.		
/s/ Brian L. Murray	/s/ David Connor	
Signature	Signature	
Brian L. Murray Chief Compliance Officer	David Connor General Counsel	
THE MACQUARIE PARTIES (LISTED ON ANNEX B HERETO)		
ATTEST BY:		
/s/ Michael Wainer	/s/ Charles Glorioso	
Signature	Signature	
Michael Wainer Associate Director	Charles Glorioso Division Director	

Annex A— Delaware FundsSM by Macquarie

- DELAWARE GROUP EQUITY FUNDS I
- DELAWARE GROUP EQUITY FUNDS II
- DELAWARE GROUP EQUITY FUNDS IV
- DELAWARE GROUP EQUITY FUNDS V
- DELAWARE GROUP INCOME FUNDS
- DELAWARE GROUP LIMITED-TERM GOVERNMENT FUNDS
- DELAWARE GROUP CASH RESERVE
- DELAWARE GROUP GOVERNMENT FUND
- DELAWARE GROUP STATE TAX-FREE INCOME TRUST
- DELAWARE GROUP TAX-FREE FUND
- DELAWARE GROUP GLOBAL & INTERNATIONAL FUNDS
- DELAWARE GROUP ADVISER FUNDS
- DELAWARE VIP TRUST
- DELAWARE POOLED TRUST
- DELAWARE GROUP FOUNDATION FUNDS
- DELAWARE INVESTMENTS DIVIDEND AND INCOME FUND, INC.
- DELAWARE ENHANCED GLOBAL DIVIDEND AND INCOME FUND
- INVESTED PORTFOLIOS
- IVY HIGH INCOME OPPORTUNITIES FUND
- **IVY FUNDS**
- IVY VARIABLE INSURANCE PORTFOLIOSSM
- VOYAGEUR INSURED FUNDS
- VOYAGEUR INTERMEDIATE TAX FREE FUNDS
- VOYAGEUR MUTUAL FUNDS
- VOYAGEUR MUTUAL FUNDS II
- VOYAGEUR MUTUAL FUNDS III
- VOYAGEUR TAX FREE FUNDS
- DELAWARE INVESTMENTS COLORADO MUNICIPAL INCOME FUND, INC.
- DELAWARE INVESTMENTS NATIONAL MUNICIPAL INCOME FUND
- DELAWARE INVESTMENTS MINNESOTA MUNICIPAL INCOME FUND II, INC.

Page 10 of 12

Annex B — the Macquarie Parties

Macquarie Group Limited
Macquarie Bank Limited
Macquarie Affiliated Managers (USA) Inc.
Macquarie Affiliated Managers Holdings (USA) Inc.
Macquarie Americas Holdings Pty Ltd.
Macquarie B.H. Pty Limited
Macquarie FG Holdings Inc.
Macquarie Funding Holdings LLC
Macquarie Investment Management Europe Limited

Page 11 of 12

EXHIBIT B

Powers of Attorney for Macquarie Group Limited and Macquarie Bank Limited incorporated by reference to 13G filings made by Macquarie Group Limited and Macquarie Bank Limited on May 25, 2021.

Page 12 of 12