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

**SCHEDULE 13G Under the Securities Exchange Act of 1934** 

(AMENDMENT NO. 1)

	QUANTERIX CORPORATION	
	(Name of Issuer)	
	COMMON STOCK, PAR VALUE \$0.001 PER SHARE	
	(Title of Class of Securities)	
	74766Q101	
	(CUSIP Number)	
	<b>DECEMBER 31, 2023</b>	
	(Date of event which requires filing of this statement)	
neck 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)		

Check

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 1	No. 7476	6Q101	SCHEDULE 13G	Page 2 of 11				
1	NAMES OF REI							
2	CHECK THE APPROPRIATE BOX IF A MEMBER OF A GROUP							
3	SEC USE ONLY							
4	CITIZENSHIP C Delaware	OR PLACE OF	ORGANIZATION					
	HINADED OF	5 SC -0-	LE VOTING POWER					
BE	IUMBER OF SHARES INEFICIALLY OWNED BY	6	IARED VOTING POWER 982,903					
EACH REPORTING PERSON WITH		7 SC -0-	DLE DISPOSITIVE POWER					
		8	IARED DISPOSITIVE POWER 082,903					
9	AGGREGATE A 2,082,903	MOUNT BEN	EFICIALLY OWNED BY EACH REPORTING PERSON					
10	CHECK BOX IF	THE AGGRE	GATE AMOUNT IN ROW (9) EXCLUDES CERTAIN SHARES					

PERCENT OF CLASS REPRESENTED BY AMOUNT IN ROW (9)

5.5%

00

12

TYPE OF REPORTING PERSON

CUSIP N	No. 74766Q101		SCHEDULE 13G	Page	3	of	11				
	NAMES OF REPORTING P	ERSONS	S								
1	Millennium Management LL	C									
			IF A MEMBER OF A GROUP								
2	(a) □	. L 2011	ii miliibliko: monoo:								
	(b)										
3	SEC USE ONLY										
	CITIZENSHIP OR PLACE OF ORGANIZATION										
4	Deleviore										
	Delaware										
		_	SOLE VOTING POWER								
		5	-0-								
	NUMBER OF		SHARED VOTING POWER								
	SHARES BENEFICIALLY	6									
	OWNED BY		2,209,566								
EACH		_	SOLE DISPOSITIVE POWER								
	REPORTING	7	-0-								
	PERSON WITH		SHARED DISPOSITIVE POWER								
		8	SILIKED DISTOSITIVE TOWER								
			2,209,566								

AGGREGATE AMOUNT BENEFICIALLY OWNED BY EACH REPORTING PERSON

PERCENT OF CLASS REPRESENTED BY AMOUNT IN ROW (9)

CHECK BOX IF THE AGGREGATE AMOUNT IN ROW (9) EXCLUDES CERTAIN SHARES

2,209,566

5.8%

00

TYPE OF REPORTING PERSON

10

11

12

CUSIP	No. 74766Q101		SCHEDULE 13G	Page 4	of 11			
.	NAMES OF REPORTING	G PERSON	IS .					
1	Millennium Group Mana	gement LLC						
	CHECK THE APPROPR	IATE BOX	IF A MEMBER OF A GROUP					
2	(a)							
	(b) 🗆							
3	SEC USE ONLY							
<b>l</b> ,	CITIZENSHIP OR PLAC	CE OF ORC	GANIZATION					
4	Delaware							
			SOLE VOTING POWER -0-					
	NUMBER OF SHARES BENEFICIALLY OWNED BY	6	SHARED VOTING POWER 2,209,566					
	EACH REPORTING PERSON WITH		SOLE DISPOSITIVE POWER -0-					
		8	SHARED DISPOSITIVE POWER 2,209,566					
	AGGREGATE AMOUN	Γ BENEFIC	CIALLY OWNED BY EACH REPORTING PERSON					
9								
	2,209,566							
	CHECK BOX IF THE AGGREGATE AMOUNT IN ROW (9) EXCLUDES CERTAIN SHARES							

PERCENT OF CLASS REPRESENTED BY AMOUNT IN ROW (9)

TYPE OF REPORTING PERSON

CUSIP N	No. 74766Q101		SCHEDULE 13G	Page [	5	of		11	
1	NAMES OF REPORTING F	ERSONS	5						
2	CHECK THE APPROPRIATE  (a) □  (b) □	re box i	IF A MEMBER OF A GROUP						
3	SEC USE ONLY								
4	CITIZENSHIP OR PLACE OF ORGANIZATION  United States								
	NAMES OF	5	SOLE VOTING POWER -0-						
	NUMBER OF SHARES BENEFICIALLY OWNED BY EACH REPORTING PERSON WITH	6	SHARED VOTING POWER 2,209,566						
		7	SOLE DISPOSITIVE POWER -0-						
		8	SHARED DISPOSITIVE POWER 2,209,566						

AGGREGATE AMOUNT BENEFICIALLY OWNED BY EACH REPORTING PERSON

PERCENT OF CLASS REPRESENTED BY AMOUNT IN ROW (9)

CHECK BOX IF THE AGGREGATE AMOUNT IN ROW (9) EXCLUDES CERTAIN SHARES

2,209,566

TYPE OF REPORTING PERSON

10

11

12

CUSIP 1	No.	74766Q101 SCHEDULE 13G Page	
Item 1.			
	(a	n) Name of Issuer:	
		Quanterix Corporation	
	(t	Address of Issuer's Principal Executive Offices:	
		900 Middlesex Turnpike Billerica, Massachusetts 01821	
Item 2.	(a (b (c	Address of Principal Business Office:	
		Integrated Core Strategies (US) LLC c/o Millennium Management LLC 399 Park Avenue New York, New York 10022 Citizenship: Delaware	
		Millennium Management LLC 399 Park Avenue New York, New York 10022 Citizenship: Delaware	
		Millennium Group Management LLC 399 Park Avenue New York, New York 10022 Citizenship: Delaware	
		Israel A. Englander c/o Millennium Management LLC 399 Park Avenue New York, New York 10022 Citizenship: United States	
	(0	d) <u>Title of Class of Securities</u> :	
		common stock, par value \$0.001 per share ("Common Stock")	
	(6	e) <u>CUSIP Number:</u>	
		74766Q101	
Item 3.	If this sta	atement is filed pursuant to Rule 13d-1(b), or 13d-2(b), check whether the person filing is a:	
(a	a) 🗆	Broker or dealer registered under section 15 of the Act (15 U.S.C. 780);	
(t	o) 🗆	Bank as defined in section 3(a)(6) of the Act (15 U.S.C. 78c);	
(0	e) 🗆	Insurance company as defined in section 3(a)(19) of the Act (15 U.S.C. 78c);	
(0	d) 🗆	Investment company registered under section 8 of the Investment Company Act of 1940 (15 U.S.C. 80a-8);	

An investment adviser in accordance with  $\S 240.13d-1(b)(1)(ii)(E);$ 

An employee benefit plan or endowment fund in accordance with §240.13d-1(b)(1)(ii)(F);

(e)

(f)

CUSIP No.		74766Q101	SCHEDULE 13G	Page 7 of 11				
(g)		A parent holding comp	any or control person in accordance with §240.13d-1(b)(1)(ii)(G);					
(h)		A savings association a	s defined in Section 3(b) of the Federal Deposit Insurance Act (12 U	J.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)		Group, in accordance v	vith §240.13d-1(b)(1)(ii)(J).					
Item 4. Own	<u>ership</u>	<u>)</u>						
Provide th	e follo	owing information regard	ling the aggregate number and percentage of the class of securities o	of the issuer identified in Item 1.				
(a) Amount	Renef	icially Owned:						

# (<u>a</u>) <u>.</u>

See response to Item 9 on each cover page.

# (b) Percent of Class:

See response to Item 11 on each cover page.

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

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

See response to Item 5 on each cover page.

CUSIP No.	74766Q101	SCHEDULE 13G	Page	8	of	11
-----------	-----------	--------------	------	---	----	----

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

See response to Item 6 on each cover page.

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

See response to Item 7 on each cover page.

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

See response to Item 8 on each cover page.

The securities disclosed herein as potentially beneficially owned by Millennium Management LLC, Millennium Group Management LLC and Mr. Englander are held by entities subject to voting control and investment discretion by Millennium Management LLC and/or other investment managers that may be controlled by Millennium Group Management LLC (the managing member of Millennium Management LLC) and Mr. Englander (the sole voting trustee of the managing member of Millennium Group Management LLC). The foregoing should not be construed in and of itself as an admission by Millennium Management LLC, Millennium Group Management LLC or Mr. Englander as to beneficial ownership of the securities held by such entities.

#### 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  $\Box$ .

## 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.

Not applicable.

#### Item 8. Identification and Classification of Members of the Group

See Exhibit I.

#### Item 9. Notice of Dissolution of Group

Not applicable.

## Item 10. Certification

By signing below each of the undersigned certifies that, to the best of its 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.

CUSIP No. 74766Q101 SCHEDULE 13G Page 9

Exhibits:

Exhibit I: Joint Filing Agreement, dated as of January 26, 2024, by and among Integrated Core Strategies (US) LLC, Millennium Management LLC, Millennium Group Management LLC and Israel A. Englander.

CUSIP No. 74766Q101 SCHEDULE 13G Page 10 of 11

#### **SIGNATURE**

After reasonable inquiry and to the best of its knowledge and belief, each of the undersigned certifies that the information with respect to it set forth in this statement is true, complete, and correct.

Dated: January 26, 2024

INTEGRATED CORE STRATEGIES (US) LLC

By: Integrated Holding Group LP, its Managing Member

By: Millennium Management LLC, its General Partner

By: /s/Gil Raviv

Title: Global General Counsel

MILLENNIUM MANAGEMENT LLC

By: /s/Gil Raviv

Name: Gil Raviv

Title: Global General Counsel

MILLENNIUM GROUP MANAGEMENT LLC

By: /s/Gil Raviv

Name: Gil Raviv

Title: Global General Counsel

/s/ Israel A. Englander

Israel A. Englander

CUSIP No.	74766Q101	SCHEDULE 13G	Page	11	of	11

### **EXHIBIT I**

## JOINT FILING AGREEMENT

This will confirm the agreement by and among the undersigned that the Schedule 13G filed with the Securities and Exchange Commission on or about the date hereof with respect to the beneficial ownership by the undersigned of the Common Stock, par value \$0.001 per share, of Quanterix Corporation will be filed on behalf of each of the persons and entities named below in accordance with Rule 13d-1(k) under the Securities Exchange Act of 1934, as amended. This Agreement may be executed in two or more counterparts, each of which shall be deemed an original, but all of which together shall constitute one and the same instrument.

Dated: January 26, 2024

INTEGRATED CORE STRATEGIES (US) LLC

By: Integrated Holding Group LP, its Managing Member

By: Millennium Management LLC, its General Partner

By: /s/ Gil Raviv

Name: Gil Raviv

Title: Global General Counsel

MILLENNIUM MANAGEMENT LLC

By: /s/Gil Raviv

Name: Gil Raviv

Title: Global General Counsel

MILLENNIUM GROUP MANAGEMENT LLC

By: /s/Gil Raviv

Name: Gil Raviv

Title: Global General Counsel

/s/ Israel A. Englander

Israel A. Englander