FORM 4

Check this box if no longer subject to Section 16. Form 4 or Form 5 obligations may continue. See Instruction 1(b).

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington,	DC	20549
vvasi ii iytori,	D.C.	20349

STATEMENT	OF CHANGES	IN BENEFICIAL	OWNERSHIP

OMB APPROVAL									
OMB Number:	3235-0287								
Estimated average burden									
hours per response:	0.5								

Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934 or Section 30(h) of the Investment Company Act of 1940

Name and Address of Reporting Person* <u>Sriram Vandana</u>						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]									all applic Directo	or		10% Owner	
(Last)	,	irst)	(Middle)			3. Date of Earliest Transaction (Month/Day/Year) 08/21/2023								X	below)	(give title Other (s below) nief Financial Officer		респу	
900 MIDDLESEX TURNPIKE				4. 1	4. If Amendment, Date of Original Filed (Month/Day/Year)								Individual or Joint/Group Filing (Check Applicable Line)						
(Street)	ICA M	ΙA	01821											X		led by More	•	orting Persor n One Repor	
(City)	(S	tate)	(Zip)		R	Rule 10b5-1(c) Transaction Indication													
			Check this box to indicate that a transaction was made pursuant to a contract, instruction or written plan that is intended to satisfy the affirmative defense conditions of Rule 10b5-1(c). See Instruction 10.																
		Tab	le I - No	n-Deriv	vativ	e Se	curities	s Ac	quired,	Dis	posed o	f, or Be	neficia	lly C	Owned				
1. Title of Security (Instr. 3) 2. Transc Date (Month/D				Execution Date,		Transaction Disposed Code (Instr. 5)		ties Acquired (A) or d Of (D) (Instr. 3, 4 an		d	5. Amour Securitie Beneficia Owned F Reported	es For ally (D) Following (I) (I		n: Direct r Indirect Instr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)				
								Code	v	Amount	Amount (A) or (D)			Transact (Instr. 3 a	ion(s)			msu. 4)	
Common Stock			08/2	1/2023				A	A 12,62		4 ⁽¹⁾ A \$		00 12,624(2)		524 ⁽²⁾		D		
Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities)																			
Derivative Conversion Date Execution Date, T Security or Exercise (Month/Day/Year) if any			ransaction of ode (Instr. Derivative		ive ies ed ed nstr.	6. Date Exercisable and Expiration Date (Month/Day/Year)			7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		De Se	Price of crivative curity str. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s (Instr. 4)	ly	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)			
					Code	v	(A)	(D)	Date Exercisal:		Expiration Date	Title	Amount or Number of Shares						
Stock Option (right to buy)	\$24.32	08/21/2023			A		29,455		(3)	C	08/21/2033	Common Stock	29,45	5	\$0.00	29,455	5	D	

Explanation of Responses:

- 1. Represents shares of stock issuable on vesting of restricted stock units. The restricted stock units vest as to one-third of shares on each of August 21, 2024, August 21, 2025 and August 21, 2026.
- 2. Includes 12,624 restricted stock units.
- 3. The option vests as to one-third of the shares on August 21, 2024, and the remaining two-thirds of the shares vest monthly thereafter over two years in 24 successive equal installments.

Remarks:

/s/ Bonnie McManus, as Attorney-in-Fact

08/22/2023

** Signature of Reporting Person

Date

Reminder: Report on a separate line for each class of securities beneficially owned directly or indirectly.

- * If the form is filed by more than one reporting person, see Instruction 4 (b)(v).
- ** Intentional misstatements or omissions of facts constitute Federal Criminal Violations See 18 U.S.C. 1001 and 15 U.S.C. 78ff(a).

Note: File three copies of this Form, one of which must be manually signed. If space is insufficient, see Instruction 6 for procedure.

Persons who respond to the collection of information contained in this form are not required to respond unless the form displays a currently valid OMB Number.

POWER OF ATTORNEY

Know all by these presents, that the undersigned hereby constitutes and appoints Laurie Churchill, Brian Keane and Bonnie McManus, each of Quanterix Corporation, and Ian McDonald of Covington & Burling LLP signing singly, with full power of substitution, the undersigned's true and lawful attorney-in-fact to:

- (1) execute for and on behalf of the undersigned, forms and authentication documents for EDGAR Filing Access;
- (2) do and perform any and all acts for and on behalf of the undersigned which may be necessary or desirable to complete and execute any such forms and authentication documents;
- (3) execute for and on behalf of the undersigned, in the undersigned's capacity as an officer, director and/or 10% shareholder of the Company, Forms 3, 4 and 5 in accordance with Section 16(a) of the Securities Exchange Act of 1934 and the rules thereunder;
- (4) do and perform any and all acts for and on behalf of the undersigned which may be necessary or desirable to complete and execute any such Form 3, 4 or 5 and timely file such form with the United States Securities and Exchange Commission and any stock exchange or similar authority; and
- (5) take any other action of any type whatsoever in connection with the foregoing which, in the opinion of such attorney-in-fact, may be of benefit to, in the best interest of, or legally required by, the undersigned, it being understood that the documents executed by such attorney-in-fact, on behalf of the undersigned pursuant to this Power of Attorney, shall be in such form and shall contain such terms and conditions as such attorney-in-fact may approve in such attorney-in-fact's discretion.

The undersigned hereby grants to each such attorney-in-fact full power and authority to do and perform any and every act and thing whatsoever requisite, necessary, or proper to be done in the exercise of any of the rights and powers herein granted, as fully to all intents and purposes as the undersigned might or could do if personally present, with full power of substitution or revocation, hereby ratifying and confirming all that such attorney-in-fact, or such attorney-in-fact's substitute or substitutes, shall lawfully do or cause to be done by virtue of this power of attorney and the rights and powers herein granted. The undersigned acknowledges that the foregoing attorneys-in-fact, in serving in such capacity at the request of the undersigned, is not assuming, nor is the Company assuming, any of the undersigned's responsibilities to comply with Section 16 of the Securities Exchange Act of 1934.

This Power of Attorney shall remain in full force and effect until the undersigned is no longer required to file Forms 3, 4 and 5 with respect to the undersigned's holdings of and transactions in securities issued by the Company, unless earlier revoked by the undersigned in a signed writing delivered to the foregoing attorneysin-fact.

IN WITNESS WHEREOF, the undersigned has caused this Power of Attorney to be executed this 21st day of August 2023.

/s/ Vandana Sriram

Vandana Sriram