FORM 4

## **UNITED STATES SECURITIES AND EXCHANGE COMMISSION**

Washington, I	D.C. 20549
---------------	------------

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

Check this box to indicate that a transaction was made pursuant to a contract, instruction or written plan for the purchase or sale of equity securities of the issuer that is intended to satisfy the affirmative

## STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

	OMB APPROVAL								
	OMB Number: 3235-0287								
l	Estimated average burden								
ı	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

	e conditions of ee Instruction 1																		
Name and Address of Reporting Person* <u>Sriram Vandana</u>				2. Issuer Name and Ticker or Trading Symbol     Quanterix Corp [ QTRX ]  3. Date of Earliest Transaction (Month/Day/Year)     09/15/2024									5. Relationship of Reporting Person(s) to Issuer (Check all applicable)  Director 10% Owner					vner	
(Last) (First) (Middle) C/O QUANTERIX CORPORATION 900 MIDDLESEX TURNPIKE													Officer (give title Other (specify below)  Chief Financial Officer						
(Street) BILLERICA MA 01821  (City) (State) (Zip)						Amend	ment, Da	ate o	f Origina	al File	d (Month/Da	y/Year)		6. Indivi ∟ine) ☑	Form	filed by On	e Rep	g (Check A orting Perso n One Repo	on
		Table	I - No	n-Deriva	tive S	Secui	rities /	Acq	uired,	Dis	posed of	, or B	enefic	cially	Own	ed			
1. Title of Security (Instr. 3)  2. Transac Date (Month/Da				Execution Date,		3. Transaction Code (Instr. 8)  4. Securitie Disposed C 5)							ies cially Following	Form (D) or	: Direct r Indirect str. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)			
									Code	v	Amount	(A) c (D)	Pric	, l	Transa	ction(s) 3 and 4)			(111501. 4)
Common Stock 09/15/				09/15/2	/2024		F		78(1)	D	\$12	2.82	.82 36,294(2)			D			
		Tal	ole II -							•	osed of, convertib			-	wne	d			
Derivative   Conversion   Date   Execution Date,   Security   or Exercise   (Month/Day/Year)   if any		4. Transa Code ( 8)			6. Date Exercisable and Expiration Date (Month/Day/Year)			7. Title and Amount of Securities Underlying Derivative Security (Inst 3 and 4)				9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s) (Instr. 4)		10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownershij (Instr. 4)				

Date Exercisable

Expiration Date

## **Explanation of Responses:**

- 1. Shares withheld by Quanterix Corporation solely to cover tax obligations upon the vesting of 263 RSUs.
- 2. Includes 33,880 restricted stock units.

## Remarks:

/s/ Brian Keane, as Attorneyin-Fact

or Shares

Title

09/17/2024

\*\* 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).

Code

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.

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.  $\slash$  Vandana Sriram

Vandana Sriram