FORM 4

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

Washington, D.C. 20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

vvaoriingtori,	D.O. 200 10	

OMB APPROVAL

OMB Number: 3235-0287

Estimated average burden

0.5

hours per response:

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

for the purchase or sale of equity securities of the issuer that is

Instruction 1(b).

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

Check this box to indicate that a transaction was made pursuant to a contract, instruction or written plan

	isfy the affirmative ions of Rule 10b5-uction 10.										
1. Name and Address of Reporting Person*  Sriram Vandana  (Last) (First) (Middle)  C/O QUANTERIX CORPORATION  900 MIDDLESEX TURNPIKE				suer Name <b>and</b> Tick anterix Corp [ substitute of Earliest Transists 5/2024	QTRX ]			5. Relationship of Reporting Person(s) to Issuer (Check all applicable)  Director 10% Owner  Officer (give title Other (specify below)  Chief Financial Officer			
(Street) BILLERICA (City)	MA (State)	01821 (Zip)	4. If <i>i</i>	Amendment, Date o	f Original Fil	ed (Month/Day.	6. Indi Line)	_			
		Table I - No	n-Derivative \$	Securities Acq	uired, Di	sposed of,	or Bei	neficially	/ Owned		
Date			2. Transaction Date (Month/Day/Year)	Execution Date, Transaction Disposed Of (D) (Instr. 3					5. Amount of Securities Beneficially Owned Following Reported Transaction(s)	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)

							,	,			
Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned											
(e.g., puts, calls, warrants, options, convertible securities)											

1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	3A. Deemed Execution Date, if any (Month/Day/Year)	4. Transaction Code (Instr. 8)		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		6. Date Exercisable and Expiration Date (Month/Day/Year)		7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		8. Price of Derivative Security (Instr. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s) (Instr. 4)	Ownership Form: Direct (D)	11. Nature of Indirect Beneficial Ownership (Instr. 4)
				Code	v	(A)	(D)	Date Exercisable	Expiration Date	Title	Amount or Number of Shares				

## Explanation of Responses:

Common Stock

- 1. Shares withheld by Quanterix Corporation solely to cover tax obligations upon the vesting of 263 RSUs.
- 2. Includes 33,617 restricted stock units.

## Remarks:

/s/ Brian Keane, as Attorneyin-Fact

10/16/2024

(Instr. 3 and 4)

\*\* Signature of Reporting Person Da

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.

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