FORM 4

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

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

Washington, D.C. 20549

OMB APPROVAL									
OMB Number: 3235-028									
Estimated average burden									
hours per response:	0.5								

Check this box if no longer subject to Section 16. Form 4 or Form 5 obligations may continue. See 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

1. Name and Address of Reporting Person*  DONNELLY WILLIAM P					2. Issuer Name <b>and</b> Ticker or Trading Symbol  Quanterix Corp [ QTRX ]										all app	licable) tor	ng Per	rson(s) to Is	vner	
(Last)	(Fir	st) (N	Middle)		3. Date of Earliest Transaction (Month/Day/Year) 07/01/2024											Officer (give title below)		Other (s below)	specify	
C/O QUANTERIX CORPORATION					4. If Amendment, Date of Original Filed (Month/Day/Year)							6. Individual or Joint/Group Filing (Check Applicable								
900 MIDDLESEX TURNPIKE													Line)  Form filed by One Reporting Person							
(Street)	*												Form filed by More than One Reporting Person							
,———	1071 1017		1021		Rule 10b5-1(c) Transaction Indication															
(City)	(Sta	ate) (Z	Zip)		Check this box to indicate that a transaction was made satisfy the affirmative defense conditions of Rule 10b5-									nt to a contract, instruction or written plan that is intended to e Instruction 10.						
		Table	I - No	on-Deriva	tive S	Secui	rities	Acc	quired	l, Dis	posed of	, or B	enefici	ially	Own	ed				
1. Title of Security (Instr. 3)  2. Transaction Date (Month/Day/					Year) Executi		eemed ition Date, h/Day/Year)		3. Transaction Code (Instr. 8)  4. Securities Ad Disposed Of (D					and 5) Secu Bene Own		cially Following	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)		7. Nature of Indirect Beneficial Ownership	
									Code	v	Amount	(A) or (D)	Price		Reported Transaction(s) (Instr. 3 and 4)				(Instr. 4)	
Common Stock 07/01/20					024						985(1)	A	\$12.6	68 <sup>(2)</sup> 33		33,583(3)		D		
	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)	rative Conversion Date Execution Date, or Exercise (Month/Day/Year) if any			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)		10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)		
					Code	v	(A)	(D)	Date Exerci	isable	Expiration Date	Title	Amount or Number of Shares							

## **Explanation of Responses:**

- 1. Consists of stock granted in lieu of cash fees for service on the Company's Board of Directors and committees thereof for the second quarter of 2024.
- 2. Closing price of the Company's common stock on the Nasdaq Global Market on July 1, 2024.
- 3. Includes 11,185 restricted stock units.

## Remarks:

/s/ Brian Keane, as Attorneyin-Fact

07/02/2024

\*\* 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).

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 each of Laurie Churchill, Brian Keane, Ian Macdonald and Bonnie McManus as the undersigned's true and lawful attorney-in-fact to:

- (1) execute for and on behalf of the undersigned, in the undersigned's capacity as a beneficial owner of Quanterix Corporation (the "Company"), Forms 3, 4, and 5 in accordance with Section 16(a) of the Securities Exchange Act of 1934 and the rules thereunder and Schedules 13D and 13G in accordance with Section 13 of the Securities Exchange Act of 1934 and the rules thereunder;
- (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 Form 3, 4, or 5 or Schedules 13D or 13G, complete and execute any amendment or amendments thereto, and timely file such form with the U.S. Securities and Exchange Commission (the "SEC") and any securities exchange or similar authority; and
- (3) 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 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 also ratifies hereby any action previously taken by each attorney-in-fact that would have been authorized by this power of attorney if it has been in effect at the time such action was taken. The undersigned acknowledges that each foregoing attorney-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 Sections 13 and 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 and Schedules 13D and 13G 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 each foregoing attorney-in-fact.

IN WITNESS WHEREOF, the undersigned has caused this Power of Attorney to be executed as of this  $20 \, \text{th}$  day of August, 2023.

/s/ William P. Donnelly

William P. Donnelly