FORM 4

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

Washington D.C. 20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

Washington, B.S. 200 10			

OMB APPROVAL								
OMB Number: 3235-02								
Estimated average	burden							
hours per response	e: 0.5							

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 defense conditions of Rule 10b5-1(c). See Instruction 10.

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*     DONNELLY WILLIAM P						2. Issuer Name <b>and</b> Ticker or Trading Symbol Quanterix Corp [ QTRX ]							5. Relationship of Reporting Person(s) to Issuer (Check all applicable)						
DONNELLY WILLIAM P						Zamilian Folk ( cont.)						1	Director			10% Ov			
(Last)	(F	irst) (f	t) (Middle) 3. Date of Earliest Transaction (Month/Day/Year)								]	belov	er (give title v)		Other (s below)	вресну -			
C/O QUANTERIX CORPORATION					10/0	1/202	.4												
900 MIDDLESEX TURNPIKE					4. If Amendment, Date of Original Filed (Month/Day/Year)						6. Individual or Joint/Group Filing (Check Applicable								
(Ott)											Line)	Line)							
(Street) BILLER	ICA N	IA 0	1821											1	_	filed by On-		•	
— DIEEER	1071 10		71021												Perso		re man	опе керс	orung
(City)	(8	tate) (2	Zip)																
Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																			
1. Title of Security (Instr. 3)  2. Transact Date (Month/Day					Execution Date,		3. Transaction Code (Instr. 8) 4. Securities Acquired Disposed Of (D) (Instr. 5)		ıired (A nstr. 3,	) or 4 and	Securi Benefi Owned	Securities For Example 1 For Example 2 For E		Direct Indirect str. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)				
								Code	v	Amount (A) or (D)		rice	Reported Transaction(s) (Instr. 3 and 4)				(instr. 4)		
Common Stock 10/01/20					2024				A		1,000(1)	A	\$	12.5(2)	34	,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)	ive Conversion Date Execution Date, or Exercise (Month/Day/Year) if any			Fransaction of Code (Instr. Derivative		6. Date Exercisable and Expiration Date (Month/Day/Year)		7. Title and Amount of Securities Underlying Derivative Security (Inst 3 and 4)		De Se (Ir	Price of erivative ecurity estr. 5)	9. Number derivative Securities Beneficially Owned Following Reported Transactio (Instr. 4)	y O D O (I)	0. Iwnership orm: irect (D) r Indirect ) (Instr. 4)	Beneficial Ownership (Instr. 4)				
					Code	v	(A)	(D)	Date Exercis	sable	Expiration Date	Title	Amor or Numi of Share	ber					

## **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 third quarter of 2024.
- 2. Closing price of the Company's common stock on the Nasdaq Global Market on October 1, 2024.
- 3. Includes 11,185 restricted stock units.

## Remarks:

/s/ Brian Keane, as Attorneyin-Fact

10/02/2024

\*\*

\*\* Signature of Reporting Person Date

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

- $^{\star}$  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