FORM 4

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

Washington, D.C. 20549

Washington, D.C. 20010	

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

## 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 OLSON LAURIE J						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [ QTRX ]									onship onship on all application of application on all applications on all applicat	,		son(s) to Issuer  10% Owner	
(Last)	(Fi	irst)	(Middle)		3. Date of Earliest Transaction (Month/Day/Year) 04/01/2024									Officer (give title below)			Other (s below)	pecify	
C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE				4. I	4. If Amendment, Date of Original Filed (Month/Day/Year)									Individual or Joint/Group Filing (Check Applicable Line)     X Form filed by One Reporting Person					
(Street) BILLER	ICA M	A	01821												Form f Persor		e than	One Repo	ting
(City)	(S	tate)	(Zip)	Ru	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 - Nor	n-Deriv	ative	e Se	curiti	ies Ac	quired	, Dis	posed o	of, or Be	eneficia	illy O	wnec	ı			
			2. Transaction Date (Month/Day/Year)			2A. Deemed Execution Date, if any (Month/Day/Year)		Code (Instr.		4. Securities Acquired (A) Disposed Of (D) (Instr. 3, 5)			4 and Securiti Benefic		es ally Following	Form (D) o	r Indirect Instr. 4)	7. Nature of Indirect Beneficial Ownership	
									Code	v	Amount	(A) o	Price	1	ransac Instr. 3	tion(s)			Instr. 4)
Common Stock			04/0	/01/2024				М		1,000	) A	\$12	.86	16,238			D		
Common Stock 04			04/0	1/2024				S <sup>(1)</sup>		1,000 D		\$23	.5	5 15,238(2)			D		
		7	able II -						,		osed of converti	,		,	/ned				
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	3A. Deeme Execution if any (Month/Da	Date,	4. Transaction Code (Instr 8)				6. Date Exercisable and Expiration Date (Month/Day/Year)			7. Title at Amount of Securitie Underlyin Derivativ (Instr. 3 a	Deri Sec (Inst	rice of vative urity tr. 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)	Beneficial Ownership (Instr. 4)	
					Code	v	(A)	(D)	Date Exercisa		Expiration Date	Title	Amount or Number of Shares	1					
Stock Option (Right to	\$12.86	04/01/2024			M			1,000	12/31/20	23	01/03/2033	Common Stock	1,000	\$(	0.00	10,124	!	D	

## Explanation of Responses:

- 1. The sale of shares reported on this Form 4 was made pursuant to a Rule 10b5-1 trading plan adopted on August 14, 2023.
- 2. Includes 4,461 restricted stock units.

## Remarks:

/s/ Brian Keane, as Attorney-

04/01/2024

in-Fact

\*\* 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 John Fry and Brian Keane, each of Quanterix Corporation, and Megan Gates, John Condon, Keunjung Cho, Amanda Mei, Anne Leland and Brenda Meyette, each of Mintz, Levin, Cohn, Ferris, Glovsky and Popeo, P.C., 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, 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; (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 and timely file such form with the United States Securities and Exchange Commission and any stock 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 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 28th day of April 2021. /s/ Laurie J. Olson
Laurie J. Olson