SEC For		4			тс						VСЦ А				SCION				
	FORM	UNITED STATES SECURITIES AND EXCHANGE COMMIS Washington, D.C. 20549														ОМВ	APPRO	VAL	
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										HIP	Estim	OMB Number: 3235-0 Estimated average burden hours per response:		3235-0287 en 0.5
1. Name and Address of Reporting Person* <u>OLSON LAURIE J</u>						2. Issuer Name and Ticker or Trading Symbol <u>Quanterix Corp</u> [QTRX]									5. Relationship of Reporting Person(s) to Issuer (Check all applicable) X Director 10% Owner Officer (give title Other (specify below) below)				
(Last) (First) (Middle) C/O QUANTERIX CORPORATION,						3. Date of Earliest Transaction (Month/Day/Year) 01/03/2022													specify
900 MIDDLESEX TURNPIKE						4. If Amendment, Date of Original Filed (Month/Day/Year)									6. Individual or Joint/Group Filing (Check Applicable Line)				
(Street) BILLERICA MA 01821															X Form filed by One Reporting Person Form filed by More than One Reporting Person				
(City) (State) (Zip)																			
		Tal	ole I - Nor	n-Deriv	/ati	ve Se	ecuritie	es Acq	uired,	Dis	posed o	of, o	or Bene	eficiall	y Owned			,	
Date				te onth/Day/Year) i		2A. Deemed Execution Date, if any (Month/Day/Year)		3. Transaction Code (Instr. 8)		4. Securities Acquired (A) Disposed Of (D) (Instr. 3, 4 5)			5. Amou Securitie Benefici Owned F	es ally 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 01/03						2022		Α		2,668	(1)	¹⁾ A \$) 7,2	7,208 ⁽²⁾		D		
			Table II -								osed of, onvertil				Owned				
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	Execution Date, Day/Year) if any		4. Transaction Code (Instr. 8)		Derivative E		6. Date Exercisable and Expiration Date (Month/Day/Year)			An Se Un De	Title and nount of ccurities nderlying erivative s istr. 3 and	Security	8. Price of Derivative Security (Instr. 5)	9. Number derivative Securitie Beneficia Owned Following Reported Transacti (Instr. 4)	e s ally g	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4	Beneficial Ownership t (Instr. 4)

/s/ Brian Keane, as Attorney-in-01/05/2022

\$<mark>0.00</mark>

4,002

D

Fact

Expiration Date

01/03/2032

Title

Common

Stock

Date Exercisable

12/31/2022

(D)

1. Represents restricted stock units that vest as to 100% of the shares on December 31, 2022 granted pursuant to the Quanterix Corporation 2018 Non-Employee Director Compensation Policy. Each restricted stock unit represents the right to receive one share of the Issuer's common stock.

v

Code

A

(A)

4,002⁽³⁾

3. Represents non-qualified stock option to purchase 4,002 shares of common stock granted pursuant to the Quanterix Corporation 2018 Non-Employee Director Compensation Policy.

** Signature of Reporting Person Date

Amount or Number of Shares

4,002

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).

01/03/2022

Stock Option (Right to

Buy)

Remarks:

\$42.41

2. Includes 7,208 restricted stock units.

Explanation of Responses:

** 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-infact'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