SEC For	rm 4																		
FORM 4 UNITED STA				TES SECURITIES AND EXCHANGE COMMISSION Washington, D.C. 20549												OMB APPROVAL			
Section 16. Form 4 or Form 5 obligations may continue. See					d purs	IT OF CHANGES IN BENEFICIAL OWNE pursuant to Section 16(a) of the Securities Exchange Act of 1934 or Section 30(h) of the Investment Company Act of 1940									SHIP	Estim	Estimated average burden		3235-0287 en 0.5
1. Name and Address of Reporting Person [*] Roskey Mark T.						2. Issuer Name and Ticker or Trading Symbol <u>Quanterix Corp</u> [QTRX]									eck all appli Directo	cable) or	, 10% Ow		
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						3. Date of Earliest Transaction (Month/Day/Year) 02/08/2021									X Officer (give title Other (specify below) below) SVP, Commercial & Accelerator				
(Street) BILLERICA MA 01821					, 4. l	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 Form filed by More than One Reporting Person 			on .	
(City)	(City) (State) (Zip)																		
		Tab	le I - Nor	n-Deriv	ative	e Se	curitie	s Ac	cquired,	Dis	sposed o	of, o	or Ber	eficial	ly Owned	d			
1. Title of Security (Instr. 3) 2. Transa Date (Month/E					ar)	2A. Deemed Execution Date, if any (Month/Day/Year		Code	Transaction Code (Instr.					Securitie Benefici Owned I	5. Amount of Securities Beneficially Owned Following		vnership 1: Direct r Indirect 1str. 4)	7. Nature of Indirect Beneficial Ownership	
									Code	v	Amount	unt (A) or (D) Pi		Price	Reported Transaction(s) (Instr. 3 and 4)				(Instr. 4)
Common Stock 02/08/					/202	2021			Α		3,373	,373 ⁽¹⁾ A S		\$ <mark>0.0</mark>	25,105 ⁽²⁾			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)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year	3A. Deeme Execution if any (Month/Da	Date,	4. Transacti Code (Ins 8)			tive ties red sed 3, 4	6. Date E: Expiration (Month/D:	n Date	e	Amount of Securities Underlying Derivative Sec (Instr. 3 and 4)		14)	8. Price of Derivative Security (Instr. 5)	9. Numbe derivative Securities Beneficia Owned Following Reported Transacti (Instr. 4)	s Ily I	10. Ownershij Form: Direct (D) or Indirect (I) (Instr. 4	Beneficial Ownership (Instr. 4)
					Code	v	(A)	(D)	Date Exercisat		Expiration Date	Title		Amount or Number of Shares					

buy) Explanation of Responses:

\$77.79

Stock Option (right to

1. Represents shares of stock issuable on vesting of restricted stock units. The restricted stock units vest as to 25% of the shares on February 8, 2022, and the remaining 75% vest in 36 equal monthly installments at the end of each month beginning on February 28, 2022.

7,870

(3)

02/08/2021

2. Includes 17,742 restricted stock units.

3. The option vests as to 25% of the shares on February 8, 2022, and the remaining 75% vest in 36 equal monthly installments at the end of each month beginning on February 28, 2022. Remarks:

> /s/ Brian Keane, Attorney-in-02/10/2021 fact

7,870

\$<mark>0.00</mark>

7,870

D

Common Stock

02/08/2031

** 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.

A

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.