FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

.....g.c., 2101 200

STATEMENT	OF CHANGES	IN BENEFICIAL	. OWNERSHIP

OMB Number: 3235-0287
Estimated average burden
hours per response: 0.5

OMB APPROVAL

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

Name and Address of Reporting Person* Roskey Mark T.						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]								(Check	all applica Director	able)	p Person(s) to Issuer 10% Owner Other (spec		/ner
	(First) (Middle) QUANTERIX CORPORATION, IIDDLESEX TURNPIKE					3. Date of Earliest Transaction (Month/Day/Year) 02/18/2020								X	Officer (give title Other (specify below) SVP, Commercial & Accelerator				·
(Street) BILLER	ICA M	IA	01821		4.	4. If Amendment, Date o				of Original Filed (Month/Day/Year)					vidual or Joint/Group Filing (Check Applicable Form filed by One Reporting Person Form filed by More than One Reporting Person				
(City)	(S	itate)	(Zip)																
Table I - Non-Deriv 1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/				action	on 2A. Deemed Execution Date,		3. 4. Securities			f, or Beneficially Acquired (A) or (D) (Instr. 3, 4 and 5)			5. Amount of Securities Beneficially Owned 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 (Ins Transaction(s) (Instr. 3 and 4)				(Instr. 4)
Common	Stock			02/18	/2020				M		5,000	A	\$	2.96	19,1	19,184 ⁽¹⁾ D		D	
Common	Stock			02/18	/2020				S		5,000 ⁽²⁾	D	\$27.	8388 ⁽³⁾) 14,184 ⁽¹⁾ D		D		
Common	Stock			02/18	/2020				A		8,357(4)	A	\$	0.00	22,541 ⁽¹⁾ D		D		
			Table I								posed of, convertib				wned				
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	ise (Month/Day/Year) /e	3A. Deemed Execution Date, if any (Month/Day/Year)		4. Transaction Code (Instr. 8)		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3 4 and 5)		Expiration I (Month/Day			7. Title and Amo of Securities Underlying Derivative Secur (Instr. 3 and 4)		curity (3. Price of Derivative Security (Instr. 5)	9. Number derivative Securities Beneficial Owned Following Reported	e s Illy	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership (Instr. 4)
				Code V (A) (D) Exercis			Expiration Date	Title	or Nu of	mber ares		Transaction(s) (Instr. 4)							
Stock Option (Right to Buy)	\$2.96	02/18/2020			М			5,000		(5)	09/23/2024	Commo Stock		,000,	\$0.00	114,45	5	D	
Stock Option (right to buy)	\$27.76	02/18/2020			A		19,500			(6)	02/18/2030	Commo	on 19	,500	\$0.00	19,500		D	

Explanation of Responses:

- 1. Includes 19.602 restricted stock units.
- 2. The sale of shares reported on this Form 4 was made pursuant to a Rule 10b5-1 trading plan adopted on December 13, 2019.
- 3. This transaction was executed in multiple trades at prices ranging from \$27.4100 to \$28.4099. The price reported reflects the weighted average sale price. The reporting person undertakes to provide upon request to the SEC staff, the issuer or any security holder of the issuer full information regarding the number of shares and prices at which the transaction was effected.
- 4. Represents shares of stock issuable on vesting of restricted stock units. The restricted stock units vest as to 25% of the shares on February 18, 2021 and vest as to an additional 2.083% of the shares per month for 36 months thereafter.
- 5. Option is fully vested.
- 6. The option vests as to 25% of the shares on February 18, 2021 and vests as to an additional 2.083% of the shares per month for 36 months thereafter.

Remarks:

/s/ Brian Keane, Attorney-infact 02/20/2020

** 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.