FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

OMB APPROVAL									
OMB Number:	3235-0287								
Estimated average burd	den								
hours per response:	0.5								

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

1. Name ar		2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]									all appli Directo	licable) tor		Person(s) to Issuer 10% Owner Other (specify					
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						Date of /18/20		est Tra	nsactio	n (Mon	nth/Day/Year)	X Officer (give title Other (specify below) below) SVP, Commercial & Accelerator					·		
(Street) BILLER (City)	ICA M	Α (01821 (Zip)		4. 1	4. If Amendment, Date of Original Filed (Month/Day/Year)									rvidual or Joint/Group Filing (Check Applicable Form filed by One Reporting Person Form filed by More than One Reporting Person				
	Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																		
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Y				ion	2A. De Execu if any (Monti	eemed	i Date,	3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 and			5. Amo Securit Benefic Owned		unt of ies ially Following	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)		7. Nature of Indirect Beneficial Ownership	
										v	Amount	(A) or (D)	Price			orted saction(s) r. 3 and 4)			(Instr. 4)
Common Stock 05/1				05/18/2	020	20			M		5,000	Α	\$2	.96	26,	26,741(1)		D	
Common Stock 05/18/202			020	20			S		4,600(2)	D	\$25.5	5479 ⁽³⁾	22,	22,141 ⁽¹⁾		D			
Common Stock 05/18/202			020)20					400(2)	D	\$26	\$26.23(4)		21,741(1)		D			
		Т	able								sposed of , converti				wned				
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	3A. Deemed Execution Date, if any (Month/Day/Year)		4. Transa Code (8)			6. Date Exerc Expiration D (Month/Day/		ate	7. Title and Amount of Securities Underlying Derivative Secur (Instr. 3 and 4)		rity	Price of rivative curity str. 5)	9. Number derivative Securities Beneficiall Owned Following Reported Transactio (Instr. 4)	y O F D oi (I)	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)	
					Code	v	(A)	(D)	Date Exerc	isable	Expiration Date	Title	Amo or Num of Shar	ber					
Stock Option (Right to	\$2.96	05/18/2020			M			5,000	(!	5)	09/23/2024	Common Stock	5,0	00	\$0.00	99,455	5	D	

Explanation of Responses:

- 1. Includes 17,051 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 \$25.19 USD to \$26.06 USD; the price reported above reflects the weighted average sale price.
- 4. This transaction was executed in multiple trades at prices ranging from \$26.22 USD to \$26.25 USD; the price reported above reflects the weighted average sale price.
- 5. Option is fully vested.

Remarks:

/s/ Brian Keane, Attorney-infact

05/20/2020

** 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.