SEC F	Form 4
-------	--------

П

FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

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

Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934 or Section 30(h) of the Investment Company Act of 1940

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

1. Name and Address of Reporting Person [*] Mattoon Dawn			2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]		ationship of Reporting Pe < all applicable) Director	10% Owner	
(Last) (First) (Middle)		()	3. Date of Earliest Transaction (Month/Day/Year) 02/18/2020	X	Officer (give title below) Sr. VP, Researc	Other (specify below) h Products	
C/O QUANTERIX CORPORATION,							
900 MIDDLESEX TURNPIKE		KE					
(7)			4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Indi Line)	vidual or Joint/Group Filing (Check Applicable		
(Street) BILLERICA	МА	01821		X	Form filed by One Re	porting Person	
	10171	01021			Form filed by More th Person	an One Reporting	
(City)	(State)	(Zip)					

Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned

1. Title of Security (Instr. 3)	2. Transaction Date (Month/Day/Year)	2A. Deemed Execution Date, if any (Month/Day/Year)			4. Securities A Disposed Of (I 5)			5. Amount of Securities Beneficially Owned Following Reported	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership
			Code	v	Amount	(A) or (D)	Price	Transaction(s) (Instr. 3 and 4)		(Instr. 4)
Common Stock	02/18/2020		A		11,143(1)	Α	\$0.00	33,247 ⁽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. Deemed Execution Date, if any (Month/Day/Year)	4. Transa Code (8)		of		6. Date Exercisable and Expiration Date (Month/Day/Year)		7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		8. Price of Derivative Security (Instr. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s) (Instr. 4)	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)
				Code	v	(A)	(D)	Date Exercisable	Expiration Date	Title	Amount or Number of Shares				
Stock Option (right to buy)	\$27.76	02/18/2020		A		26,000		(3)	02/18/2030	Common Stock	26,000	\$0.00	26,000	D	

Explanation of Responses:

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 18, 2021 and vest as to an additional 2.083% of the shares per month for 36 months thereafter.

2. Includes 27,665 restricted stock units.

3. 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, as Attorney-in-02/20/2020

<u>Fact</u>

** 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.