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

# FORM 4

Check this box if no longer subject

UNITED STATES SECURITIES AND EXCHANGE COMMISSION Washington, D.C. 20549

OMB APPROVAL

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

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		Estimated average burden hours per response: 0.5									
	or Section 30(h) of the Investment Company Act of 1940	1										
1. Name and Address of Reporting Person <sup>*</sup> Mattoon Dawn	2. Issuer Name <b>and</b> Ticker or Trading Symbol <u>Quanterix Corp</u> [ QTRX ]	(Check all applicab Director	10% Owner									
(Last) (First) (Middle) C/O QUANTERIX CORPORATION,	3. Date of Earliest Transaction (Month/Day/Year) 03/13/2020	• X Officer (giv below) Sr. VP,	ve title Other (specify below) , Research Products									
900 MIDDLESEX TURNPIKE	4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Individual or Joint/Group Filing (Check Applicable										
(Street) BILLERICA MA 01821		1	by One Reporting Person by More than One Reporting									
(City) (State) (Zip)		Person										
Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned												

1. Title of Security (Instr. 3)	Date (Month/Day/Year)	2A. Deemed Execution Date, if any (Month/Day/Year)	Transaction Code (Instr.		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 and 5)				Form: Direct (D) or Indirect (I)	7. Nature of Indirect Beneficial Ownership
			Code	v	Amount	(A) or (D)	Price	Reported Transaction(s) (Instr. 3 and 4)	(Instr. 4)	(Instr. 4)
Common Stock	03/13/2020		S		566(1)	D	\$16.0434 <sup>(2)</sup>	32,566 <sup>(3)</sup>	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)		5. Number 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				

#### Explanation of Responses:

1. The sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 1,807 RSUs.

2. This transaction was executed in multiple trades at prices ranging from \$15.88 to \$16.25. 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. 3. Includes 25,520 restricted stock units.

### Remarks:

#### /s/ Brian Keane, as Attorney-03/17/2020 in-Fact

\*\* Signature of Reporting Person Date

Reminder: Report on a separate line for each class of securities beneficially owned directly or indirectly.

 $^{\ast}$  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.

## STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP