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 bu	rden									

Estimated average burden	

1. Name and Address of Reporting Person <sup>*</sup> <u>Mattoon Dawn</u> (Last) (First) (Middle) C/O QUANTERIX CORPORATION,			2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [ QTRX ] 3. Date of Earliest Transaction (Month/Day/Year) 06/04/2019	(Check a	5. Relationship of Reporting Person(s) to Issuer (Check all applicable) Director 10% Owner X Officer (give title Other (specify below) below) Sr. VP, Research Products				
900 MIDDLESEX TURNPIKE			4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Individual or Joint/Group Filing (Check Applicable Line)					
Street) BILLERICA MA 01821		01821		X	Form filed by One Reporti Form filed by More than O Person	0			
(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)	Transaction Code (Instr.					5. Amount of Securities Beneficially Owned Following Reported	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
			Code		Amount	(A) or (D)	Price	Transaction(s) (Instr. 3 and 4)		
Common Stock	06/04/2019		S		64(1)	D	\$25.81	23,090	D	
Common Stock	06/04/2019		S		<b>49</b> <sup>(1)</sup>	D	\$25.81	23,041 <sup>(2)</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)		of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		6. Date Exerc Expiration Da (Month/Day/)	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			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 208 RSUs and 160 RSUs, respectively, and were effected pursuant to a Rule 10b5-1 trading plan adopted on December 14, 2018.

2. Includes 19,496 restricted stock units.

**Remarks:** 

/s/ Brian Keane, as Attorney-

06/05/2019

\*\* Signature of Reporting Person

in-Fact

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.

Ī