FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

STATEMENT	OF CHANGES	IN BENEFICIAL	OWNERSHIP
_		_	

OWR APPRO	JVAL			
OMB Number:	3235-0287			
Estimated average burd	en			
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 and Address of Reporting Person* Mattoon Dawn						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]								(Check	all app	olicable)	Person(s) to Issuer 10% Owner Other (specify	
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE					3. Date of Earliest Transaction (Month/Day/Year) 08/01/2019								X	belov	w) ``	below arch Products)``	
(Street) BILLERICA MA 01821 (City) (State) (Zip)				_ 4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)							Individual or Joint/Group Filing (Check Applicable Line) X 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/Ye			.	Execution Date,		3. Transaction Code (Instr. 8) 4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 and				nd 5) Secu Bend Own		ficially d Following	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership				
									Code	v	Amount	(A) or (D)	Price			action(s) 3 and 4)		(Instr. 4)
Common Stock 08/01/201					019	19		S		49(1)	D	\$31.61	.53 ⁽²⁾ 2		22,880	D		
Common Stock 08/01/201					019	19		S		64(1)	D	\$31.61	153 ⁽²⁾	2:	2,816 ⁽³⁾	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	e (Month/Day/Year) if any (Month/	eemed 4. Transa Code h/Day/Year) 8)				Expiration Date (Month/Day/Year)		eate Year)	7. Title and Amount of Securities Underlying Derivative Security (Instr. and 4)		Deriv Secu (Instr	vative urity tr. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s (Instr. 4)	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	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 160 RSUs and 208 RSUs, respectively, and were effected pursuant to a Rule 10b5-1 trading plan adopted on December 14, 2018.
- 2. This transaction was executed in multiple trades at prices ranging from \$31.55 to \$31.795. 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 18,760 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorneyin-Fact

08/02/2019

** 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.