FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

STATEMENT	OF CHANGES	IN BENEFICIAL	OWNERSHIP

l	OMB APPRO	DVAL
	OMB Number:	3235-0287
l	Estimated average burd	en
l	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

Name and Address of Reporting Person* Mattoon Dawn					2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]								Check	all app	olicable)	g Person(s) to Is		
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 113 HARTWELL AVENUE						3. Date of Earliest Transaction (Month/Day/Year) 05/01/2019								X	belov	w) ``	below arch Products)``
(Street) LEXINGTON MA 02421					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				
(City)	(St	ate) (Zip)											Form filed by More than One Reporting Person				orting
		Tabl	e I - N	Non-Deriv	ative	Seci	uritie	s Ac	quire	ed, Di	isposed o	f, or E	Benefici	ally (Owne	ed		
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Ye						Execution Date,		·	3. Transaction Code (Instr. 8) 4. Securities Acquired (A Disposed Of (D) (Instr. 3,				and 5)		ount of ities icially d Following rted	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
									Code	v	Amount	(A) or (D)	Price		Trans	action(s) 3 and 4)		(111511.4)
Common Stock 05/01/201)19	19			S		65(1)	D	\$22.59	25 ⁽²⁾	2	3,806(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	3. Transaction Date (Month/Day/Year)	Execu	eemed Ition Date, h/Day/Year)	4. Transa Code (8)		5. Numof Derive Securion Acquired (A) or Disposof (D) (Instrand 5	ative rities ired sed	Expir (Mon	te Exer ration D th/Day/	Year) Expiration	7. Title Amour Securi Underl Deriva Securi and 4)	nt of ties ying			9. Number of derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)

Explanation of Responses:

- 1. The sale of shares reported on this Form 4 was made solely to cover tax obligations upon the vesting of 208 RSUs and was 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 \$22.7812 to \$22.40. 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 21,784 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorneyin-Fact 05/02/2019

** 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.