FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL										
OMB Number:	3235-0287									
Estimated average t	ourden									

0.5

hours per response:

Check this box if no longer subject to
Section 16. Form 4 or Form 5
obligations may continue. See
Instruction 1(h)

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*					2. Is	2. Issuer Name and Ticker or Trading Symbol									5. Relationship of Reporting Person(s) to Issuer (Check all applicable)				
Mattoon Dawn							Quanterix Corp [QTRX]								CHECK	Direc	ctor		Owner
(I cot)	3 D	Date of Earliest Transaction (Month/Day/Year)								X	Offic belov	er (give title w)	Other below	(specify)					
(Last) (First) (Middle) C/O OUANTERIX CORPORATION,							02/26/2019								Sr VP, Stgc Mrktg & Assay Tech				
113 HARTWELL AVENUE																			
						. 4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Individual or Joint/Group Filing (Check Applicable Line)				
(Street)	TON	N # A		12.421										l,	X	Forn	n filed by One	e Reporting Per	son
LEXING	TON	MA)2421											Form filed by More than One Reporting Person				
(City)	((Stat	e) (a	Zip)											F 613	OH			
			Tabl	e I - N	Non-Deriv	ative	Sec	uritie	s Ac	quire	ed, Di	sposed 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 A Disposed Of (d 5) Sec Ben Owr		ount of ities 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		Reported Transaction(s) (Instr. 3 and 4)		action(s)		(Instr. 4)
Common Stock 02/26/201)19	.9			S		890(1)	D	\$23.72	35 ⁽²⁾	16,790(3)		D		
			Та	ble II	- Derivat (e.g., p	ive S uts, c	ecuri alls,	ties warr	Acqı ants	uired , opti	, Disp ions,	osed of, convertib	or Be	neficial curities)	ly Ov	vned			
1. Title of Derivative Security (Instr. 3)	rivative Conversion Date Execution Date, curity or Exercise (Month/Day/Year) if any				ransaction ode (Instr.		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		6. Date Exercisable and Expiration Date (Month/Day/Year)			and nt of ties ying tive ty (Instr. 3			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)		
					•	Code	v	(A)	(D)	Date Exerc	cisable	Expiration Date	Title	Amount or Number of Shares					

Explanation of Responses:

- 1. The sale of shares reported on this Form 4 was made solely to cover tax obligations upon the vesting of 2,500 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 \$23.56 to \$23.751. 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 15,180 shares of restricted stock units.

Remarks:

/s/ Brian Keane, as Attorneyin-Fact 02/27/2019

** 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.