FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

STATEMENT	OF CHANGES	IN BENEFICIAL	OWNERSHIP

OMB APPROVAL									
OMB Number:	3235-0287								
Estimated average burden									
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	tionship of Reportin all applicable) Director Officer (give title		10% C			
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						3. Date of Earliest Transaction (Month/Day/Year) 04/01/2020									X	below) Sr. VP, Research Products				
(Street) BILLERICA MA 01821 (City) (State) (Zip)					4. If	Amend	ment, [Date (of Ori	iginal F	iled (Month	n/Da	y/Year		i. Indiv. ine) X	Form	Joint/Group filed by One filed by Mor on	e Rep	orting Pers	son
		Table	1 - 1	Non-Deriva	tive	Secui	rities	Acc	quir	ed, D	isposed	of	, or E	Benefic	ially	Own	ed			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Ye			Execution Date,		, t	3. Transaction Code (Instr. 8) 4. Securities A Disposed Of (I			es Ad Of (D	cquired)) (Insti	d (A) or c. 3, 4 and	nd 5) Secui Benet		cially d Following	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)		7. Nature of Indirect Beneficial Ownership (Instr. 4)			
								c	ode	v	Amount	(1	A) or D)	Price		Transa	action(s) 3 and 4)	(11130	+,	(111301. 4)
Common	Stock			04/01/202	0				S		97(1)		D	\$17.44	32,469 ⁽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 (Instr. 3) 3. Transaction Date Execution Date, if any (Month/Day/Year)		Code 8)	Transaction of Code (Instr. Derivative			Expiration Date (Month/Day/Year)			7. Title and Amount of Securities Underlying Derivative Security (Insti 3 and 4) Amoun or Numbe of Title Shares		unt ber		9. Number derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	y	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership (Instr. 4)				

Explanation of Responses:

- 1. The sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 518 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$17.12 to \$17.76. 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 25,002 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorneyin-Fact

04/01/2020

** 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.