FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington D.C. 20549

wasinington,	D.C.	20343	

OMB APPROVAL

OMB Number: 3235-0287

Estimated average burden

0.5

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

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

Name and Address of Reporting Person* Mattoon Dawn					2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]									Check	ionship of Reportin all applicable) Director Officer (give title		ng Person(s) to Is 10% Ov Other (s		wner
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE					3. Date of Earliest Transaction (Month/Day/Year) 10/01/2020									X	below) Sr. VP, Research			below)	Specify
(Street) BILLER (City)	BILLERICA MA 01821				4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Indiv Line) X	ividual or Joint/Group Filing (Check Applicable Form filed by One Reporting Person Form filed by More than One Reporting Person					
		Table	1 - 1	Non-Deriva	tive	Secui	rities	Acc	quire	ed, D	isposed c	of, or E	Benefic	cially	Own	ed			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Yea			ear) i	2A. Deemed Execution Date, if any (Month/Day/Year)		Ti				Acquired (A) or (D) (Instr. 3, 4 ar		nd 5) Secur Benef		rities F ficially (ed Following I		n: Direct or rect (I)	7. Nature of Indirect Beneficial Ownership		
								ode	v	Amount	(A) or (D)	Price		Transa	ansaction(s) estr. 3 and 4)		tr. 4)	(Instr. 4)	
Common Stock 10/01/20			10/01/202	.0			S		159(1)	D	\$34.12	4.1224 ⁽²⁾		31,503(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)	Derivative Conversion Date Execution Date, Gecurity or Exercise (Month/Day/Year) if any				Transaction of Code (Instr. Derivat		tive ties ed sed	Expiration Date (Month/Day/Year)			Amor Secu Unde Deriv Secu	7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		rice of vative urity tr. 5)	9. Number derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	y	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)	
					Code	v	(A)	(D)	Date Exe	e rcisabl	Expiratior e Date	ı 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 518 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$33.881 to \$34.51. 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 21,894 restricted stock units.

Remarks:

<u>/s/ Brian Keane, as Attorney-in-Fact</u>

10/02/2020

** 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.