Instruction 1(b).

FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Check this box if no longer subject	STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP
to Section 16. Form 4 or Form 5	
obligations may continue. See	

OMB APPROVAL 3235-0287 Estimated average burden hours per response: 0.5

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]									ionship of Reportinall applicable) Director Officer (give title		ng Person(s) to Is 10% O Other (wner
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						3. Date of Earliest Transaction (Month/Day/Year) 01/05/2021									below below Sr. VP, Research Products				
(Street) BILLERICA MA 01821 (City) (State) (Zip)				1	4. If Amendment, Date of Original Filed (Month/Day/Year)									i. Indiv ine) X					
		Table	I - I	Non-Deriva	tive	Secui	rities	Acc	quir	ed, Di	isposed o	f, or E	3enefic	ially	Own	ed			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Yea				ear) E	2A. Deem Execution if any (Month/D		ָּרְ זַׁ כ	3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 ar			nd 5) Sec Bei Ow		Amount of curities neficially ned Following ported		wnership n: Direct or rect (I) tr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
								c	Code	v	Amount	(A) or (D)	Price		Transa	nsaction(s) str. 3 and 4)		u. 4)	(11150.4)
Common Stock 01/05/2023					1	1			S		170(1)	D	\$44.13	\$44.1317 ⁽²⁾		31,016 ⁽³⁾		D	
		Tal	ble	II - Derivati (e.g., pu						,	posed of, convertil			•)wne	d			
1. Title of Derivative Security (Instr. 3)	vative Conversion Date Execution Date, Irity or Exercise (Month/Day/Year) if any				action (Instr.				Expiration Date (Month/Day/Year)			le and unt of rities rrlying vative rity (Instr. i 4) Amount or Number of	nt		9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s) (Instr. 4)		10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)	
					Code	l v	(A)	(D)		rcisable		Title							

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 \$43.91 to \$44.62. 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 20,340 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorneyin-Fact

01/06/2021

** 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.