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									
OMB Number:	3235-0287								
Estimated average burden									
hours per response	e: 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 <b>and</b> Ticker or Trading Symbol Quanterix Corp [ QTRX ]										all app Direc		ng Per	rson(s) to Is 10% O Other (	wner	
(Last) (First) (Middle) C/O QUANTERIX CORPORATION,						3. Date of Earliest Transaction (Month/Day/Year) 02/02/2022										below)  Sr. VP, Di		iagno	below)	
900 MIDDLESEX TURNPIKE  (Street)  BILLERICA MA 01821				4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)									6. Individual or Joint/Group Filing (Check Applicable Line)  X Form filed by One Reporting Person Form filed by More than One Reporting Person					on	
(City)	(St		Zip)	lon Deriva	tive Securities Acquired, Disposed of, or Beneficially Owned															
1. Title of Security (Instr. 3)  2. Transaction Date (Month/Day/Yea				ear) i	2A. Deemed Execution Date,		3 T C	3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 an			d (A) or	5. Am Secur Benef Owne		ount of ties cially d Following	Forn (D) o Indir	n: Direct or rect (I)	7. Nature of Indirect Beneficial Ownership	
								c	Code	v	Amount		(A) or (D)	Price		Reported Transaction(s) (Instr. 3 and 4)		(Instr. 4)		(Instr. 4)
Common Stock 02/02/20			02/02/202	2				S		196(1)		D	\$31.20	66(2)	37,793(3)			D		
Common Stock 02/02/2022			2			S		71 <sup>(1)</sup> D \$		\$32.	2.15 3		37,722 <sup>(3)</sup>		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)	Conversion or Exercise Price of Derivative Security  Date (Month/Day/Year)   Execution Date, if any (Month/Day/Year)   Execution Date, if any (Month/Day/Year)   Security   Sec					5. Num of Deriva Securi Acquir (A) or Dispos of (D) (Instr.: and 5)	tive ties ed sed	Exp	ate Ex iration nth/Da	y/Year) Securities Underlying Derivative Security (Inst 3 and 4)		unt of rities rlying rative rity (Instr.	Deri Sec (Ins	rice of ivative urity tr. 5)	ive derivative Securities		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	Expiration Date Ti		Title	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 762 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$30.84 to \$31.83. 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 11,704 restricted stock units.

## Remarks:

/s/ Brian Keane, as Attorney-

in-Fact

02/04/2022

\*\* 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.