FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

wasiiiigton,	D.C. 20549	

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

OMB APPROVAL									
OMB Number: 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]									Check	all app	o of Reportir dicable) tor er (give title	ng Pe	erson(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) 02/09/2022									X	below) Sr. VP, Diagr		iagn	below)	.,,
(Street) BILLERICA MA 01821 (City) (State) (Zip)				4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Indiv Line) X	′						
		Table	1 - 1	Non-Deriva	tive S	Secui	rities	Ac	quir	ed, D	isposed (of, or	Benefic	ially	Own	ed			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Ye.				2A. Deemed Execution Date, if any (Month/Day/Year		,	3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 ar			nd 5) Secu Bene		icially d Following	Fori (D) (Indi	Ownership m: Direct or irect (I) str. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)		
								Code	v	Amount	(A) or (D)	Price		Transaction(s) (Instr. 3 and 4)		(3		(111301.4)	
Common Stock 02/09/2023				.2			s 318 ⁽¹⁾		D	D \$34.9686 ⁽²⁾		37,404 ⁽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	3. Transaction Date (Month/Day/Year)	Execution Date, T			action (Instr.				Expiration Date (Month/Day/Year)			tle and unt of urities erlying vative urity (Instr. d 4) Amount or Number of	nt				10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficia Ownership (Instr. 4)
					Code V (A) (D			(D)		ercisable		Title		1					

Explanation of Responses:

- 1. The sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 919 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$34.9435 to \$34.985. 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 10,785 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorney-** Signature of Reporting Person

in-Fact

02/09/2022

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.