Instruction 1(b).

FORM 4

Check this box if no longer subject

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington,	D.C.	20549
-------------	------	-------

STATEMENT	OF CHAI	NGES IN E	BENEFICIAL

OMB APPROVAL								
OMB Number: 3235-0287								
Estimated average burden								
hours per response: 0.								

to Section 16. Form 4 or Form 5 obligations may continue. See 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	,	ng Pei	10% O	
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						3. Date of Earliest Transaction (Month/Day/Year) 10/01/2021								X		below) Sr. VP, Diagno			,
(Street) BILLER (City)	ICA M.	A 0	1821 Zip)		4. If A	4. If Amendment, Date of Original Filed (Month/Day/Year)								. Indiv ine) X	Form Form	Form filed by One Reporting Person Form filed by More than One Reporting Person			
		Table	I - No	on-Deriva	tive S	Secur	rities	Acc	quired	d, Dis	sposed of	, or E	Benefic	ially	Own	ed			
1. Title of Security (Instr. 3) 2. Transactio Date (Month/Day/N				(Year)	Execution Date,			3. Transaction Code (Instr. 8) 4. Securities Acquire Disposed Of (D) (Inst						5. Amount of Securities Beneficially Owned Following Reported		wnership m: Direct or Indirect nstr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)		
									Code	v	Amount	(A) or (D)	Price		Transa	action(s) 3 and 4)			(IIISU. 4)
Common	Stock			10/01/20)21			S		342(1)	D	\$51.2	1.2006		36,848(2)		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)	vative Conversion Date Execution Date, urity or Exercise (Month/Day/Year) if any			5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		Expiration Date			7. Title and Amount of Securities Underlying Derivative Security (Inst 3 and 4)		Deri Sec	rice of ivative urity tr. 5)	9. Number derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	Ownersh Form: Direct (D) or Indirect (I) (Instr.		Beneficial Ownership t (Instr. 4)			
					Code	v	(A)	(D)	Date Exerc	isable	Expiration Date	Title	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 750 RSUs.
- 2. Includes 14,716 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorneyin-Fact

OWNERSHIP

10/04/2021

** 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.