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 resp	onse: 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

1. Name and Address of Reporting Person* <u>Hlavinka Sarah E.</u>						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]									ationship all app Direc	,	ng Per	son(s) to Is	
(Last)	(Fi	st) (N	Middle)			3. Date of Earliest Transaction (Month/Day/Year) 10/02/2023									Office	er (give title v)		Other (s below)	specify
C/O QUANTERIX CORPORATION 900 MIDDLESEX TURNPIKE			4. If A	4. If Amendment, Date of Original Filed (Month/Day/Year)								Indiv ne)	-/						
(Street) BILLER	ICA M	A 0	01821												Form filed by More than One Reporting Person				
(City)	(St	ate) (2	Zip)		$ _{\Box}$	Rule 10b5-1(c) Transaction Indication Check this box to indicate that a transaction was made pursuant to a contract, instruction or written plan that is intended to satisfy the affirmative defense conditions of Rule 10b5-1(c). See Instruction 10.													
		Table	I - No	on-Deriva	tive S	Secui	rities	Acc	quirec	d, Dis	sposed of	, or E	Benefici	ially	Own	ed			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/				Execu (Year) if any		eemed Ition Date, h/Day/Year)		3. Transaction Code (Instr. 8)			Acquii f (D) (In:	red (A) or str. 3, 4 and	nd 5) Securit Benefic Owned		ities Foicially (Did Following (I)		n: Direct or Indirect ostr. 4)	7. Nature of Indirect Beneficial Ownership	
								Code	v	Amount	(A) oi (D)	Price	Reported Transaction(s) (Instr. 3 and 4)				(Instr. 4)		
Common Stock 10/02/20					023				A		561(1)	A	\$26.7	1 ⁽²⁾	39,490 ⁽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)	rative Conversion Date Execution Date, or Exercise (Month/Day/Year) if any		4. Transaction Code (Instr. 8)		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		6. Date Exercisable and Expiration Date (Month/Day/Year)			7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		8. Price of Derivative Security (Instr. 5)		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	v	(A)	(D)	Date Exerc	isable	Expiration Date	Title	Amount or Number of Shares						

Explanation of Responses:

- 1. Consists of stock granted in lieu of cash fees for service on the Company's Board of Directors and committees thereof for the third quarter of 2023.
- 2. Closing price of the Company's common stock on the Nasdaq Global Market on October 2, 2023.
- 3. Includes 8,083 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorney-10/04/2023

** Signature of Reporting Person

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.