SEC Form 4	
FORM	4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

Check this box if no longer subject to Section 16. Form 4 or Form 5 obligations may continue. See Instruction 1(b)
Instruction 1(b).

Check this box to indicate that a transaction was made pursuant to a contract, instruction or written plan for the purchase or sale of equity securities of the issuer that is intended to satisfy the affirmative defense conditions of Rule 10b5-1(c). See Instruction 10.

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934 or Section 30(h) of the Investment Company Act of 1940

OMB APPROVAL							
OMB Number:	3235-0287						
Estimated average	hurden						

Estimated average burden hours per response: 0.5

		on [*]	2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]		tionship of Reporting Pers all applicable) Director	Person(s) to Issuer 10% Owner Other (specify below)	
(Last) (First) (Middle) C/O QUANTERIX CORPORATION,		,	3. Date of Earliest Transaction (Month/Day/Year) 08/20/2024		Officer (give title below)		
900 MIDDLESEX TURNPIKE (Street) BILLERICA MA 01821		01821	4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Indiv Line)	idual or Joint/Group Filing Form filed by One Repo Form filed by More than Person	orting Person	
(City)	(State)	(Zip)					

Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned

1. Title of Security (Instr. 3)	2. Transaction Date (Month/Day/Year)	2A. Deemed Execution Date, if any (Month/Day/Year)		A. Securities Acquir Disposed Of (D) (Ins ode (Instr.				5. Amount of Securities Beneficially Owned Following Reported	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
		Code V Amount (A) or Price		Price	Transaction(s) (Instr. 3 and 4)		(1130. 4)			
Common Stock	10/12/2022		G		1,000	D	\$0.00	1,441,342	D	
Common Stock	03/09/2023		G		1,000	D	\$0.00	1,440,342	D	
Common Stock	08/20/2024		Р		47,000	A	\$13.29 ⁽¹⁾	1,487,342 ⁽²⁾	D	

Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities)

			(0) 1	,	,			• •				,			
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	3A. Deemed Execution Date, if any (Month/Day/Year)	4. Transa Code (8)		of		6. Date Exercisable and Expiration Date (Month/Day/Year)		nd 7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		mount of ecurities Security nderlying (Instr. 5) erivative ecurity (Instr.		10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)
				Code	v	(A)	(D)	Date Exercisable	Expiration Date	Title	Amount or Number of Shares				

Explanation of Responses:

1. This transaction was executed in multiple trades at prices ranging from \$13.03 to \$13.4045. 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.

2. Includes 2,947 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorneyin-Fact 09/2

09/17/2024

** 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.