SEC Form 4

FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION Washington, D.C. 20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL

OMB Number:	3235-0287								
Estimated average burden									
hours per response:	0.5								

to Section 16.	t if no longer subje Form 4 or Form 5 y continue. <i>See</i>).	_	Filed pursuant to Section 16(a) of the Securities Exchange Act of 19 or Section 30(h) of the Investment Company Act of 1940	_	OMB Number Estimated ave hours per res	erage burden
1. Name and Addu Hlavinka Sa (Last) C/O QUANTE	(First) (RIX CORPO	(Middle)	2. Issuer Name and Ticker or Trading Symbol <u>Quanterix Corp</u> [QTRX] 3. Date of Earliest Transaction (Month/Day/Year) 04/01/2022	5. Relationship of I (Check all applicat X Director Officer (g below)	ole)	on(s) to Issuer 10% Owner Other (specify below)
900 MIDDLESEX TURNPIKE (Street) BILLERICA MA		KE 01821 (Zip)	4. If Amendment, Date of Original Filed (Month/Day/Year)	Line) X Form filed	d by One Repo	(Check Applicable rting Person One Reporting
		Table I - Non-De	rivative Securities Acquired, Disposed of, or Ben	eficially Owned		

1. Title of Security (Instr. 3)	2. Transaction Date (Month/Day/Year)	2A. Deemed Execution Date, if any (Month/Day/Year)	3. Transa Code (8)		4. Securities Disposed Of			Securities Beneficially	Form: Direct	7. Nature of Indirect Beneficial Ownership (Instr. 4)
			Code	v	Amount	(A) or (D)	Price			(1130.4)
Common Stock	04/01/2022		Α		451 ⁽¹⁾	Α	\$30.47 ⁽²⁾	15,291(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. 3. Transaction 3A. Deemed 4.			5. Number of Derivative Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)			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 Exercisable	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 first quarter of 2022.

2. Closing price of the Company's common stock on the Nasdaq Global Market on April 1, 2022.

3. Includes 2,668 restricted stock units

Remarks:

/s/ Brian Keane, as Attorney-

in-Fact

04/05/2022

** 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.