FORM 4

Check this box if no longer subject

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							

		Table I - Non-De	rivative Securities Acquired. Disposed of, or Ben	eficially Owned			
(City)	(State)	(Zip)					
(Street) BILLERICA	MA	01821			l by One Repo l by More thar	orting Person n One Reporting	
900 MIDDLES	SEX TURNP	IKE	4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Individual or Join Line)	nt/Group Filin	g (Check Applicable	-
(Last) C/O QUANTE			3. Date of Earliest Transaction (Month/Day/Year) 07/01/2021	below)	ive uue	below)	y
1. Name and Addr Hlavinka Sa		g Person [*]	2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]	5. Relationship of f (Check all applicat X Director Officer (gi	ble)	son(s) to Issuer 10% Owner Other (specify	
to Section 16. F	Form 4 or Form 5 y continue. See		Filed pursuant to Section 16(a) of the Securities Exchange Act of 19: or Section 30(h) of the Investment Company Act of 1940	34	Estimated av hours per res	verage burden sponse: 0.5	;

1. Title of Security (Instr. 3)		2A. Deemed Execution Date, if any (Month/Day/Year)	3. Transa Code (8)	ction	4. Securities Disposed Of			Securities Beneficially	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
			Code	v	Amount	(A) or (D)	Price	Transaction(s) (Instr. 3 and 4)		(1150.4)
Common Stock	07/01/2021		Α		199 ⁽¹⁾	Α	\$59.48 ⁽²⁾	11,660 ⁽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)	3A. Deemed Execution Date, if any (Month/Day/Year)	4. Transa Code (8)		5. Number of		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 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 second quarter of 2021.

2. Closing price of the Company's common stock on the Nasdaq Global Market on July 1, 2021.

3. Includes 5,297 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorney-

in-Fact

07/06/2021

** 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.