FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

	OMB APPROVAL									
	OMB Number:	3235-0287								
- 1	Fatimated average I	aurdon								

0.5

hours per response:

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

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* Hlavinka Sarah E. (Last) (First) (Middle) C/O QUANTERIX CORPORATION					<u>Qu</u>	2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX] 3. Date of Earliest Transaction (Month/Day/Year) 10/01/2019									heck all X [I appli Directo	icable) or r (give title	g Person(s) to I: 10% C Other below	Owner (specify
900 MIDDLESEX TURNPIKE (Street) BILLERICA MA 01821 (City) (State) (Zip)					4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)									6. Individual or Joint/Group Filing (Check Applicable Line) X Form filed by One Reporting Person Form filed by More than One Reporting Person				
		Tabl	e I - No	n-Deriv	ative	Sec	uritie	s Ac	quired	d, Dis	sposed o	f, or	Ben	eficia	lly O	wne	d		
1. Title of Security (Instr. 3) 2. Transacti Date (Month/Day					Execution Date,		Date,	3. Transa Code (8)			4. Securities Acquired (A) o Disposed Of (D) (Instr. 3, 4 a			and 5) See Bei Ow		ially Following	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
						Code	v	Amount	(A) (D)	or	Price	Tr	Reported Transaction(s) (Instr. 3 and 4)			(111341. 4)			
Common Stock 10/01/2						2019					320(1)	A \$21.		\$21.41	4,860 ⁽³⁾		860(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)	vative Conversion Date Execution Date urity or Exercise (Month/Day/Year) if any		on Date,	4. Transaction Code (Instr. 8)		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)				te	or		nount mber	8. Price Derivat Securit (Instr. 5	tive (1)	9. Number o derivative Securities Beneficially Owned Following Reported Transaction((Instr. 4)	Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)	

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 2019, pursuant to an election made in July 2019.
- 2. Closing price of the Company's common stock on the Nasdaq Global Market on October 1, 2019.
- 3. Includes 4,540 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorney-

in-Fact

10/03/2019

** 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.