SEC Form 4	
------------	--

П

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).

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934 30(h) of t

OMB APPROVAL								
OMB Number:	3235-0287							
Estimated average burden								
h	0.5							

1. Name and Address of Reporting I Desjourdy Amee	Person*	2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]	5. Relationship of Reporting Person(s) to Issuer (Check all applicable) Director 10% Owner
(Last) (First) C/O QUANTERIX CORPOR 900 MIDDLESEX TURNPIK	· · · · · · · · · · · · · · · · · · ·	3. Date of Earliest Transaction (Month/Day/Year) 08/19/2019	Chief Human Resource Officer
(Street) BILLERICA MA (City) (State)	01821 (Zip)	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

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)	Transaction Code (Instr.		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 and 5)			5. Amount of Securities Beneficially Owned Following Reported		7. Nature of Indirect Beneficial Ownership (Instr. 4)
			Code	v	Amount	(A) or (D)	Price	Transaction(s) (Instr. 3 and 4)		(11501 4)
Common Stock	08/19/2019		Α		11,834(1)	A	\$ <mark>0</mark>	11,834	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. Numb of Derivati Securiti Acquire (A) or Dispose of (D) (II 3, 4 and	ve es d ed nstr.	6. Date Exerc Expiration Da (Month/Day/N	te of Securities		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				
Stock Option (right to buy)	\$27.14	08/19/2019		A		17,354		(2)	08/19/2029	Common Stock	17,354	\$0	17,354	D	

Explanation of Responses:

1. Represents shares of common stock issuable upon vesting of restricted stock units. The restricted stock units vest as to 25% of the shares on August 19, 2020 and vest as to an additional 2.083% of the shares per month at the end of each month for 36 months thereafter beginning August 31, 2020, subject to certain accelerated vesting provisions upon a sale of the company within 18 months of the grant date. 2. The option vests as to 25% of the shares on August 19, 2020 and vests as to an additional 2.083% of the shares per month at the end of each month for 36 months thereafter beginning August 31, 2020, subject to certain accelerated vesting provisions upon a sale of the company within 18 months of the grant date.

/s/ Brian Keane, Attorney-infact

08/21/2019

Date

** 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.

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