FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington D.C. 20549

wasiiiigton,	D.C. 20549

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

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL							
OMB Number:	3235-0287						
Estimated average burden							
haura nar raananaa							

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

Name and Address of Reporting Person* Mattoon Dawn						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]									tionship of Reporting all applicable) Director Officer (give title		10% (
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE					3. Date of Earliest Transaction (Month/Day/Year) 12/02/2019									belov	N) .	below arch Products)	
(Street) BILLERI (City)	ICA M	A ()1821 Zip)		4. If	4. If Amendment, Date of Original Filed (Month/Day/Year)								5. Indiv _ine) X	<i>'</i>			
		Tabl	e I - N	Non-Deriv	ative	Seci	uritie	s Ac	quire	ed, D	isposed o	f, or E	Benefici	ially	Owne	ed		
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Yo				Execution Date,		3. Transaction Code (Instr. 8)					d 5) Sed Bed Ow		ount of ities ficially d Following	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	7. Nature of Indirect Beneficial Ownership			
									Code	v	Amount	(A) or (D)	Price		Reported Transaction(s) (Instr. 3 and 4)			(Instr. 4)
Common Stock 12/02/20				019	19			S		115(1)	D	\$25.0629(2)		22,356 ⁽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)	Derivative decurity lenstr. 3) Date (Month/Day/Year) Date (Month/Day/Year) Fixed of Derivative Security Date (Month/Day/Year) Execution Date, if any (Month/Day/Year) Month/Day/Year) System of Date (Month/Day/Year)		4. Transa Code (8)		5. Num of Derive Secum Acqui (A) or Dispo of (D) (Instr. and 5	ative rities ired osed	6. Date Exercisable and Expiration Date (Month/Day/Year) Date Expiration Exercisable Date			7. Title and Amount of Securities Underlying Derivative Security (Instr. and 4) Amoun or Numbe of Title Shares		nt er		9. Number of 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. The sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 368 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$24.99 to \$25.13. 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.
- 3. Includes 17,288 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorney-

<u>in-Fact</u>

** Signature of Reporting Person Date

12/03/2019

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.