FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 2004

neck this box if no longer subject to STATE action 16. Form 4 or Form 5	STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP
obligations may continue. See	
Instruction 1(b).	Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934

OMB APPROVAL								
OMB Number:	3235-0287							
Estimated average burden								
haura nar raananaa.	٥٦							

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 below) Sr. VP, Resea		10% (
`						3. Date of Earliest Transaction (Month/Day/Year) 03/02/2020											below arch Products)``
(Street) BILLER (City)	ICA M	Α (01821 (Zip)		4. 11	f Amen	Amendment, Date of Original Filed (Month/Day/Year)								ividual or Joint/Group Filing (Check Applicable Form filed by One Reporting Person Form filed by More than One Reporting Person			
		Tabl	le I - N	lon-Deriv	ative	Sec	uritie	s Ac	quire	ed, D	isposed c	f, or E	Benefici	ally	Owne	ed		
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Ye					Execution Date,		···/	3. Transaction Code (Instr. 8) 4. Securities A Disposed Of (I			Acquire (D) (Inst	d (A) or r. 3, 4 and	Secur Benef Owne		icially d Following	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		Reported Transaction(s) (Instr. 3 and 4)			(111511.4)
Common	ommon Stock 03/02/20				020	20		S		115(1)	D	\$22.92	54 ⁽²⁾	3	3,132(3)	D		
		Та	able II							,	posed of, convertib			•	vned			
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	if any	ion Date, Tra	Transa Code (5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		6. Date Exe Expiration I (Month/Day			7. Title and Amount of Securities Underlying Derivative Security (Instr and 4)			vative ırity r. 5)	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 \$22.90 to \$22.99. 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 27,327 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorney-03/04/2020 in-Fact

** 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.