FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

washington,	D.C. 20549

Check this box if no longer subject to	STATEMENT OF CHA
Section 16. Form 4 or Form 5	
obligations may continue. See	
Instruction 1(b).	Filed pursuant to Sect

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL

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

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 (Last) (First) (Middle)							Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX] 3. Date of Earliest Transaction (Month/Day/Year) 02/08/2021									ationship of Reporting Pe k all applicable) Director Officer (give title below) Sr. VP, Research			vner specify
C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE						V2.00, 2021													
(Street) BILLER (City)			01821 (Zip)		_ 4. lt	f Ame	ndment,	Date of Original Filed (Month/Day/Year)							e) X Form	ual or Joint/Group Filing (Check Applicable Form filed by One Reporting Person Form filed by More than One Reporting Person			
		Tab	le I - No	n-Deriv	vative	e Se	curitie	s Ac	quired,	Dis	posed o	of, or	Ben	eficial	ly Owne	ł			
1. Title of Security (Instr. 3) 2. Transac Date (Month/Date)						ar) E	A. Deemed xecution Date, any Month/Day/Year)		Code (Instr.		4. Securities Acquired (A) Disposed Of (D) (Instr. 3, 4 5)				Benefic	es ally Following	Form (D) o	n: Direct r Indirect istr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)
						Code	v	Amount	nount (A		Price	Transac (Instr. 3	tion(s)			(1130.4)			
Common Stock 02/08/						2021			A		3,679	^{'9(1)} A \$		\$0.00	34,	34,513 ⁽²⁾		D	
		T	able II -						uired, C s, optior						Owned				
	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	3A. Deemed Execution E if any (Month/Day	Date,	4. Transa Code (8)		of I		Expiration	s. Date Exercisal Expiration Date Month/Day/Year		Amour Securi Under Deriva	7. Title and Amount of Securities Underlying Derivative Securit (Instr. 3 and 4)		8. Price of Derivative Security (Instr. 5)	9. Number derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4)	e S Illy	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)
					Code	v	(A) (D)		Date Exercisab		expiration Date	Title	0 0	Amount or Number of Shares					
Stock Option (right to	\$77.79	02/08/2021			A		8,585		(3)	0	2/08/2031	Comm		8,585	\$0.00	8,585	5	D	

Explanation of Responses:

- 1. Represents shares of stock issuable on vesting of restricted stock units. The restricted stock units vest as to 25% of the shares on February 8, 2022, and the remaining 75% vest in 36 equal monthly installments at the end of each month beginning on February 28, 2022.
- 2. Includes 23,501 restricted stock units.
- 3. The option vests as to 25% of the shares on February 8, 2022, and the remaining 75% vest in 36 equal monthly installments at the end of each month beginning on February 28, 2022.

Remarks:

<u>/s/ Brian Keane, as Attorney-in-Fact</u>

02/10/2021

** 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.