FORM 4

Check this box if no longer subject to Section 16. Form 4 or Form 5

UNITED STATES SECURITIES AND EXCHANGE COMMISSION Washington, D.C. 20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL

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

Instruction 1(b)	ly continue. See).		Filed pur	rsuant to Section 16	hours per response:		0.5				
	,.					Company Act of 1940	-				
1. Name and Addr Mattoon Da		ng Person*		Issuer Name and Juanterix Cor		0 ,		ationship of R k all applicabl Director Officer (giv	le)	ng Person(s) to 10% C Other	
(Last) C/O QUANTE 900 MIDDLES				Date of Earliest Tra 6/01/2021	ansaction (Mo	nth/Day/Year)		below) Sr. VP, Di		below) biagnostics)
			4.	If Amendment, Da	te of Original I	Filed (Month/Day/Year)	6. Individual or Joint/Group Filing (Check Applicable Line)				
(Street) BILLERICA MA 018		01821					X			e Reporting Per re than One Rep	
(City)	(State)	(Zip)									
		Table I - Non-De	erivativ	e Securities A	cquired, D	isposed of, or Bene	ficially	v Owned			
				an Desmand				E	- 6	A Aumanahim	7. 1

1. Title of Security (Instr. 3)	2. Transaction Date (Month/Day/Year)	2A. Deemed Execution Date, if any (Month/Day/Year)	3. Transa Code (8)				5. Amount of Securities Beneficially Owned Following Reported	Form: Direct (D) or Indirect (I)	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
			Code	v	Amount	(A) or (D)	Price	Transaction(s) (Instr. 3 and 4)	(Instr. 4)	(instr. 4)
Common Stock	06/01/2021		S		243(1)	D	\$48.3181 ⁽²⁾	32,733 ⁽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)	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)		n of		6. Date Exercisable and Expiration Date (Month/Day/Year)		7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		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				

Explanation of Responses:

1. The sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 750 RSUs.

2. This transaction was executed in multiple trades at prices ranging from \$48.00 to \$48.75. 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 17,716 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorney-06/02/2021 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.