

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

**STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP**

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

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

1. Name and Address of Reporting Person* <u>DEKKERS MARIJN E</u>  (Last) (First) (Middle) <u>C/O QUANTERIX CORPORATION,</u> <u>900 MIDDLESEX TURNPIKE</u>  (Street) <u>BILLERICA MA 01821</u>  (City) (State) (Zip)	2. Issuer Name and Ticker or Trading Symbol <u>Quanterix Corp [ QTRX ]</u>	5. Relationship of Reporting Person(s) to Issuer (Check all applicable) <input checked="" type="checkbox"/> Director 10% Owner Officer (give title below) Other (specify below)
	3. Date of Earliest Transaction (Month/Day/Year) <u>01/11/2021</u>	
	4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Individual or Joint/Group Filing (Check Applicable Line) <input checked="" type="checkbox"/> 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)	3. Transaction Code (Instr. 8)		4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 and 5)			5. Amount of Securities Beneficially Owned Following Reported Transaction(s) (Instr. 3 and 4)	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			
Common Stock	06/04/2021		M		1,951	A	\$8.16	11,031 <sup>(1)</sup>	D	
Common Stock	06/04/2021		M		3,168	A	\$23.12	14,199 <sup>(1)</sup>	D	
Common Stock	06/04/2021		s <sup>(2)</sup>		7,610	D	\$55.0483 <sup>(3)</sup>	6,589 <sup>(1)</sup>	D	
Common Stock	06/07/2021		M		4,732	A	\$23.12	11,321 <sup>(1)</sup>	D	
Common Stock	06/07/2021		s <sup>(2)</sup>		4,360	D	\$55.5768 <sup>(4)</sup>	6,961 <sup>(1)</sup>	D	
Common Stock	06/07/2021		s <sup>(2)</sup>		4,691	D	\$56.2449 <sup>(5)</sup>	2,270 <sup>(1)</sup>	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. Transaction Code (Instr. 8)		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		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
Stock Option (right to buy)	\$8.16	06/04/2021		M		1,951		<sup>(6)</sup>	03/31/2027	Common Stock	1,951	\$0.00	0.00	D		
Stock Option (Right to Buy)	\$23.12	06/04/2021		M		3,168			12/31/2020	01/02/2030	Common Stock	3,168	\$0.00	4,732	D	
Stock Option (Right to Buy)	\$23.12	06/07/2021		M		4,732			12/31/2020	01/02/2030	Common Stock	4,732	\$0.00	0.00	D	

**Explanation of Responses:**

- Includes 2,270 restricted stock units.
- The sale of shares was effected pursuant to a Rule 10b5-1 trading plan adopted on March 12, 2021.
- This transaction was executed in multiple trades at prices ranging from \$55.00 to \$55.11. 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.
- This transaction was executed in multiple trades at prices ranging from \$55.17 to \$56.16. 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.
- This transaction was executed in multiple trades at prices ranging from \$56.17 to \$56.41. 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.
- The option vests as to 25% of the shares on March 13, 2018 and vests as to an additional 2.083% of the shares per month thereafter until February 28, 2021.

**Remarks:**

/s/ Brian Keane, as Attorney-in-Fact 06/08/2021

\*\* 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.

