SEC Form 4	
------------	--

П

FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

OMB APPROVAL
OMB Number: 3235-0287

0.5

Estimated average burden

hours per response:

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

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

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

	of Section 30(n) of the investment Company Act of 1940						
1. Name and Address of Reporting Person*	2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]	5. Relationship of Reporting Person(s) to Issuer (Check all applicable)					
DEKKERS MARIJN E	Stamiterus cosh (duan)	X Director 10% Owner					
(Last) (First) (Middle) C/O QUANTERIX CORPORATION,	3. Date of Earliest Transaction (Month/Day/Year) 01/11/2021	Officer (give title Other (specify below) below)					
900 MIDDLESEX TURNPIKE	4. If Amendment, Date of Original Filed (Month/Day/Year)	6. Individual or Joint/Group Filing (Check Applicable Line)					
(Street)		X Form filed by One Reporting Person					
BILLERICA MA 01821		Form filed by More than One Reporting Person					
(City) (State) (Zip)							

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)					5. Amount of Securities Beneficially Owned 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	01/11/2021		М		7,900	A	\$18.24	16,980	D	
Common Stock	01/11/2021		М		7,900	A	\$14.8	24,880	D	
Common Stock	01/11/2021		М		29,162	A	\$8.16	54,042	D	
Common Stock	01/11/2021		S ⁽¹⁾		7,200	D	\$47.1656(2)	46,842	D	
Common Stock	01/11/2021		S ⁽¹⁾		22,830	D	\$47.8651 ⁽³⁾	24,012	D	
Common Stock	01/11/2021		S ⁽¹⁾		2,333	D	\$49.1564(4)	21,679	D	
Common Stock	01/11/2021		S ⁽¹⁾		4,400	D	\$50.3767(5)	17,279	D	
Common Stock	01/11/2021		S ⁽¹⁾		8,199	D	\$51.0821(6)	9,080 ⁽⁷⁾	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)		of Deri Sec Acq (A) (Disp of (E	umber vative urities uired or oosed 0) (Instr. and 5)	6. Date Exerc Expiration Da (Month/Day/Y	ate	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)	\$18.24	01/11/2021		М			7,900	12/31/2019	01/02/2029	Common Stock	7,900	\$0.00	0.00	D	
Stock Option (Right to Buy)	\$14.8	01/11/2021		М			7,900	12/31/2018	08/07/2028	Common Stock	7,900	\$0.00	0.00	D	
Stock Option (right to buy)	\$8.16	01/11/2021		М			29,162	(8)	03/31/2027	Common Stock	29,162	\$0.00	1,951	D	

Explanation of Responses:

1. The sale of shares was effected pursuant to a Rule 10b5-1 trading plan adopted on December 9, 2020.

2. This transaction was executed in multiple trades at prices ranging from \$46.47 to \$47.46. 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. This transaction was executed in multiple trades at prices ranging from \$47.47 to \$48.43. 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.

4. This transaction was executed in multiple trades at prices ranging from \$48.74 to \$49.70. 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.

5. This transaction was executed in multiple trades at prices ranging from \$49.79 to \$50.78. 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.

6. This transaction was executed in multiple trades at prices ranging from \$50.79 to \$51.64. 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.

7. Includes 2,270 restricted stock units.

8. 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:

** 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.