FORM 4

## **UNITED STATES SECURITIES AND EXCHANGE COMMISSION**

Vashington,	D.C.	20549	

## STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

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

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

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

1. Name and Address of Reporting Person*  DEKKERS MARIJN E					2. Issuer Name <b>and</b> Ticker or Trading Symbol Quanterix Corp [ QTRX ]								Relationship heck all appli X Direct	cable) or	g Pers	10% Ov	vner	
(Last)	`	irst)	(Middle) N,			3. Date of Earliest Transaction (Month/Day/Year) 01/04/2021								Office below	(give title		Other (s below)	pecify
900 MIDDLESEX TURNPIKE					4. 1	If Amendment, Date of Original Filed (Month/Day/Year)							6. Individual or Joint/Group Filing (Check Applicable					
(Street)	ICA M	Α	01821										Form	Form filed by One Reporting Person Form filed by More than One Reporting Person				
(City)	(S	tate)	(Zip)															
	Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																	
Date					action 2A. Deemed Execution Date, if any (Month/Day/Year)		Code (Instr.   5)			Benefic	es Form ally (D) of following (I) (Ir		: Direct r Indirect str. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)				
							Code	v	Amount	(A) or (D)	Price	Tropos	tion(s)			(msu. 4)		
Common Stock 01/04			4/202	21			A 2,270 <sup>(1)</sup> A \$		\$0.	00 9,	9,080(2)		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 D if any (Month/Day/	ate, 1		ansaction Derivative E		6. Date Exercisable and Expiration Date (Month/Day/Year)  (Month/Day/Year)  T. Title and Amount of Securities Underlying Derivative Sec (Instr. 3 and 4)			of s ng e Securit	8. Price of Derivative Security (Instr. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s)	e S Illy	Ownership Form: Direct (D)	11. Nature of Indirect Beneficial Ownership (Instr. 4)		
					Code	v	(A)	(D)	Date Exercisal		Expiration Date	Title	Amour or Number of Shares	r	(Instr. 4)			
Stock Option (Right to Buy)	\$44.1	01/04/2021			A		7,900 <sup>(3)</sup>		12/31/20	21	01/04/2031	Common Stock	7,90	\$0.00	7,900		D	

## **Explanation of Responses:**

- 1. Represents restricted stock units that vest as to 100% of the shares on December 31, 2021 granted pursuant to the Quanterix Corporation 2018 Non-Employee Director Compensation Policy. Each restricted stock unit represents the right to receive one share of the Issuer's common stock.
- 2. Includes 2,270 restricted stock units.
- 3. Represents non-qualified stock option to purchase 7,900 shares of common stock granted pursuant to the Quanterix Corporation 2018 Non-Employee Director Compensation Policy.

## Remarks:

/s/ Brian Keane, as Attorney-in-01/05/2021

**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.