FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington,	D.C.	20549	

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

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

				_		_		_		_				_				_	
Name and Address of Reporting Person* Madaus Martin D				2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]								5. Relationship of Reporting Person(s) to Issuer (Check all applicable)							
Widdaus Wartin D													X Director			10% Ow			
(Last) (First) (Middle) C/O QUANTERIX CORPORATION,					3. Date of Earliest Transaction (Month/Day/Year) 11/04/2020									Officer below)	(give title	•	Other below	(specify)	
-		ΓURNPIKE	,		\vdash														
					_ 4. If	f Amer	ndmer	nt, Date	of Origin	al File	ed (Month/D	ay/Year)		Indiv ne)	ridual or 3	Joint/Grou	up Filing	g (Check A	pplicable
(Street)														X	Form f	iled by Or	ne Repo	orting Pers	son
BILLER	ICA M	ÍΑ	01821														ore thar	n One Rep	orting
(City)	(S	tate)	(Zip)		-										Persor	1			
(0.0)	()			on Dori	vative	. 500	uviti	ioc Ao	auiroc		sposed o	of or Bo	noficia	ally (Ownoo				
1. Title of	Security (Ins		16 1 - 140	2. Transa		_	Deem		3.	i, Di	·	es Acquired		Ť	5. Amoun		6. Own	nership	7. Nature of
1. The of security (man. 5)		Date (Month/Day/Ye) if a	Execution Date, if any (Month/Day/Year)		Transaction Code (Instr. 8)		Disposed Of (D) (Inst		. 3, 4 and		Securities Beneficially Owned Following	lly	Form: Direct (D) or Indirect (I) (Instr. 4)	Direct Indirect tr. 4)	Indirect Beneficial Ownership		
									Code	v	Amount	(A) or (D)	Price	1	Reported Fransactio (Instr. 3 au				(Instr. 4)
Common	Stock			11/04	/2020				M		1,600	A	\$2.2	8	13,31	16 ⁽¹⁾	D		
Common	Stock			11/04	/2020				S		1,600(2) D	\$39.	4	11,716(1)]	D	
Common	Stock			11/05	/2020				M		1,372	A	\$2.2	8	13,08	38(1)	D		
Common	Stock			11/05	/2020				S		1,372(2) D	\$42.4	1 5	11,71	16 ⁽¹⁾	D		
Common	Stock														164,	966	I See Footnot		See Footnote ⁽³⁾
		Т	able II					-	-		oosed of converti			-	wned				
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	3A. Deer Execution if any (Month/I	med	4. Transa Code (8)	ction	5. No of Deriv Secu Acqu (A) of Disp of (D	vative urities uired or oosed o)	· •	Exerci	sable and te	7. Title an Amount o Securities Underlyin Derivative (Instr. 3 au	d f s g Security	8. De Se (In	Price of erivative ecurity estr. 5)	9. Number derivative Securities Beneficial Owned Followin Reported Transact (Instr. 4)	re es ally g	10. Ownershi Form: Direct (D) or Indirec (I) (Instr. 4	Beneficial Ownership t (Instr. 4)
					Code	v	(A)	(D)	Date Exercisa	ıble	Expiration Date	Title	Amount or Number of Shares						
Stock Option (Right to Buy)	\$2.28	11/04/2020			М			1,600	(4)		10/31/2021	Common Stock	1,600		\$0.00	1,37	72	D	
Stock Option (Right to	\$2.28	11/05/2020			M			1,372	(4)		10/31/2021	Common Stock	1,372		\$0.00	0.00	0	D	

Explanation of Responses:

- 1. Includes 2,270 Restricted Stock Units.
- 2. All sales reported in this Form 4 were effected pursuant to a Rule 10b5-1 trading plan adopted by the Reporting Person on September 9, 2019.
- 3. Held by a trust for which the Reporting Person is a trust advisor who shares voting and investment control.
- 4. All options are vested and exercisable.

Remarks:

/s/ Brian Keane, Attorney-in-

11/06/2020

Fact

** Signature of Reporting Person

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.