FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

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

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* Hlavinka Sarah E.							Name and terix Co				Symbol	(Ch	eck all applic	all applicable) Director		p Person(s) to Issuer 10% Owner		
`	ANTERIX	irst) CORPORATIOΩ ΓURNPIKE	(Middle) V			3. Date of Earliest Transaction (Month/Day/Year) 01/03/2022								Officer below)	Officer (give title below)		Other (s below)	specify
(Street) BILLER (City)	ICA M	IA	01821 (Zip)		4. 1	If Ame	endment, C	ate o	f Origina	al Filed	d (Month/Day	Line	ndividual or Joint/Group Filing (Check Applicable e) X Form filed by One Reporting Person Form filed by More than One Reporting Person					
		Tal	ole I - No	n-Deri	vativ	e Se	curities	Ac	quired	l, Dis	sposed of	f, or Ber	neficiall	y Owned				
1. Title of Security (Instr. 3) 2. Transac Date (Month/Da					Executy/Year) if any		A. Deemed Execution Date, f any Month/Day/Year)		ction Instr.	4. Securities Acquired (A) of Disposed Of (D) (Instr. 3, 4 a			Benefici	es ally Following	Form (D) o	: Direct r Indirect str. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
									Code	v	Amount	(A) or (D)	Price	Transac (Instr. 3	tion(s)			
Common Stock				01/03	3/2022	2			A		2,668(1)	A	\$0.00	14,5	560 ⁽²⁾		D	
Common Stock 01/0				01/03	3/2022	2022			A		280(3)	A	\$42.41	14,8	14,840(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)	Derivative Conversion Date Security or Exercise (Month/Day/Ye		3A. Deemed Execution Date, if any (Month/Day/Year)		4. Transactio Code (Inst 8)				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 derivative Securitie: Beneficia Owned Following Reported Transacti	e s ally	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership t (Instr. 4)
					Code	v	(A)	(D)	Date Exercis	able	Expiration Date	Title	Amount or Number of Shares		(Instr. 4)			
Stock Option (Right to Buy)	\$42.41	01/03/2022					4,002 ⁽⁵⁾		12/31/2	2022	01/03/2032	Common Stock	4,002	\$0.00	4,002	2	D	

Explanation of Responses:

- 1. Represents restricted stock units that vest as to 100% of the shares on December 31, 2022 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 2,668 restricted stock units.
- 3. Consists of stock granted in lieu of cash fees for service on the Company's Board of Directors and committees thereof for the fourth quarter of 2021.
- 4. Closing price of the Company's common stock on the Nasdaq Global Market on January 3, 2022.
- 5. Represents non-qualified stock option to purchase 4,002 shares of common stock granted pursuant to the Quanterix Corporation 2018 Non-Employee Director Compensation Policy.

Remarks:

/s/ Brian Keane, as Attorney-in-Fact 01/05/2022

** 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.