FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington,	D.C.	20549
-------------	------	-------

Check this box if no longer subject to Section 16. Form 4 or Form 5	STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP
obligations may continue. See	
Instruction 1(b).	Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934

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

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* Flynn Karen					2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]										tionship of Reportir all applicable) Director		10% Ov		vner	
_ `		CORPORATION	Middle)			3. Date of Earliest Transaction (Month/Day/Year) 07/01/2022									Officer (give title below)			Other (below)		
900 MIDDLESEX TURNPIKE					4. If Amendment, Date of Original Filed (Month/Day/Year)									6. Individual or Joint/Group Filing (Check Applicable Line)						
(Street) BILLER	ICA M	A 0	1821												X Form filed by One Reporting Person Form filed by More than One Reporting Person					
(City)	(St	ate) (2	Zip)																	
	Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																			
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/				/Year) Executio		cution Date,		3. Transaction Code (Instr. 8)				and 5) Securi Benefi		cially I Following	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	Trans		action(s) 3 and 4)			(msu. 4)		
Common Stock 07/01/2				022			A		164(1)	A	\$16.	6.79(2)		13,391(3)		D				
		Tal	ble II								osed of, convertib				Owne	d				
1. Title of Derivative Security (Instr. 3)	erivative Conversion Date Execution Date, ecurity or Exercise (Month/Day/Year) if any					ansaction of De Se Ac (A) Dis		osed) r. 3, 4	Expiration I (Month/Day		ate	7. Title and Amount of Securities Underlying Derivative Security (Inst 3 and 4)		Der Sec (Ins	Price of ivative curity str. 5)	9. Number derivative Securities Beneficiall Owned Following Reported Transactio (Instr. 4)	y C	Ownershi Form: Direct (D) or Indirect (I) (Instr. 4	Beneficial Ownership (Instr. 4)	
					Code	v	(A)	(D)	Date Exerc	isable	Expiration Date	Title	or Number of Shares							

Explanation of Responses:

- 1. Consists of stock granted in lieu of cash fees for service on the Company's Board of Directors and committees thereof for the second quarter of 2022.
- 2. Closing price of the Company's common stock on the Nasdaq Global Market on July 1, 2022.
- 3. Includes 13,227 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorney-

07/06/2022

in-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.