## 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 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

Name and Address of Reporting Person*     WALT DAVID R			2. Issuer Name <b>and</b> Ticker or Trading Symbol Quanterix Corp [ QTRX ]							eck all applic  X Director	10% Owner		ner						
(Last)	`	irst)	(Middle) N,		3. Date of Earliest Transaction (Month/Day/Year) 01/03/2023							Officer below)	(give title		Other (s below)	pecify			
900 MIDDLESEX TURNPIKE					4. If /	4. If Amendment, Date of Original Filed (Month/Day/Year)							6. Individual or Joint/Group Filing (Check Applicable Line)						
(Street) BILLER	ICA M	ΙA	01821										Lim	X Form fi	,	•	rting Person One Report		
(City)	(S	tate)	(Zip)																
Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																			
Date		2. Transa Date (Month/E	h/Day/Year)		2A. Deemed Execution Date, if any (Month/Day/Year)		3. Transaction Code (Instr. 8)  4. Securities Acquired (A Disposed Of (D) (Instr. 3, 1)			5. Amour Securitie Beneficia Owned F Reported	s Form ally (D) o ollowing (I) (In		: Direct Indirect Istr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)					
								Code	v	Amount	(A) or (D)	Price	Transact (Instr. 3 a	on(s)			iiisu. 4)		
Common Stock 01/03			01/03	3/2023	3			A		8,083(1	) A	\$0.0	0 1,439	1,439,395(2)		D			
Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities)																			
1. Title of Derivative Conversion Date December December Of Execution Date, Of Execution		Co	ansaction Derivative		6. Date Exercisable and Expiration Date (Month/Day/Year)  7. Title and A of Securities Underlying Derivative Se (Instr. 3 and 4		ties ig e Security	Derivative Security	9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s)	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)								
				Co	ode V	,	(A)	(D)	Date Exercisa	ble	Expiration Date	Title	Amount or Number of Shares		(Instr. 4)	J.11(3)			
Stock Option (Right to Buy)	\$12.86	01/03/2023		1	A		12,124 <sup>(3)</sup>		12/31/20	)23	01/03/2033	Common Stock	12,12	\$0.00	12,124	4	D		

## Explanation of Responses:

- 1. Represents restricted stock units that vest as to 100% of the shares on December 31, 2023 granted pursuant to the Quanterix Corporation Amended and Restated 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 8,083 restricted stock units.
- 3. Represents non-qualified stock option to purchase 12,124 shares of common stock granted pursuant to the Quanterix Corporation Amended and Restated 2018 Non-Employee Director Compensation Policy.

## Remarks:

/s/ Brian Keane, as Attorney-in-

**Fact** 

\*\* Signature of Reporting Person Date

01/04/2023

Reminder: Report on a separate line for each class of securities beneficially owned directly or indirectly.

- $^{\star}$  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.