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

1. Name and Address of Reporting Person*  WALT DAVID R					2. Issuer Name <b>and</b> Ticker or Trading Symbol Quanterix Corp [ QTRX ]							(Che	elationship o eck all applic	able)	g Perso	on(s) to Issu 10% Ow			
(Last) (First) (Middle) C/O QUANTERIX CORPORATION,						3. Date of Earliest Transaction (Month/Day/Year) 01/03/2022							Officer below)	er (give title V)		Other (s below)	pecify		
900 MIDDLESEX TURNPIKE					4. 1	4. If Amendment, Date of Original Filed (Month/Day/Year)								6. Individual or Joint/Group Filing (Check Applicable					
(Street)	et) .LERICA MA 01821											Line							
(City)	(S	tate)	(Zip)																
		Tak	ole I - Non	-Deriv	ativ	e Se	curities	Acc	quired,	Dis	posed of	f, or Ber	eficiall	y Owned					
1. Title of Security (Instr. 3)  2. Transa Date (Month/D					Execution Date		Date,	3. Transaction Code (Instr. 8)  4. Securities Acquired Disposed Of (D) (Instr. 5)		d (A) or r. 3, 4 and	5. Amour Securitie Beneficia Owned F Reported	es ally Following		: Direct I r Indirect E str. 4) (	7. Nature of Indirect Beneficial Ownership				
									Code	v	Amount	(A) or (D)	Price	Transact (Instr. 3 a	ion(s)			Instr. 4)	
Common Stock 01/03				3/2022		A		2,668(1	1) <b>A</b>	\$0.00	1,281	1,281,312(2)		D					
			Table II - I								osed of, onvertib			Owned					
1. Title of Derivative Security (Instr. 3)	rivative Conversion Date Execution Date, curity or Exercise (Month/Day/Year) if any		ate, T	ransaction Derivative Securities		e s I (A) sed str.	Expiration Date (Month/Day/Year)		е	7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		8. Price of Derivative Security (Instr. 5)	9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s)		10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership (Instr. 4)			
				c	Code	v	(A)	(D)	Date Exercisal:		Expiration Date	Title	Amount or Number of Shares		(Instr. 4)	on(s)			
Stock Option (Right to Buy)	\$42.41	01/03/2022			A		4,002 <sup>(3)</sup>		12/31/202	22	01/03/2032	Common Stock	4,002	\$0.00	4,002	!	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. 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-01/05/2022

**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.