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								

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* CRANDELL KEITH					2. Issuer Name and Ticker or Trading Symbol Ouanterix Corp [OTRX]								Relationship of Reporting Person(s) to Issuer (Check all applicable)					
CRANDELL KEITH					-`									C Director	r	10% C		ner
(Last) 8755 WE	,	irst) NS ROAD, SUI	(Middle) TE 1025		3. Date of Earliest Transa 01/03/2023				action (Month/Day/Year)					below)			Other (s below)	
(Street)	O IL		60631		_ 4.								Line) 【 Form fi	ual or Joint/Group Filing (Check Applicable Form filed by One Reporting Person Form filed by More than One Reporting			
(City)	(S	tate)	(Zip)											Person		e triari	One Report	irig
	Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																	
Date			Date	nsaction h/Day/Year)		2A. Deemed Execution Date, if any (Month/Day/Year)		3. Transaction Code (Instr. 8)		4. Securities Acquired (A) o Disposed Of (D) (Instr. 3, 4 a		(A) or 3, 4 and 5)	5. Amour Securitie Beneficia Owned F Reported	es Formally (D) (Sollowing (I) (I		: Direct Indirect Str. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)	
									Code	v	Amount	(A) or (D)	Price	Transact (Instr. 3 a	ion(s)			(111501.4)
Common Stock 01/03				3/202	3			A		8,083(1)	A	\$0.00	129,	051(2)		D		
Common Stock 01/03/				3/202	2023		A		1,015(3)	A	\$12.86	⁴⁾ 130,	066(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)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	3A. Deemed Execution D if any (Month/Day/	Date,	4. Transa Code (8)				6. Date Exercisable and Expiration Date (Month/Day/Year)		ate	7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4)		8. Price of Derivative Security (Instr. 5)	9. Number derivative Securities Beneficial Owned Following Reported Transactio	e S Illy	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	Beneficial Ownership (Instr. 4)
					Code	v	(A)	(D)	Date Exercis	sable	Expiration Date	Title	Amount or Number of Shares	er (Insti		J.1(J)		
Stock Option (Right to Buy)	\$12.86	01/03/2023			A		12,124 ⁽⁵⁾		12/31/2	2023	01/03/2033	Common Stock	12,124	\$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. 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 2022.
- 4. Closing price of the Company's common stock on the Nasdaq Global Market on January 3, 2023.
- 5. 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-01/04/2023 Fact for Keith L. Crandell

** 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.