FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

Washington, D.C. 20049	OMB APPROVAL			
STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP	OMB Number:	3235-0287		
THE THE TOTAL OF T	Estimated average bur	den		
Filed pursuant to Section 16(a) of the Securities Eychange Act of 1034	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* <u>CRANDELL KEITH</u>				2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]									all app Direc	tor	ng Per X	10% O	wner	
(Last) (First) (Middle) 8755 WEST HIGGINS ROAD, SUITE 1025			3. Date of Earliest Transaction (Month/Day/Year) 07/01/2022										Office below	er (give title		Other (below)	specify	
(Street)			0631		4. If #	4. If Amendment, Date of Original Filed (Month/Day/Year)							i. Indivine)	-/				
(City)	(Sta		Zip)															
Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned																		
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/N				Execution Date,		,	3. Transaction Code (Instr. 8) 4. Securities Acquired (A) Disposed Of (D) (Instr. 3, 4)				and 5) Securiti Benefic		ties For cially (D) Following (I) (wnership n: Direct or Indirect nstr. 4)	7. Nature of Indirect Beneficial Ownership (Instr. 4)		
								Code	v	Amount	(A) or (D)	Price	Transa		ction(s) 3 and 4)			(IIISU. 4)
Common Stock 07/01/2			07/01/20)22			A		893(1)	A	\$16.7	79 ⁽²⁾ 119,		,642(3)		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 Execution Date, Security or Exercise (Month/Day/Year) if any		4. Transa Code (8)	(Instr.	nstr. Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		6. Date Exercisable and Expiration Date (Month/Day/Year) Date Expiration		7. Title and Amount of Securities Underlying Derivative Security (Inst: 3 and 4) Amour or Numbe of		Derivative d Security S B C F F T (I		9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s) (Instr. 4)		10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)		

Explanation of Responses:

- 1. Consists of stock granted in lieu of eash 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 2,668 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorneyin-Fact for Keith L. Crandell

07/06/2022

0.5

** 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.