FORM 4

Check this box if no longer subject to Section 16. Form 4 or Form 5 obligations may continue. See

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									

	tion 1(b).	nuc. See		Filed							ities Exchanç ompany Act o		of 1934			nours	per re	esponse:	0.5	
1. Name and Address of Reporting Person* HRUSOVSKY E KEVIN					2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]									5. Relationship of Reporting Person(s) to Issuer (Check all applicable) X Director 10% Owner						
(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE					3. Date of Earliest Transaction (Month/Day/Year) 06/23/2020									X	Office	er (give title	Other (specify below)			
(Street) BILLER (City)	ICA M	A 0	1821 Zip)		4. If Amendment, Date of Original Filed (Month/Day/Year)								Indivi ne) X	lividual or Joint/Group Filing (Check Applicable Form filed by One Reporting Person Form filed by More than One Reporting Person						
		Table	I - N	on-Deriva	tive :	Secu	rities	Ac	quire	d, Di	sposed of	f, or E	Benefic	ially	Own	ed				
1. Title of Security (Instr. 3) 2. Transactio Date (Month/Day/N					Execution Date			´				rities Acquired (A) o ed Of (D) (Instr. 3, 4		and 5) S B O		Securities Beneficially		n: Direct or Indirect nstr. 4)	7. Nature of Indirect Beneficial Ownership	
								Ī	Code	v	Amount	(A) or (D)	Price	Transa		action(s) 3 and 4)			(Instr. 4)	
Common Stock 06/23/20					20				S ⁽¹⁾		5,000	D	\$27.06	1,074,637 ⁽³⁾		74,637 ⁽³⁾		D		
		Tal	ble II	- Derivati (e.g., pເ							oosed of, convertib				wne	d				
1. Title of Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	Exec if any	Deemed ution Date, / uth/Day/Year)	4. Transaction Code (Instr. 8)		5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5)		Expir	te Exer ration C th/Day		and 7. Title and Amount of Securities Underlying Derivative Security (Ins 3 and 4)		Deriv Secu	Price of rivative curity str. 5)	9. Number derivative Securities Beneficiall Owned Following Reported Transaction (Instr. 4)	у	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirec Beneficial Ownershi (Instr. 4)	
		Code	v	(A)	(D)	Date Exerc	cisable	Expiration Date	Title	Amount or Number of Shares										

Explanation of Responses:

- 1. The sale of shares was effected pursuant to a Rule 10b5-1 trading plan adopted on December 5, 2019.
- 2. This transaction was executed in multiple trades at prices ranging from \$26.87 to \$27.48. The price reported reflects the weighted average sale price. The reporting person undertakes to provide upon request to the SEC staff, the issuer or any security holder of the issuer full information regarding the number of shares and prices at which the transaction was effected.
- 3. Includes 185,255 restricted stock units and 39,815 shares of common stock that are subject to performance based vesting,

Remarks:

/s/ Brian Keane, as Attorney-06/25/2020 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.