FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APP	OMB APPROVAL										
OMB Number:	3235-028										

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

OMB Number:	3235-0287
Estimated average bure	den
hours per response:	0.5

Name and Address of Reporting Person* Duffy David C.					Quanterix Corp [QTRX]									(Che	eck all appli Directo	cable)	g Person(s) to Issu 10% Ow Other (sp		ner
(Last) C/O QUA		3. Date of Earliest Transaction (Month/Day/Year) 09/16/2019										below) below) SVP R&D and CTO							
(Street) LEXING	GTON M	IA .	02421 (Zip)		4. If	Ame	nendment, Date of Original Filed (Month/Day/Year)							Line) K Form	or Joint/Group Filing (Check App m filed by One Reporting Person m filed by More than One Repor son			n
		Tab	le I - No	n-Deriv	ative	Sec	curiti	ies Ac	quired	Dis	posed o	of, or	Bene	ficiall	y Owned	t k			
1. Title of Security (Instr. 3)			2. Transaction Date (Month/Day/Year)		Execution Date,		, Transaction Dis Code (Instr. 5)		Disposed	I. Securities Acquired (A) Disposed Of (D) (Instr. 3, 4			5. Amou Securiti Benefic Owned Reporte	es ially Following	6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)		7. Nature of Indirect Beneficial Ownership		
									Code	v	Amount	A) (1)	A) or D)	Price	Transac (Instr. 3	ction(s)			(Instr. 4)
Common	Stock			09/16	/2019				М		1,667	7	A	\$1.93	46	46,427 D			
Common	Stock			09/16	/2019				S		1,667	(1)	D	\$25.6	8 44,760 ⁽²⁾ D				
		7	able II -								osed of converti				Owned				
Derivative Security (Instr. 3)	2. Conversion or Exercise Price of Derivative Security	3. Transaction Date (Month/Day/Year)	3A. Deem Executior if any (Month/Da	Date,		ransaction ode (Instr.		n of E		6. Date Exercisa Expiration Date (Month/Day/Yea		7. Title and Amount of Securities Underlying Derivative Sec (Instr. 3 and 4)			3. Price of Derivative Security Instr. 5)	9. Number derivative Securities Beneficial Owned Following Reported Transactio (Instr. 4)	ly	10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4)	11. Nature of Indirect Beneficial Ownership (Instr. 4)
					Code	v	(A)	(D)	Date Exercisal		Expiration Date	Title	or No of	umber					
Stock Option (Right to	\$1.93	09/16/2019			M			1,667	(3)	(07/06/2021	Comm		.,667	\$0.00	6,942		D	

Explanation of Responses:

- 1. The sale of shares reported on this Form 4 was made pursuant to a Rule 10b5-1 trading plan adopted on May 16, 2019.
- 2. Includes 9,291 restricted stock units.
- 3. Option is fully vested.

Remarks:

Buy)

/s/ Brian Keane, Attorney-in-

09/17/2019

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.