Instruction 1(b)

FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

D.C. 20549
D.C. 20549

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

OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL 3235-0287 Estimated average burden hours per response: 0.5

Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934 or Section 30(h) of the Investment Company Act of 1940

(Last) (First) (Middle) C/O QUANTERIX CORPORATION, 900 MIDDLESEX TURNPIKE (Street) BILLERICA MA 01821 Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned Following (Month/Day/Year) (Month/Day/Year) (Month/Day/Year) (Month/Day/Year) (Month/Day/Year) (State) (Zip) Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned Following (Month/Day/Year) (Month/Day/Year)	Name and Address of Reporting Person* Mattoon Dawn						2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX]								Check	all app	o of Reportir dicable) tor er (give title	ng Pe	erson(s) to Is 10% O Other (wner
City	C/O QUANTERIX CORPORATION,														X	below) below)			opeoy	
1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Year) 2. A. Deemed Execution Date, if any (Month/Day/Year) 3. Transaction (Code (Instr. 4)) 4. Amount (A) or Disposed Of (D) (Instr. 3, 4 and 5) 5. Amount of Securities Deneficially Owned Following Reported Transaction(s) (Instr. 4) 4. Deemed Execution Date, if any (Month/Day/Year) 5. Amount of Securities Deneficially Owned (Instr. 4) 5. Amount of Disposed Of (D) (Instr. 3, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 3, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 3, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 3, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 3, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 3, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 3, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 3, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 3, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 3, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 3, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 3, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 4, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 4, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 4, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 4, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 4, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 4, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 4, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 4, 4 and 5) 5. Amount of Disposed Of (D) (Instr. 4, 4 and 5) 5. Amount of Disposed Of (D) (D) (Instr. 4, 4 and 5) 5. Amount of Disposed Of (D) (D) (Instr. 4, 4 and 5) 5. Amount of Disposed Of (D) (D) (Instr. 4, 4 and 5) 5. Amount of Disposed Of (D)	(Street) BILLERICA MA 01821						4. If Amendment, Date of Original Filed (Month/Day/Year)								.ine)	X Form filed by One Reporting Person Form filed by More than One Reporting				
Date (Month/Day/Year) Fixed times (Month/Day/Year) Month/Day/Year) Date (Month/Day/Year) Date (Month/Day/Year) Disposed of (D) (Instr. 3, 4 and 5) Disposed of (D) (Instr. 3, 4 and 5) Securities (D) or Indirect (D) or Indir			Table	I - I	Non-Deriva	tive S	Secui	rities	Ac	quir	ed, D	isposed (of, or	Benefic	ially	Own	ed			
Common Stock O1/03/2022 S 355(1) D \$41.9066(2) 37,989(3) D	Date			Execution Date, ar) if any		· [Transaction Code (Instr.		Disposed Of (D) (Instr. 3, 4 a			and 5) Secu Bene Owne		rities F eficially (ed Following I		Form: Direct (D) or Indirect (I)	of Indirect Beneficial Ownership			
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) 1. Title of Derivative Security (Instr. 3) 2. Conversion or Exercise Price of Derivative Security (Instr. 3) 1. Title of Derivative Security (Instr. 3) 2. Conversion or Exercise (Month/Day/Year) 3. Transaction Date (Execution Date, if any (Month/Day/Year) 4. Transaction Code (Instr. 8) 5. Number of Exercisable and Expiration Date (Month/Day/Year) 6. Date Exercisable and Expiration Date (Month/Day/Year) 7. Title and Amount of Securities Security (Instr. 5) 9. Number of derivative Security (Instr. 5) 9. Ownership Form: Direct (D) Ownership Pormic (Instr. 4) 11. Nature of Indirect (I) (Instr. 4) 12. Ownership Form: Direct (D) Ownership Security (Instr. 4) 13. Transaction Date (Month/Day/Year) 14. Transaction Date (Month/Day/Year) 15. Number of Derivative Security (Instr. 3) 16. Date Exercisable and Expiration Date (Month/Day/Year) 17. Title and Amount of Securities Underlying Derivative Security (Instr. 3) 18. Price of Derivative Security (Instr. 3) 19. Number of Derivative Security (Instr. 4) 10. Ownership Form: Direct (D) Ownership Form: Dire									١	Code	v	Amount	(A) or (D)	Price		Transa	action(s)	(1113		(111341.4)
(e.g., puts, calls, warrants, options, convertible securities) 1. Title of Derivative Security (Instr. 3) 1. Title of Derivative Security (Instr. 3) 1. Title of Derivative Security (Instr. 3) 2. Conversion or Exercise (Month/Day/Year) 2. Conversion On Exercise (Month/Day/Year) 3. Transaction Date (Execution Date, if any (Month/Day/Year) 4. Transaction Code (Instr. 8) 4. Transaction Date (Month/Day/Year) 5. Number of Expiration Date (Month/Day/Year) 6. Date Exercisable and Expiration Date (Month/Day/Year) 6. Date Exercisable and Expiration Date (Month/Day/Year) 8. Price of Derivative Security (Instr. 5) 9. Number of derivative Security (Instr. 5) 9. Number of Derivative Security (Instr. 4) 10. Ownership Form: Direct (D) Ownership Porman Direct (D) Ownership Porman Date (Instr. 4) 11. Nature (Month/Day/Year) 12. Amount of Securities (Instr. 4) 13. Transaction Date (Month/Day/Year) 14. Transaction Date (Month/Day/Year) 15. Number of Derivative Security (Instr. 3) 16. Date Exercisable and Expiration Date (Month/Day/Year) 16. Date Exercisable and Expiration Date (Month/Day/Year) 17. Title and Amount of Securities Underlying Derivative Security (Instr. 3) 18. Price of Derivative Security (Instr. 3) 19. Number of derivative Security (Instr. 3) 10. Ownership Form: Direct (D) Own	Common Stock 01/03/202				2	s 355 ⁽¹⁾		D	\$41.9066 ⁽²⁾		37,989 ⁽³⁾			D						
Derivative Security (Instr. 3) Price of Derivative Security Security Security Price of Derivative Security Secu																				
ı ı ı ı ı lılılıpate lexpiration lot i l l	Derivative Security	Conversion or Exercise Price of Derivative	Date	Exe if ar	cution Date, ny	Transa Code (Transaction Code (Instr. 8) Berivativ Securitie Acquirer (A) or Dispose of (D) (Instr. 3,			Expiration Date (Month/Day/Year) sid			Amo Secu Unde Deriv Secu 3 and	unt of prities erlying vative prity (Instr. d 4)	Deriva Securi (Instr.		derivative Securities Beneficially Owned Following Reported Transaction(s)		Ownership Form: Direct (D) or Indirect	11. Nature of Indirect Beneficial Ownership (Instr. 4)

Explanation of Responses:

- 1. The sales of shares reported on this Form 4 were made solely to cover tax obligations upon the vesting of 750 RSUs.
- 2. This transaction was executed in multiple trades at prices ranging from \$41.71 to \$42.16. 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 12,466 restricted stock units.

Remarks:

/s/ Brian Keane, as Attorneyin-Fact

01/05/2022

** 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.