Check

FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

| Washington, | D.C. 20549 | |
|-------------|------------|--|
|-------------|------------|--|

| Check this box if no longer subject |
|-------------------------------------|
| to Section 16. Form 4 or Form 5 |
| obligations may continue. See |
| Instruction 1(h) |

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

| | OMB APPROVAL | | | | | | | | | |
|---|--------------------------|-----|--|--|--|--|--|--|--|--|
| | OMB Number: 3235-0287 | | | | | | | | | |
| | Estimated average burden | | | | | | | | | |
| ı | 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

| Name and Address of Reporting Person* Mattoon Dawn | | | | | 2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX] | | | | | | | | (Check | all app | ionship of Reportir all applicable) Director Officer (give title | | rson(s) to Is 10% O Other (| wner | | |
|---|--|--|--------------|---|--|-------------------|--|---|-------------------------------------|--------------|--|---------------|----------------------------|---|--|---|-----------------------------------|--|---|--|
| (Last) (First) (Middle) C/O QUANTERIX CORPORATION, | | | | | 3. Date of Earliest Transaction (Month/Day/Year) 11/02/2021 | | | | | | | | | X | below | | iagn | below) | | |
| 900 MIDDLESEX TURNPIKE | | | | 4. If . | 4. If Amendment, Date of Original Filed (Month/Day/Year) | | | | | | | | | 6. Individual or Joint/Group Filing (Check Applicable Line) | | | | | | |
| (Street) BILLERICA MA 01821 | | | | | | | | | | | | ľ | X | | | | | | | |
| (City) | (Sta | ate) (Z | Zip) | | | | | | | | | | | | 1 0130 | , , , , , , , , , , , , , , , , , , , | | | | |
| | Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned | | | | | | | | | | | | | | | | | | | |
| 1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/Yea | | | | ear) if | 2A. Deemed Execution Date if any (Month/Day/Ye | | Code (In | | | | | | | nd 5) Secui Benet Owne | | rities Fo ficially (D ed Following Inc | | m: Direct or rect (I) | 7. Nature of Indirect Beneficial Ownership (Instr. 4) | |
| | | | | | | | | Code | v | Amoun | t | (A) or (D) | Price | | Transa | Reported Transaction(s) (Instr. 3 and 4) | | u. 4) | (IIISU. 4) | |
| Common Stock 11/02 | | | | 11/02/202 | 21 | | | | S | | 240 | [1) | D | \$52.39 | 74(2) | 36,608(3) | | | D | |
| Common | Stock | | | 11/02/202 | 1 | | | S | | 93(| 1) | D | D \$53.1449 ⁽⁴⁾ | | 36,515 | | | 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) | Exe if an | Deemed cution Date, ry nth/Day/Year) | | action (Instr. | 5. Nu of Deriv Secu Acqu (A) o Dispo of (D) (Insti | vative irities uired r osed) r. 3, 4 | Expiration Date (Month/Day/Year) | | 7. Title and Amount of Securities Underlying Derivative Security (Ins 3 and 4) | | Deri Sec (Ins | Price of rivative curity str. 5) | 9. Number derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4) | Owners Form: Direct (I or Indire (I) (Instr | Ownership | 11. Nature of Indirect Beneficial Ownership (Instr. 4) | | |
| | | | | | Code | v | (A) | (D) | Dat Exe | e ercisab | | iration | Title | or Number of Shares | r | | | | | |

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 \$51.86 to \$52.85. 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 13,966 restricted stock units.
- 4. This transaction was executed in multiple trades at prices ranging from \$52.96 to \$53.78. 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

Remarks:

/s/ Brian Keane, as Attorneyin-Fact

11/04/2021

** 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.