FORM 4

Check this box if no longer subject

to Section 16. Form 4 or Form 5 obligations may continue. See

Instruction 1(b).

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

| STATEMENT | OF | CHANGES | IN BEI | NEFICIAL | OWNERS | HIP |
|-----------|----|----------------|--------|----------|--------|-----|
| | | | | | | |

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

| 1. Name and Address of Reporting Person* MEISTER PAUL M | | | | | 2. Issuer Name and Ticker or Trading Symbol Quanterix Corp [QTRX] | | | | | | | | | | all app | | ng Pers | 10% O | wner |
|--|---|------------|-----------------------------|--|---|---|--------|---|--------------------------|--|--------------------|--|---|-----------|-----------------------|--|---------|--|--|
| · ` | | CORPORATIO | Middle) <mark>N</mark> , | | | ate of E 01/202 | | Trans | saction (Month/Day/Year) | | | | | | Office below | er (give title v) | | Other (below) | specify |
| 900 MIDDLESEX TURNPIKE | | | | 4. If Amendment, Date of Original Filed (Month/Day/Year) | | | | | | | | | 6. Individual or Joint/Group Filing (Check Applicable Line) | | | | | | |
| (Street) BILLER | ICA M | A 0 | 1821 | | | | | | | | | | | X | | | | | |
| (City) | (St | ate) (2 | Zip) | | | | | | | | | | | | | | | | |
| | | Table | I - No | on-Deriva | tive S | Secui | rities | Acc | quirec | d, Dis | sposed of | , or E | Benefic | ially | Own | ed | | | |
| 1. Title of Security (Instr. 3) 2. Transaction Date (Month/Day/ | | | | - 1 | Execution Da | | | 3. Transaction Code (Instr. 8) | | 4. Securities Acquired (A) Disposed Of (D) (Instr. 3, 4 | | | and 5) Secur Benef | | cially I Following | 6. Ownership Form: Direct (D) or Indirect (I) (Instr. 4) | | 7. Nature of Indirect Beneficial Ownership (Instr. 4) | |
| | | | | | | | | Code | v | Amount | (A) or (D) | Price | Trans | | action(s) 3 and 4) | | | (111341.4) | |
| Common Stock 10/01/2 | | | 10/01/20 |)21 | | | A | | 269(1) | A | \$51. | 1.18(2) | | 13,930(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 Execution Date Execution Date | | | ition Date, | | Transaction Code (Instr. 8) Sec Acq (A) Disport | | osed) r. 3, 4 | Expiration Da | | ate | e and 7. Title and Amount of Securities Underlying Derivative Security (Ins 3 and 4) | | | | 9. Number derivative Securities Beneficially Owned Following Reported Transaction (Instr. 4) | y G | 10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4) | 11. Nature of Indirect Beneficial Ownership (Instr. 4) |
| | | | | | Code | v | (A) | (D) | Date Exerc | isable | Expiration Date | Title | or Number of Shares | | | | | | |

Explanation of Responses:

- 1. Consists of stock granted in lieu of cash fees for service on the Company's Board of Directors and committees thereof for the third quarter of 2021.
- 2. Closing price of the Company's common stock on the Nasdaq Global Market on October 1, 2021.
- 3. Includes 2,270 restricted stock units.

Remarks:

/s/ Brian Keane, Attorney-in-

<u>fact</u>

10/05/2021 ** 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.