| SEC Form 4 |
|------------|
|------------|

FORM 4

UNITED STATES SECURITIES AND EXCHANGE COMMISSION Washington, D.C. 20549

STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

OMB APPROVAL

| OMB Number: | 3235-0287 |
|-----------------------|-----------|
| Estimated average but | rden |
| hours per response: | 0.5 |

| obligations may continue. See Instruction 1(b). | Filed pursuant to Section 16(a) of the Securities Exchange Act of 193 | hours per response: | 0.5 | | |
|--|---|-----------------------------------|------------------------------------|--------|--|
| | 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 <u>Quanterix Corp</u> [QTRX] | (Check all applicat X Director | 10% Owr | ner | |
| (Last) (First) (Middle) C/O QUANTERIX CORPORATION, | 3. Date of Earliest Transaction (Month/Day/Year) 07/01/2020 | - Officer (g below) | ive title Other (sp below) | Decify | |
| 900 MIDDLESEX TURNPIKE | 4. If Amendment, Date of Original Filed (Month/Day/Year) | 6. Individual or Joi Line) | int/Group Filing (Check Applicable | | |
| (Street) | | X Form filed | Form filed by One Reporting Perso | | |
| BILLERICA MA 01821 | | Form filed Person | d by More than One Repor | rting | |
| (City) (State) (Zip) | | | | | |

Table I - Non-Derivative Securities Acquired, Disposed of, or Beneficially Owned

| 1. Title of Security (Instr. 3) | 2. Transaction Date (Month/Day/Year) | 3. Transa Code (8) | ction | | | | Securities Beneficially | Form: Direct (D) or Indirect | 7. Nature of Indirect Beneficial Ownership (Instr. 4) |
|---------------------------------|--|------------------------------|-------|---------------------------|---------------|-----------------------|------------------------------------|---------------------------------|---|
| | | Code | v | Amount | (A) or (D) | Price | Transaction(s) (Instr. 3 and 4) | | (1150.4) |
| Common Stock | 07/01/2020 | Α | | 494 ⁽¹⁾ | Α | \$27.8 ⁽²⁾ | 10,230 ⁽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) | 2. Conversion or Exercise Price of Derivative Security | 3. Transaction Date (Month/Day/Year) | 3A. Deemed Execution Date, if any (Month/Day/Year) | 4. Transa Code (8) | | of Deriv | r osed) 7. 3, 4 | 6. Date Exercisable and Expiration Date (Month/Day/Year) | | 7. Title and Amount of Securities Underlying Derivative Security (Instr. 3 and 4) | | 8. Price of Derivative Security (Instr. 5) | 9. Number of derivative Securities Beneficially Owned Following Reported Transaction(s) (Instr. 4) | 10. Ownership Form: Direct (D) or Indirect (I) (Instr. 4) | 11. Nature of Indirect Beneficial Ownership (Instr. 4) |
| | | | | Code | v | (A) | (D) | Date Exercisable | Expiration Date | Title | Amount 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 second quarter of 2020, pursuant to an election made in December 2019. 2. Closing price of the Company's common stock on the Nasdaq Global Market on July 1, 2020.

3. Includes 2,270 restricted stock units.

Remarks:

/s/ Brian Keane, Attorney-in-

<u>fact</u>

07/06/2020

** 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.

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