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

 $\square$ 

## FORM 4

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

OMB APPROVAL

I

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

## STATEMENT OF CHANGES IN BENEFICIAL OWNERSHIP

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 burden |     |  |  |  |  |  |  |  |  |
| hours per response:      | 0.5 |  |  |  |  |  |  |  |  |

| 1. Name and Address of Reporting Person <sup>*</sup><br>Orticerio Ernest     |               |                | 2. Issuer Name <b>and</b> Ticker or Trading Symbol<br><u>Quanterix Corp</u> [ QTRX ] |                        | tionship of Reporting Pe<br>all applicable)<br>Director                                 | rson(s) to Issuer<br>10% Owner           |
|--|---------------|----------------|--|------------------------|---|--|
| (Last) (First) (Middle)<br>C/O QUANTERIX CORPORATION,<br>113 HARTWELL AVENUE |               | ( )            | 3. Date of Earliest Transaction (Month/Day/Year)<br>12/13/2017                       |                        | Officer (give title<br>below)<br>SVP Finance and C                                      | Other (specify<br>below)<br>orporate Dev |
| (Street)<br>LEXINGTON<br>(City)  | MA<br>(State) | 02421<br>(Zip) | 4. If Amendment, Date of Original Filed (Month/Day/Year)                             | 6. Indiv<br>Line)<br>X | idual or Joint/Group Filin<br>Form filed by One Rep<br>Form filed by More tha<br>Person | orting Person                            |

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

| 1. Title of Security (Instr. 3) | 2. Transaction<br>Date<br>(Month/Day/Year) | 2A. Deemed<br>Execution Date,<br>if any<br>(Month/Day/Year) |   |        | 4. Securities /<br>Disposed Of (<br>5) |       |                                    | Securities<br>Beneficially | (I) (Instr. 4) | 7. Nature<br>of Indirect<br>Beneficial<br>Ownership<br>(Instr. 4) |
|---------------------------------|--|---|---|--------|--|-------|------------------------------------|----------------------------|----------------|---|
|                                 | Code                                       |   | v | Amount | (A) or<br>(D)                          | Price | Transaction(s)<br>(Instr. 3 and 4) |                            | (1130.4)       |   |
| Common Stock                    | 12/13/2017                                 |   | М |        | 46,671                                 | Α     | \$2.28                             | 46,671                     | D              |   |

## Table II - Derivative Securities Acquired, Disposed of, or Beneficially Owned (e.g., puts, calls, warrants, options, convertible securities)

|   |   |  | 1   |                              |   |     |        |  |                    | -   |  |   |  |  |  |
|---|---|--|---|------------------------------|---|-----|--------|--|--------------------|---|--|---|--|--|--|
| 1. Title of<br>Derivative<br>Security<br>(Instr. 3) | 2.<br>Conversion<br>or Exercise<br>Price of<br>Derivative<br>Security | 3. Transaction<br>Date<br>(Month/Day/Year) | 3A. Deemed<br>Execution Date,<br>if any<br>(Month/Day/Year) | 4.<br>Transa<br>Code (<br>8) |   | of  |        | 6. Date Exercisable and<br>Expiration Date<br>(Month/Day/Year) |                    | 7. Title and Amount<br>of Securities<br>Underlying<br>Derivative Security<br>(Instr. 3 and 4) |  | 8. Price of<br>Derivative<br>Security<br>(Instr. 5) | 9. Number of<br>derivative<br>Securities<br>Beneficially<br>Owned<br>Following<br>Reported<br>Transaction(s)<br>(Instr. 4) | 10.<br>Ownership<br>Form:<br>Direct (D)<br>or Indirect<br>(I) (Instr. 4) | 11. Nature<br>of Indirect<br>Beneficial<br>Ownership<br>(Instr. 4) |
|   |   |  |   | Code                         | v | (A) | (D)    | Date<br>Exercisable  | Expiration<br>Date | Title   | Amount<br>or<br>Number<br>of<br>Shares |   |  |  |  |
| Stock<br>Option<br>(right to<br>buy)                | \$2.28  | 12/13/2017                                 |   | М                            |   |     | 46,671 | (1)  | 02/03/2022         | Common<br>Stock   | 46,671                                 | \$0   | 46,670   | D  |  |

Explanation of Responses:

1. All shares underlying this option have vested.

/s/ Brian Keane, as Attorney-in-12/18/2017

\*\* Signature of Reporting Person Date

Fact

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.