# SEC Form 4

Instruction 1(b).

# FORM 4

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

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

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

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

| 1. Name and Address of Reporting Person <sup>*</sup><br>CRANDELL KEITH           |         | g Person <sup>*</sup> | 2. Issuer Name <b>and</b> Ticker or Trading Symbol <u>Quanterix Corp</u> [ QTRX ]  |                  | 5. Relationship of Reporting Person(s) to Issuer<br>(Check all applicable)<br>X Director 10% Owner |                       |  |  |  |
|--|---------|-----------------------|--|------------------|--|-----------------------|--|--|--|
| (Last)   | (First) | (Middle)              | 3. Date of Earliest Transaction (Month/Day/Year)<br>04/03/2023   |                  | Officer (give title below)   | Other (specify below) |  |  |  |
| 8755 WEST HIGGINS ROAD, SUITE 1025   |         | AD, SUITE 1025        | 4. If Amendment, Date of Original Filed (Month/Day/Year)   | 6. Indi<br>Line) | 6. Individual or Joint/Group Filing (Check Applicat  |                       |  |  |  |
| (Street)   |         |                       |  | X                | Form filed by One Re   | eporting Person       |  |  |  |
| CHICAGO  | IL      | 60631                 |  |                  | Form filed by More th<br>Person  | an One Reporting      |  |  |  |
| (City)   | (State) | (Zip)                 | Rule 10b5-1(c) Transaction Indication  |                  |  |                       |  |  |  |
|  |         |                       | Check this box to indicate that a transaction was made pursuant to a contract, instruction or written plan that is intended to satisfy the affirmative defense conditions of Rule 10b5-1(c). See Instruction 10. |                  |  |                       |  |  |  |
| 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) | Code ( |   | 4. Securities<br>Disposed Of |               |                               | Beneficially<br>Owned Following                | Form: Direct<br>(D) or Indirect | 7. Nature<br>of Indirect<br>Beneficial<br>Ownership<br>(Instr. 4) |
|---------------------------------|--|---|--------|---|------------------------------|---------------|-------------------------------|--|---------------------------------|---|
|                                 |  |   | Code   | v | Amount                       | (A) or<br>(D) | Price                         | Reported<br>Transaction(s)<br>(Instr. 3 and 4) |                                 | (11150. 4)  |
| Common Stock                    | 04/03/2023                                 |   | Α      |   | 1,170(1)                     | A             | <b>\$11.16</b> <sup>(2)</sup> | 131,236 <sup>(3)</sup>                         | D                               |   |

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

|   |   |  | (* 5 )  **  | ,                            | , |   | ,   |  |   |       |   | ,  |  |                                       |  |
|---|---|--|---|------------------------------|---|---|-----|--|---|-------|---|--|--|---------------------------------------|--|
| 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) |   | 5. Number<br>of<br>Derivative<br>Securities<br>Acquired<br>(A) or<br>Disposed<br>of (D)<br>(Instr. 3, 4<br>and 5) |     | 6. Date Exerc<br>Expiration Da<br>(Month/Day/) | 7. Title and<br>Amount of<br>Securities<br>Underlying<br>Derivative<br>Security (Instr.<br>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) | 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              |  |  |                                       |  |

### 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 first quarter of 2023.

2. Closing price of the Company's common stock on the Nasdaq Global Market on April 3, 2023.

3. Includes 8,083 restricted stock units.

**Remarks:** 

#### /s/ Brian Keane, as Attorney-

in-Fact for Keith L

04/04/2023

\*\* Signature of Reporting Person Date

Crandell

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.