

| 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

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

|  |   |  |
|--|---|--|
| 1. Name and Address of Reporting Person*<br><u>HRUSOVSKY E KEVIN</u><br><br>(Last) (First) (Middle)<br><u>C/O QUANTERIX CORPORATION,</u><br><u>900 MIDDLESEX TURNPIKE</u><br><br>(Street)<br><u>BILLERICA MA 01821</u><br><br>(City) (State) (Zip) | 2. Issuer Name and Ticker or Trading Symbol<br><u>Quanterix Corp [ QTRX ]</u> | 5. Relationship of Reporting Person(s) to Issuer<br>(Check all applicable)<br><br><input checked="" type="checkbox"/> Director 10% Owner<br><input checked="" type="checkbox"/> Officer (give title below) Other (specify below)<br><u>Chairman, President &amp; CEO</u> |
|  | 3. Date of Earliest Transaction (Month/Day/Year)<br><u>02/08/2021</u>         |  |
|  | 4. If Amendment, Date of Original Filed (Month/Day/Year)                      | 6. Individual or Joint/Group Filing (Check Applicable Line)<br><br><input checked="" type="checkbox"/> Form filed by One Reporting Person<br><input type="checkbox"/> Form filed by More than One Reporting Person   |

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

| 1. Title of Security (Instr. 3) | 2. Transaction Date (Month/Day/Year) | 2A. Deemed Execution Date, if any (Month/Day/Year) | 3. Transaction Code (Instr. 8) |   | 4. Securities Acquired (A) or Disposed Of (D) (Instr. 3, 4 and 5) |            |                          | 5. Amount of Securities Beneficially Owned Following Reported Transaction(s) (Instr. 3 and 4) | 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                    |   |  |   |
| Common Stock                    | 02/08/2021                           |  | A                              |   | 18,397 <sup>(1)</sup>   | A          | \$0.00                   | 1,015,936 <sup>(2)</sup>  | D  |   |
| Common Stock                    | 02/09/2021                           |  | s <sup>(3)</sup>               |   | 700   | D          | \$76.8077 <sup>(4)</sup> | 1,015,236 <sup>(2)</sup>  | D  |   |
| Common Stock                    | 02/09/2021                           |  | s <sup>(3)</sup>               |   | 3,100   | D          | \$78.2426 <sup>(5)</sup> | 1,012,136 <sup>(2)</sup>  | D  |   |
| Common Stock                    | 02/09/2021                           |  | s <sup>(3)</sup>               |   | 1,200   | D          | \$78.8079 <sup>(6)</sup> | 1,010,936 <sup>(2)</sup>  | 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. Transaction Code (Instr. 8) |   | 5. Number of Derivative Securities Acquired (A) or Disposed of (D) (Instr. 3, 4 and 5) |     | 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 |
| Stock Option (right to buy)                | \$77.79  | 02/08/2021                           |  | A                              |   | 42,925   |     | (7)  | 02/08/2031      | Common Stock  | 42,925                                     | \$0.00   | 42,925  | D  |       |

**Explanation of Responses:**

- Represents shares of stock issuable on vesting of restricted stock units. The restricted stock units vest as to 25% of the shares on February 8, 2022, and the remaining 75% vest in 36 equal monthly installments at the end of each month beginning on February 28, 2022.
- Includes 146,668 restricted stock units and 39,815 shares of common stock that are subject to performance based vesting.
- The sale of shares was effected pursuant to a Rule 10b5-1 trading plan adopted on December 14, 2020.
- This transaction was executed in multiple trades at prices ranging from \$76.53 to \$77.30. 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.
- This transaction was executed in multiple trades at prices ranging from \$77.57 to \$78.55. 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.
- This transaction was executed in multiple trades at prices ranging from \$78.64 to \$79.16. 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.
- The option vests as to 25% of the shares on February 8, 2022, and the remaining 75% vests in 36 equal monthly installments at the end of each month beginning on February 28, 2022.

**Remarks:**

/s/ Brian Keane, as Attorney-in-Fact 02/10/2021

\*\* 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.