

**UNITED STATES  
SECURITIES AND EXCHANGE COMMISSION**

Washington, D.C. 20549

**FORM 8-K**

**CURRENT REPORT  
Pursuant to Section 13 or 15(d)  
of the Securities Exchange Act of 1934**

Date of Report (Date of earliest event reported): **August 29, 2018**

**QUANTERIX CORPORATION**

(Exact name of registrant as specified in its charter)

**Delaware**  
(State or other jurisdiction  
of incorporation)

**001-38319**  
(Commission File Number)

**20-8957988**  
(IRS Employer  
Identification No.)

**113 Hartwell Avenue**  
**Lexington, MA**  
(Address of principal executive offices)

**02421**  
(zip code)

Registrant's telephone number, including area code: **(617) 301-9400**

Check the appropriate box below if the Form 8-K filing is intended to simultaneously satisfy the filing obligation of the registrant under any of the following provisions:

- Written communications pursuant to Rule 425 under the Securities Act (17 CFR 230.425)
- Soliciting material pursuant to Rule 14a-12 under the Exchange Act (17 CFR 240.14a-12)
- Pre-commencement communications pursuant to Rule 14d-2(b) under the Exchange Act (17 CFR 240.14d-2(b))
- Pre-commencement communications pursuant to Rule 13e-4(c) under the Exchange Act (17 CFR 240.13e-4(c))

Indicate by check mark whether the registrant is an emerging growth company as defined in Rule 405 of the Securities Act of 1933 (17 CFR §230.405) or Rule 12b-2 of the Securities Exchange Act of 1934 (17 CFR §240.12b-2).

Emerging Growth Company

If an emerging growth company, indicate by check mark if the registrant has elected not to use the extended transition period for complying with any new or revised financial accounting standards provided pursuant to Section 13(a) of the Exchange Act.

---

**Item 1.01 Entry into a Material Definitive Agreement**

On August 30, 2018, Quanterix Corporation (the "Company") entered into an amendment to the Loan and Security Agreement, dated April 14, 2014, as amended, with Hercules Capital, Inc. (the "Loan Agreement"). The amendment extends the interest only payment period through March 1, 2020 and also extends the loan maturity date to March 1, 2020. As of June 30, 2018, there was approximately \$7.7 million outstanding under the Loan Agreement.

**Item 1.02 Termination of a Material Definitive Agreement.**

On August 29, 2018, the Company exercised an option to terminate the Lease Agreement, dated September 24, 2007, by and between RAR2 Boston Industrial QRS-MA, Inc. and Aushon Biosystems, Inc. (a wholly owned subsidiary of the Company), as amended (the "Lease"), for approximately 21,500 rentable square feet of space at 43 Manning Road, Billerica, Massachusetts. The termination of the Lease will be effective as of September 1, 2019. In accordance with the terms of the Lease, the Company is required to pay \$75,933 no later than July 1, 2019 in consideration for the early termination.

---

**SIGNATURES**

Pursuant to the requirements of the Securities Exchange Act of 1934, the Registrant has duly caused this report to be signed on its behalf by the undersigned thereunto duly authorized.

**QUANTERIX CORPORATION**

By: /s/ Joseph Driscoll  
Joseph Driscoll  
Chief Financial Officer

Date: September 5, 2018