



Quanterix Announces Closing of Public Offering Including Exercise of Underwriters' Option to Purchase Additional Shares

August 13, 2019

BILLERICA, Mass.--(BUSINESS WIRE)--Aug. 13, 2019-- Quanterix Corporation (Nasdaq: QTRX), a company digitizing biomarker analysis with the goal of advancing the science of precision health, today announced the closing of its previously announced underwritten public offering of 2,732,673 shares of its common stock at a public offering price of \$25.25 per share, including 356,435 shares sold pursuant to the full exercise of the underwriters' option to purchase additional shares. Gross proceeds from the sale of the shares, before deducting underwriting discounts and commissions and offering expenses, were approximately \$69.0 million.

J.P. Morgan Securities LLC and SVB Leerink LLC acted as joint book-running managers for the offering. Canaccord Genuity LLC acted as co-manager for the offering. Perella Weinberg Partners acted as independent capital markets advisor to Quanterix for the offering.

The public offering was made pursuant to a shelf registration statement on Form S-3 that was previously filed with and declared effective by the Securities and Exchange Commission ("SEC"). The final prospectus supplement and the accompanying prospectus relating to this offering has been filed with the SEC and is available on the SEC's website located at www.sec.gov, copies of which can be obtained from J.P. Morgan Securities LLC, c/o Broadridge Financial Solutions, 1155 Long Island Avenue, Edgewood, NY 11717, telephone: (866) 803-9204; or SVB Leerink LLC, Attention: Syndicate Department, One Federal Street, 37th Floor, Boston, MA, 02110, by telephone at (800) 808-7525, ext. 6132 or by e-mail at syndicate@svbleerink.com.

This press release shall not constitute an offer to sell, or a solicitation of an offer to buy, nor will there be any sale of these securities in any state or other jurisdiction in which such an offer, solicitation or sale would be unlawful prior to the registration or qualification under the securities laws of any such state or other jurisdiction.

About Quanterix

Quanterix is a company that's digitizing biomarker analysis with the goal of advancing the science of precision health. The company's digital health solution, Simoa, has the potential to change the way in which healthcare is provided today by giving researchers the ability to closely examine the continuum from health to disease. Quanterix' technology is designed to enable much earlier disease detection, better prognoses and enhanced treatment methods to improve the quality of life and longevity of the population for generations to come. The technology is currently being used for research applications in several therapeutic areas, including oncology, neurology, cardiology, inflammation and infectious disease. The company was established in 2007 and is located in Billerica, Massachusetts.

View source version on businesswire.com: <https://www.businesswire.com/news/home/20190813005721/en/>

Source: Quanterix Corporation

Quanterix IR Contact:

Amol Chaubal, CFO, 617-301-9495
achaubal@quanterix.com

Quanterix PR Contact:

PAN Communications
Erin Smith, 617-502-4371
quanterix@pancomm.com