



Quanterix to Participate in the Canaccord Genuity 42nd Annual Growth Conference

August 3, 2022

BILLERICA, Mass.--(BUSINESS WIRE)--Aug. 3, 2022-- [Quanterix Corporation](#) (NASDAQ: QTRX), a company expanding the limits of exploration with ultrasensitive biomarker detection, today announced that President and Chief Executive Officer Masoud Toloue will present at the [Canaccord Genuity 42nd Annual Growth Conference](#) on Wednesday, August 10, 2022 at 1:30 p.m. ET in Boston, MA. Toloue will also host one-on-one and group meetings with institutional investors that day.

A [live webcast](#) of the conversation will be available on the Investors section of the Quanterix website at <https://ir.quanterix.com/>. Replays of the webcast will be available on the Quanterix website for 90 days following the conference.

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. For additional information, please visit <https://www.quanterix.com>.

Forward-Looking Statements

This press release contains forward-looking statements within the meaning of the Private Securities Litigation Reform Act of 1995. Words such as "may," "will," "expect," "plan," "anticipate," "estimate," "intend" and similar expressions (as well as other words or expressions referencing future events, conditions or circumstances) are intended to identify forward-looking statements. Forward-looking statements in this news release are based on Quanterix' expectations and assumptions as of the date of this press release. Each of these forward-looking statements involves risks and uncertainties. Factors that may cause Quanterix' actual results to differ from those expressed or implied in the forward-looking statements in this press release are discussed in Quanterix' filings with the U.S. Securities and Exchange Commission, including the "Risk Factors" sections contained therein. Except as required by law, Quanterix assumes no obligation to update any forward-looking statements contained herein to reflect any change in expectations, even as new information becomes available.

View source version on [businesswire.com](https://www.businesswire.com/news/home/20220803005859/en/): <https://www.businesswire.com/news/home/20220803005859/en/>

Media:

PAN Communications
Lauren Force
(610) 574-0565
pan.quanterix@pancomm.com

Investor Relations:

Stephen Hrusovsky
(774) 278-0496
ir@quanterix.com

Source: Quanterix Corporation