



Discovery Fueled by Ultra-Sensitivity

Quanterix Presenting at Canaccord Genuity 44th Annual Growth Conference

August 6, 2024

BILLERICA, Mass.--(BUSINESS WIRE)--Aug. 6, 2024-- [Quanterix Corporation](#) (NASDAQ: QTRX), a company fueling scientific discovery through ultra-sensitive biomarker detection, today announced that Chief Executive Officer Masoud Toloue and Chief Financial Officer Vandana Sriram will speak in a fireside chat at the [Canaccord Genuity 44th Annual Growth Conference](#) on August 13, 2024, at 3:00 p.m., ET.

The presentation will be available virtually and participants can register using [this link](#). The presentation will be available following the conference on Quanterix's website [here](#).

To learn more about Quanterix, visit www.quanterix.com/about. To learn more about Quanterix's Simoa® technology, visit: <https://www.quanterix.com/simoa-technology/>.

About Quanterix

From discovery to diagnostics, Quanterix's ultra-sensitive biomarker detection is fueling breakthroughs only made possible through its unparalleled sensitivity and flexibility. The Company's Simoa® technology has delivered the gold standard for earlier biomarker detection in blood, serum or plasma, with the ability to quantify proteins that are far lower than the Level of Quantification (LoQ). Its industry-leading precision instruments, digital immunoassay technology and CLIA-certified Accelerator laboratory have supported research that advances disease understanding and management in neurology, oncology, immunology, cardiology and infectious disease. Quanterix has been a trusted partner of the scientific community for nearly two decades, powering research published in more than 2,900 peer-reviewed journals. Find additional information about the Billerica, Massachusetts-based company at <https://www.quanterix.com> or follow us on [Twitter](#) and [LinkedIn](#).

View source version on [businesswire.com](https://www.businesswire.com/news/home/20240806815882/en/): <https://www.businesswire.com/news/home/20240806815882/en/>

Media:

PAN Communications
Maya Nimnicht
510-334-6273
quanterix@pancomm.com

Investor Relations:

Quanterix
Francis Pruell
(508)-789-1725
ir@quanterix.com

Source: Quanterix Corporation