



Discovery Fueled by Ultra-Sensitivity

## Quanterix to Participate at the Morgan Stanley 21st Annual Global Healthcare Conference

Sep 5, 2023

BILLERICA, Mass.--(BUSINESS WIRE)--Sep. 5, 2023-- [Quanterix Corporation](#) (NASDAQ: QTRX), a company fueling scientific discovery through ultrasensitive biomarker detection, today announced that Chief Executive Officer Masoud Toloue will take part in a fireside chat at the Morgan Stanley 21<sup>st</sup> Annual Global Healthcare Conference on September 13, 2023 at 3:35 p.m. ET. The presentation will be available virtually and participants can register [here](#). The presentation will be available for viewing following the conference on Quanterix's website here: [ir.quenterix.com](#).

To learn more about Quanterix, visit [www.quenterix.com/company/](#) To learn more about Quanterix's Simoa® technology, visit: [www.quenterix.com/simoa-technology/](#).

### About Quanterix

From discovery to diagnostics, Quanterix's ultrasensitive 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 Limit of Quantification (LoQ) of conventional analog methods. 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 2,000 peer-reviewed journals. Find additional information about the Billerica, Massachusetts-based company at <https://www.quenterix.com> or follow us on [Twitter](#) and [LinkedIn](#).

View source version on [businesswire.com](#): <https://www.businesswire.com/news/home/20230905628089/en/>

### Media Contact:

PAN Communications  
Maya Nimnicht  
510-334-6273  
[quenterix@pancomm.com](mailto:quenterix@pancomm.com)

### Investor Relations Contact:

Ed Joyce  
610-306-9917  
[ir@quenterix.com](mailto:ir@quenterix.com)

Source: Quanterix Corporation