



Discovery Fueled by Ultra-Sensitivity

Quanterix to Present at 45th Annual Goldman Sachs Global Healthcare Conference

June 5, 2024

BILLERICA, Mass.--(BUSINESS WIRE)--Jun. 5, 2024-- [Quanterix Corporation](#) (NASDAQ: QTRX), a company fueling scientific discovery through ultrasensitive biomarker detection, today announced that Chief Executive Officer Masoud Toloue will speak at the Goldman Sachs 45th Annual Global Healthcare Conference. Toloue will speak alongside Chief Financial Officer Vandana Sriram in a fireside chat presentation on Wednesday, June 12, 2024, at 2:40 p.m. EST. The session will be made available to attendees and the general public via webcast.

Webcast Information

To access the live webcast of Quanterix's fireside chat presentation at the conference, please visit [this link](#).

Replays of the presentation will be available for a limited period following the conference. The presentation will also be made available through the [Investor Relations](#) section of the company's website.

To learn more about Quanterix, visit www.quanterix.com/company. To learn more about Quanterix's Simoa® technology, visit www.quanterix.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 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): <https://www.businesswire.com/news/home/20240605682576/en/>

Media Contact:

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

Investor Relations Contact:

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

Source: Quanterix Corporation