



Discovery Fueled by Ultra-Sensitivity

Quanterix to Present at Leerink Partners Healthcare Crossroads Conference 2024

May 22, 2024

BILLERICA, Mass.--(BUSINESS WIRE)--May 22, 2024-- [Quanterix Corporation](#) (NASDAQ: QTRX), a company fueling scientific discovery through ultrasensitive biomarker detection, today announced that Chief Executive Officer Masoud Toloue will speak in a fireside chat at the 2024 Leerink Partners Healthcare Crossroads Conference. Quanterix's presentation is scheduled for Wednesday, May 29, 2024, at 11:40 a.m. ET. The fireside chat will be made available to attendees and to the public via webcast.

Webcast Information

The live webcast of Quanterix's presentation can be accessed [here](#), and from the Investor section of the company's website at www.Quanterix.com. Replays of the presentation will be available for a limited time following the conference.

To learn more about Quanterix, visit www.quanterix.com/about. 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/20240522008584/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