



Discovery Fueled by Ultra-Sensitivity

Quanterix To Report First Quarter 2024 Financial Results

April 29, 2024

BILLERICA, Mass.--(BUSINESS WIRE)--Apr. 29, 2024-- [Quanterix Corporation](#) (NASDAQ: QTRX), a company fueling scientific discovery through ultrasensitive biomarker detection, today announced that it will host a conference call on Wednesday, May 8, 2024 at 8:30 a.m. E.T., to discuss its first quarter 2024 financial results.

Quanterix will issue a press release regarding the first quarter 2024 earnings prior to the conference call on Tuesday, May 7, 2024 after the market closes. The press release will be posted on the Quanterix website at <https://www.quanterix.com/>.

To pre-register for the conference call with a webcast link [click here](#) and for an audio only link [click here](#) to obtain your dial-in number and passcode. Interested investors can also access the live webcast from the [News & Events](#) page within the Investors section of the Quanterix website at <http://www.quanterix.com>. An archived webcast replay will be available on the Company's website for one year.

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 2,900 peer-reviewed journals. Find additional information about the Billerica, Massachusetts-based company at <https://www.quanterix.com> or follow us on Twitter and LinkedIn.

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/20240429829499/en/): <https://www.businesswire.com/news/home/20240429829499/en/>

Media:

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

Investor Relations:

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

Source: Quanterix Corporation