

Quanterix™

Quanterix Corporation CEO to Present at MassBio: Possible Talk Series

June 26, 2020

BILLERICA, Mass.--(BUSINESS WIRE)--Jun. 26, 2020-- [Quanterix Corporation](#) (NASDAQ: QTRX), a company digitizing biomarker analysis to advance the science of precision health, today announced that Kevin Hrusovsky, Chairman and Chief Executive Officer of Quanterix and Founder of the non-profit [Powering Precision Health](#) (PPH), will participate in the [MassBio: Possible Talk](#) series on June 26, 2020 at 10 a.m. EDT. In this Possible Talk, "Building A Scientific/Political Ecosystem to Get America Safely Back On Its Feet," Hrusovsky explores the intersection of these three elements against the backdrop of COVID-19. He'll detail how the PPH syndicate is harnessing cutting-edge biomarker technology to fuel critical advancements in early detection, enable deadly cytokine storms prediction, and deliver empirical insights into immunity that could help overcome testing challenges and set the stage for reliable, widespread homecare testing.

A recording will be available at www.quanterix.com.

About Quanterix

Quanterix is a company that's digitizing biomarker analysis with the goal of advancing the science of precision health. The company's digital health solution, Simoa, has the potential to change the way in which healthcare is provided today by giving researchers the ability to closely examine the continuum from health to disease. Quanterix' technology is designed to enable much earlier disease detection, better prognoses and enhanced treatment methods to improve the quality of life and longevity of the population for generations to come. The technology is currently being used for research applications in several therapeutic areas, including oncology, neurology, cardiology, inflammation and infectious disease. The company was established in 2007 and is located in Billerica, Massachusetts. For additional information, please visit <https://www.quanterix.com>.

About Powering Precision Health

Powering Precision Health is the world's first independent, non-profit organization dedicated to bringing the world's leading physicians, scientists, innovators, investors and patient advocates together to unveil their latest research on new biomarkers that are revolutionizing precision health. Founded by Kevin Hrusovsky, a widely acclaimed thought leader and visionary in life sciences and personalized medicine, Powering Precision Health is a movement that represents the intersection of new technological capabilities with the latest medical research. It's rooted in the science of precision medicine, which shows personalized treatments to be an increasingly more effective way to maximize drug efficacy and minimize toxicity. In addition to the impact environmental and lifestyle factors can have on minimizing disease triggers, precision health marks an evolution in the way we approach disease and aims to inspire a full healthcare transformation, from philosophy to approach to outcome. In an industry often plagued by skepticism and marred by false promises, PPH puts science first and brings together stakeholders that span from fundamental research to clinical practice, investors, policy makers, patient advocacy groups, and anyone who embraces the vision of Powering Precision Health. Featuring a distinguished keynote lineup of dignitaries, the Summit unveils groundbreaking approaches to prevention, early diagnosis, and next-generation treatments. Powering Precision Health is supported thanks to the generous contributions of sponsors from a wide range of companies and organizations committed to advancing precision health.

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/20200626005267/en/): <https://www.businesswire.com/news/home/20200626005267/en/>

Media:

PAN Communications
Staci Didner, (407) 734-7325
quanterix@pancomm.com

Investor Relations:

Stephen Hrusovsky
(774) 278-0496
shrusovsky@quanterix.com

Source: Quanterix Corporation