

Quanterix

Quanterix Named One of the World's Most Innovative Companies in 2016 by Fast Company *Innovative Technology Detects Concussions with Single Finger Prick*

Lexington, Mass. – March 1, 2016 — [Quanterix Corporation](#), a leader in the transformation of healthcare through its ultrasensitive single molecule testing capabilities, today announced that it has been selected by *Fast Company* as one of the world's top 10 most innovative companies in 2016 in [Sports](#) – the only medical technology company to be recognized in this category. *Fast Company's* annual ranking of the world's [Most Innovative Companies](#) honors leading enterprises and rising newcomers that exemplify the best in nimble business and impactful innovation.

“We are honored to be recognized by *Fast Company* for our ground-breaking single molecule detection technology and its ability to detect even the mildest of concussions through a simple blood test,” said Kevin Hrusovsky, CEO and Executive Chairman, Quanterix. “Our technology is truly revolutionary and we’re excited for what’s to come in 2016 as we continue to grow and apply the science to other therapeutic areas, such as oncology, to redefine remission and transform today’s sick care to true, preventative health care.”

With proven applications across a wide spectrum of therapeutic areas, including oncology, cardiology, neurology and infectious disease, Quanterix’ ultra-sensitive diagnostic platform, Simoa, is capable of measuring individual proteins at concentrations 1000 times lower levels than the best immunoassays available today. For example, this sensitivity enables the direct measurement of a cancer biomarker in a single cell. The Company’s annual growth has exceeded 100 percent quarter over quarter in the past 18 months, and is expected to continue that momentum by further accelerating the vast potential of their technology in 2016.

Quanterix has received significant industry recognition in the past year alone, including several prominent industry award wins, such as the [GE – NFL Head Health Challenge](#), and the technology has been featured in over 50 distinguished scientific publications, including *Nature*, *JAMA Neurology* and *Clinical Chemistry*. Hrusovsky recently delivered a dynamic presentation at the JP Morgan Healthcare Conference outlining how Quanterix is powering a revolution in healthcare: <http://info.quanterix.com/watch-quanterix-jp-morgan-healthcare-presentation>.

“We’re excited by Quanterix’ technology and the potential it has to transform public health,” said Jeff Miller, NFL senior vice president of health and safety policy. “We look forward to continuing to support their work and that of all Head Health Challenge winners that will soon benefit athletes, the military and society in general.”

The “Most Innovative Companies” list is one of *Fast Company's* most significant and highly anticipated editorial efforts of the year. To produce the 2016 list, the magazine's editors spent a full year gathering and analyzing data on companies around the globe. For the first time, *Fast Company* enlisted machine intelligence to help inform the list: Quid, a San Francisco-based startup, used its software to create topographical maps of innovation across more than 40 sectors of the economy to visually organize the year’s most relevant and prominent trends.

About Quanterix

Quanterix is a developer of ground-breaking tools in high definition diagnostics. Its Simoa platform uses single molecule measurements to access previously undetectable proteins. With this unprecedented

sensitivity and full automation, Simoa offers significant benefits to both research and clinical testing applications. Quanterix was established in 2007 and is located in Lexington, Massachusetts. To learn more about Quanterix and Simoa, please visit: www.quanterix.com.

###

Contacts:

Caitlyn Keating

PAN Communications

617-502-4300

quanterix@pancomm.com