



One Kendall Square, Building 1400
Cambridge, MA 02139

617.301.9400 877.786.8749

Quanterix Develops Microfluidic Consumable That Will Enable Next Generation Molecular Diagnostic Systems Based on Single Molecule Array Technology

Lab on a Chip Publication Highlights Performance of Integrated Arrays

CAMBRIDGE, MA – December 22, 2011 – Quanterix Corporation, a company enabling a new generation of molecular diagnostic tests based on its revolutionary Single Molecule Array (SiMoA™) technology, has reported a method for detecting individual proteins within microfabricated polymer arrays produced by Sony DADC. Quanterix recently entered into a collaboration with Sony DADC for the development and manufacture of “smart consumables” based on an optical disc format. This method was published online in *Lab on a Chip*.

“We have developed a device that enables the fluidic sealing and isolation of single molecules in arrays containing hundreds of thousands of miniaturized wells. These consumables are manufactured using high precision and high volume manufacturing processes, and will be at the heart of the SiMoA instrument that we are currently developing. By using arrays enclosed in microfluidic channels we are able to perform SiMoA in a fully automated fashion.” said David Duffy, Ph.D., Vice President of Research at Quanterix and corresponding author of the publication. “This represents a significant milestone towards our goal of developing an automated platform that enables single molecule sensitivity while offering significant advantages for the manufacture of high quality, low-cost consumables. These devices will facilitate high throughput and sensitive detection of biomarkers, and address important applications in life science research and in vitro diagnostics.”

Martin Madaus, Ph.D., Quanterix Chairman and CEO added, “This achievement is an important step towards our goal to commercialize the SiMoA technology. Working closely with our strategic partners, Quanterix plans to develop and manufacture a fully automated instrument for its SiMoA technology that will utilize these consumables to support a menu of diagnostic tests. Quanterix expects to launch the life science instrument in 2013, followed by an IVD platform to be commercialized in 2014.”

About Quanterix

Quanterix Corporation is developing its proprietary Single Molecule Array (SiMoA™) technology for the in vitro diagnostics and life science research markets. The digital nature of SiMoA yields unprecedented assay performance, stemming from a 1,000-fold improvement in sensitivity compared with today’s analog only technology. SiMoA will enable researchers in life science to validate novel, low abundance biomolecules from a single droplet of blood, leading to greater insight into disease detection, diagnosis, therapy selection and disease monitoring. Automated systems based on SiMoA will provide diagnostic test information to healthcare practitioners faster, with greater reliability, unprecedented range and increased cost effectiveness. Founded in 2007, the privately held Cambridge, Massachusetts-based company is backed by leading life science investors including ARCH Venture Partners, Bain Capital Ventures, and Flagship Ventures. For additional information, please visit www.quanterix.com.

Media Contact:

David Hanlon

Quanterix Corporation

Tel: 617-301-9400

Email: media@quanterix.com