

Quanterix

Quanterix Selected as a Red Herring Top 100 North America Winner

Lexington, MA – May 28, 2014 – Quanterix, a leader in high definition diagnostics, delivering ultrasensitive single molecule measurement for the benefit of human health, announced today it has been selected as a Red Herring Top 100 North America award winner. The Top 100 award winners are selected in recognition of the leading private companies from the Americas, celebrating these startups’ innovations and technologies across their respective industries.

Red Herring Top 100 America enlists outstanding entrepreneurs and promising companies. It selects 100 award winners from among the approximately 3,000 tech startups financed each year in the U.S. and Canada. Since 1996, Red Herring has kept tabs on these up-and-comers. Red Herring editors were among the first to recognize that companies such as Facebook, Twitter, Google, Yahoo, Skype, Salesforce.com, YouTube, Marin Software, Palo Alto Networks and eBay would change the way we live and work.

“We are honored to be recognized by Red Herring and proud to be included in this prestigious group of premier tech companies in North America,” said Paul Chapman, President and CEO at Quanterix. “The fact that Quanterix was one of only a few Life Sciences companies recognized this year further illustrates the importance and potential benefit that our Simoa technology will have on human health”

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.