

# Quanterix

## Quanterix one of 33 Recipients of Massachusetts Life Sciences Center Tax Incentive Award

**Lexington, Mass. – February 20, 2014** — [Quanterix Corporation](#), a leader in high definition diagnostics, delivering ultrasensitive single molecule measurement for the benefit of human health, today announced that it was one of the 33 recipients of the [Massachusetts Life Sciences Center tax](#) awards. The Life Sciences Tax Incentive Program authorizes up to \$25 million in tax incentives each year for companies engaged in life science research and development, commercialization and manufacturing, and is designed to incentivize Massachusetts life science companies to create new jobs.

“We are honored to be a repeat winner of the Mass Life Sciences Center awards,” said Paul Chapman, President and CEO of Quanterix. “This award validates the exciting work Quanterix is doing with our ground-breaking ultrasensitive Simoa technology, and we are very appreciative of the MLSC support for the opportunity it provides us to further accelerate our strategy execution. It is particularly timely considering we are now in full commercial launch phase and are expanding our operations accordingly.”

For more information on the Massachusetts Life Sciences Center awards, please visit: <http://www.masslifesciences.com/>.

### 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/simoa-is-here](http://www.quanterix.com/simoa-is-here).

###

### Contacts:

Nikki Festa & Caitlyn Keating  
PAN Communications  
617-502-4300  
[quanterix@pancomm.com](mailto:quanterix@pancomm.com)