

# APPLIED CLINICAL TRIALS

December 16, 2014

(UVPM: 22,328)

## **Quanterix Announces the Simoa Accelerator**

Quanterix Corporation, a provider of high definition diagnostics, delivering ultrasensitive single molecule measurement, announced the opening of The Simoa Accelerator, a dedicated laboratory environment for custom assay development and clinical sample testing. Designed to provide access to the revolutionary Single Molecule Array (Simoa) HD-1 Analyzer, this state-of-the-art facility allows life science researchers to explore biomarkers and pathways with greater sensitivity and precision, paving the way for the development of new diagnostic tests and improved treatment of diseases in oncology, neurology, cardiology and other fields of medicine.