

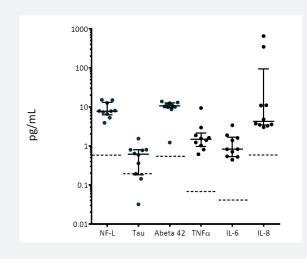
ULTRA-SENSITIVE BIOMARKER DETECTION

For the benchtop





- Multiplexed detection without compromising sensitivity or specificity
- Simple 2-step or 3-step workflow with minimal user intervention
- Compact instrument with built-in touch-screen control and comprehensive data analysis tools
- · No daily or monthly user maintenance
- Broad menu of commercially available kits
- Simple development of custom homebrew Simoa assays



Multiplex measurement of low abundance circulating proteins

Quantitative detection of normal endogenous neurodegeneration biomarkers and inflammatory cytokine levels using the Simoa bead-based assay kits on the SR-X. Distribution of sample readings are shown along with LoQ of each assay (dashed lines).



Sample to answer for 96 tests in as little as 4 hours with minimal hands-on time!

Intuitive User-guided Run Setup and Integrated Data Analysis



Initiate experiment from home screen



Define plate lavout



Analyze results and generate report

Visit quanterix.com/SR-X for more information

The SR-X Ultra-Sensitive Biomarker Detection System is the latest instrument from Quanterix powered by Simoa technology, offering researchers access to ultra-sensitive protein detection capabilities in a compact and affordable system. The SR-X is designed to support multiplexed detection of up to four biomarkers per sample, with low volume requirements to increase throughput and productivity while conserving precious samples.

A menu of over 25 Simoa assay kits have been validated on the SR-X with an additional 60 Simoa bead-based assay kits are available to researchers to measure critical biomarkers with 1000X higher sensitivity than standard immunoassay methods, enabling detection of both normal and acute biomarker levels with high precision across a range of sample types.



Quanterix Corporation 900 Middlsex Turnpike, Building 1 Billerica, MA 01821

