

simOa<sup>®</sup>



SR-X<sup>™</sup>

Ultrasensitive Biomarker Detection System

Quanterix<sup>®</sup>

# Ultrasensitive Biomarker Detection

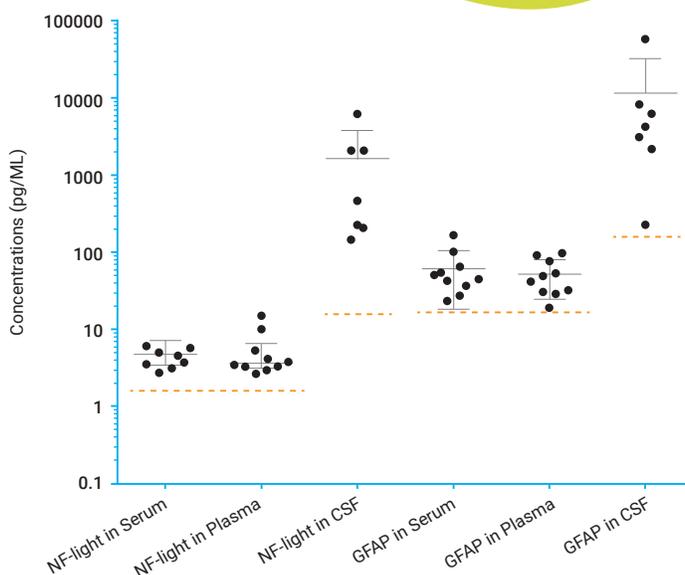
benchtop instrument and workflows supporting flexible lab settings

Simoa SR-X™ empowers scientists to discover, detect and treat, faster and with deeper sensitivity than ever before.



## SR-X™ unique capabilities

- Proprietary Simoa® Discs allow single bead isolation into femto-sized wells and single molecule detection
- Simple 2-step or 3-step workflow with minimal user intervention
- Compact instrument with built-in touch-screen control and comprehensive data analysis tools
- Broad menu of commercially available kits
- Simple development of custom home brew Simoa® assays
- Multiplexed detection without compromising sensitivity or specificity



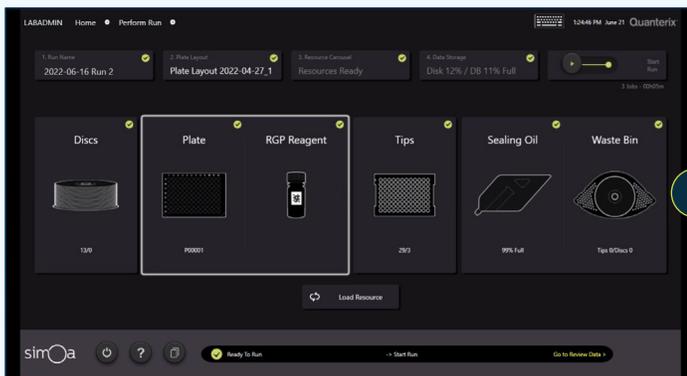
## Multiplex measurement of low abundance circulating proteins

Multiplex measurement of neurodegeneration biomarkers using the Simoa® bead-based assay on the SR-X™. Distribution of healthy donor sample readings are shown along with the LLOQ (dashed lines) of each assay in each validated matrix.

## The value of The SR-X System

The Quanterix SR-X<sup>®</sup> Ultrasensitive Biomarker Detection System is the latest benchtop instrument from Quanterix powered by Simoa<sup>®</sup> technology, offering researchers access to ultra-sensitive protein detection capabilities in a compact and affordable system. The SR-X<sup>™</sup> 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.

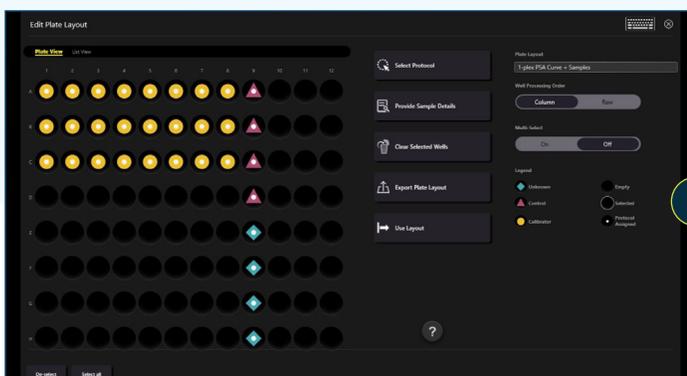
A menu of transferable kits between all bead-based assay instruments are available to researchers to measure critical biomarkers with higher sensitivity than standard immunoassay methods, enabling the detection of both normal and acute biomarker levels with high precision across a range of sample types.



1

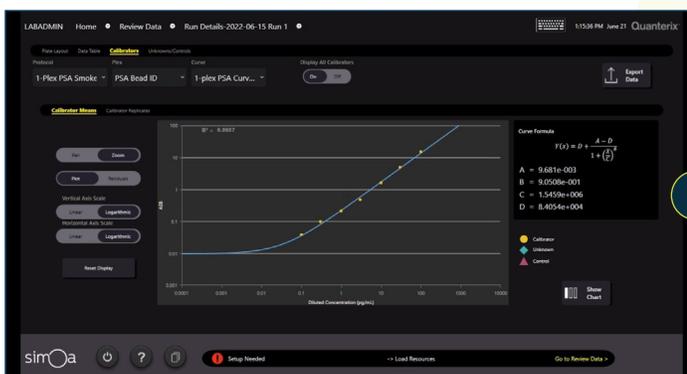
### Intuitive user-guided run setup

1. Initiate experiment from home screen
2. Define plate layout
3. Analyze results and generate report



2

Speed to insight in as little as 4 hours with minimal user intervention.



3

# Quanterix SR-X<sup>®</sup> workflow: simple, scalable, flexible.



1

Add reagents to assay plate



2

Simoa Microplate Shaker

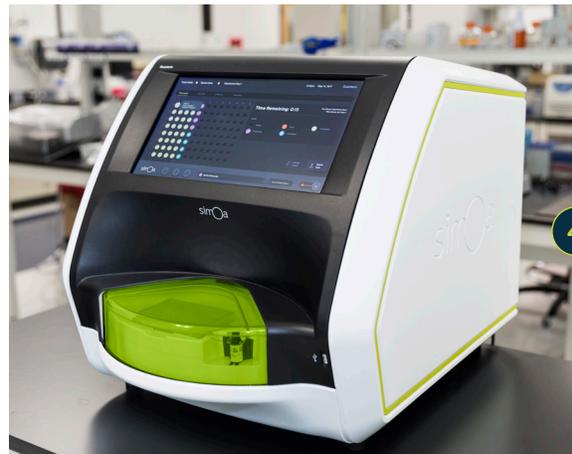
Re-suspend Simoa magnetic beads with automated shaking and incubation



3

Simoa Microplate Washer

Wash beads with automated magnetic separation



4

Load assay plate onto SR-X for single molecule detection on Simoa discs

Visit [quanterix.com/SR-X](https://www.quanterix.com/SR-X) for more information



[www.quanterix.com](https://www.quanterix.com) | 900 Middlesex Turnpike, Building One | Billerica, MA 01821 USA

©2023 Quanterix Corporation. Simoa<sup>®</sup> is a registered trademark of Quanterix Corporation for research use only. Not for use in diagnostic procedures.