

Which Instrument Is Right for You?

Quanterix offers three instruments for clinical and diagnostic ultra-sensitive biomarker research. Publications utilizing Quanterix instruments have spanned diverse fields such as neurology, oncology, cardiology, immunology, infectious disease, and beyond. Each instrument enables ultra-sensitive protein quantitation, surpassing traditional automated immunoassay analyzers by up to 1000 times. In addition, these instruments offer flexibility in assay options, supporting [Simoa® immunoassay kits](#) as well as [Homebrew Custom Assays](#).



HD-X ANALYZER™



SR-X™ BIOMARKER DETECTION SYSTEM



SP-X IMAGING AND ANALYSIS SYSTEM™

Simoa® Technology	Bead	Bead	Planar
Multiplex capacity	Up to 4	Up to 4	Up to 10
Automation	Fully automated	Semi-automated	Semi-automated
Throughput	3 x 96-well plates / 8 hr	1 x 96 well plate / 3-4 hr	1 x 96-well plate / 3-4 hr
Sample Prep	Integrated automated onboard	Offboard	Offboard
Footprint	Floor-standing	Benchtop	Benchtop
Dimensions	141.4 x 89.7 x 161.2 cm (55.66 x 35.3 x 63.5 in)	55.9 x 72.9 x 55.2 cm (22.0 x 28.7 x 21.7 in)	28 x 61 x 36.5 cm (11 x 24 x 14.4 in)



For more information, contact your Quanterix representative.



Scan the QR code to be directed to an archive of publications.