

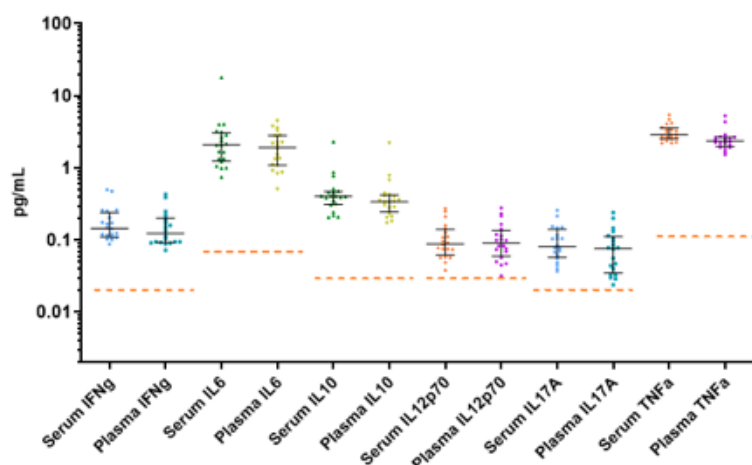
Simoa® Cytokine 6-Plex

Measures IFN γ , IL-6, IL-10, IL-12p70,
IL-17A, TNF α

The Simoa Cytokine 6-Plex Panel 1 is the most sensitive multiplex cytokine assay available to quantitatively measure 6 key immunomodulatory biomarkers in serum and plasma at normal baseline levels.

- Available for use on the HD-X and SR-X platforms
- Measures all six analytes in serum and EDTA plasma using only 50 μ l sample
- Pre-mixed ready-to-use reagents for simple assay setup and processing workflow

Simoa Human Cytokine Panel 1 Samples Reading (HD-X)



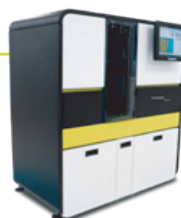
Ultra-sensitive multiplexing
without compromise

Healthy donor matched EDTA plasma (n=20), and serum (n=20) were measured. Bars depict median with interquartile range.

Comprehensive assay validation report and data sheets available demonstrating analytical performance exceeding any competitive multiplex assay capable of measuring up to 6 analytes per test.

The Quanterix SR-X™:

The first benchtop instrument to offer true multiplex detection at both acute and baseline levels.



The Simoa HD-X Analyzer™:

A fully automated instrument capable of running up to 6-plex assays.

Order your 6-Plex Today! | sales@quanterix.com

For research use only. Not for use in diagnostic procedures.

Product Code 102958