

Ultra-Sensitive, Fully Automated Simoa® Cytokine 4-Plex Advantage PLUS Assays

Quanterix®
Discovery Fueled by Ultra-Sensitivity

Three panels, 11 essential cytokine biomarkers, femtogram sensitivity

With Simoa® ultra-sensitive protein quantification, the cutting-edge Advantage PLUS assay platform, and full automation of the HD-X Analyzer, the Simoa® Cytokine 4-Plex Advantage PLUS Assays will transform cytokine biomarker analysis, delivering deeper insights and scalable results like never before.



Ultra-sensitivity

Uncover critical insights into early biological and pathological processes from low-abundance cytokines at femtogram levels.



Multiplex

Quantify four cytokines in a single assay with 25 µL input, conserving samples and resources.



Full Automation

Enhance efficiency and reproducibility with walk-away automation on the HD-X.



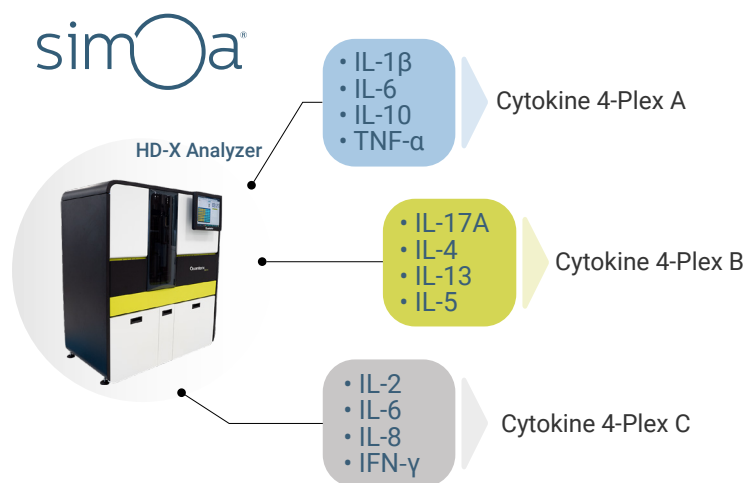
Scalability

Adapt the automated workflow for fast time-to-results or high-throughput, generating data from one 96-well plate in under 3 hours or up to 6 plates daily.



Lot-to-Lot Consistency

Improve reliability for longitudinal and large-scale studies.



Unlock crucial insights early

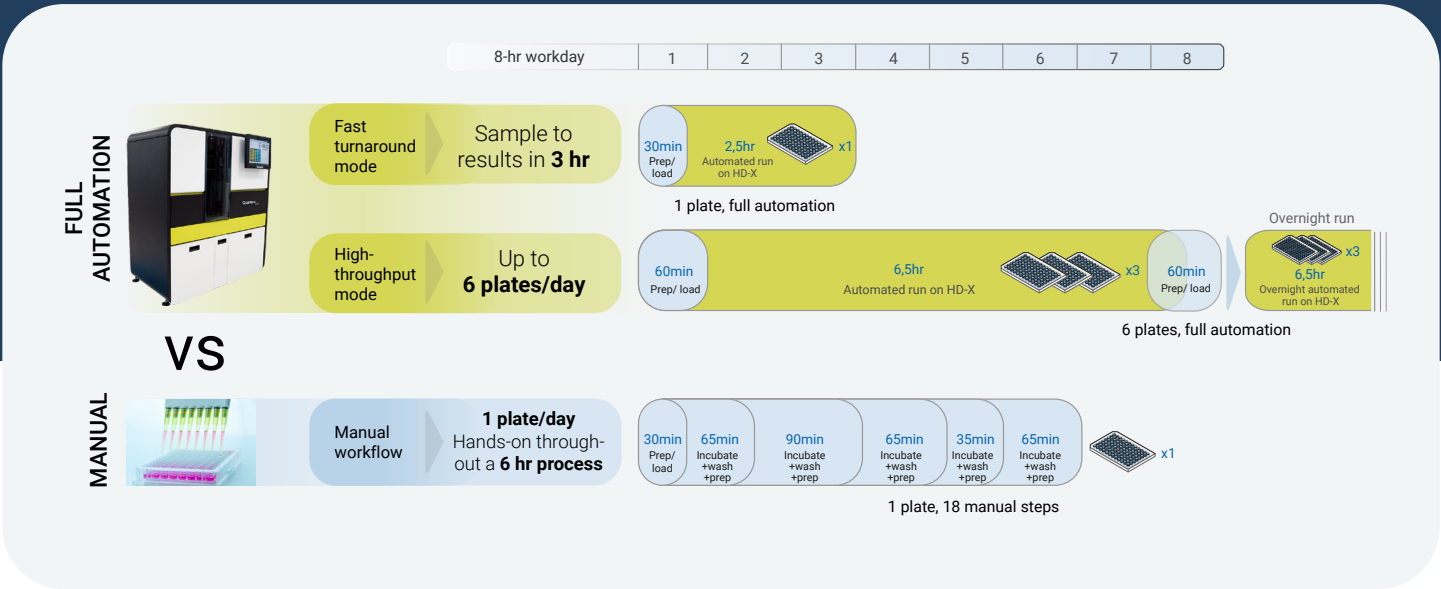
The Simoa® Cytokine 4-Plex Advantage PLUS Assays deliver sensitivity down to fg/mL levels, reliably measuring low-abundance cytokines at baseline in healthy individuals and capturing subtle yet critical signals early in biological and pathological processes.

Experience robust consistency

Powered by the Simoa® Advantage PLUS assay technology, the Simoa® Cytokine 4-Plex Advantage PLUS assays are designed to deliver consistent quality and performance lot after lot - a key to the success of large-scale and longitudinal studies. Full automation on the HD-X further enhances reproducibility by minimizing variables associated with manual handling.

Accomplish a week's work in a single day

With the Simoa® Cytokine 4-Plex Advantage PLUS Assays fully automated on the HD-X Analyzer, you can accomplish a week's worth of work in just an 8-hour day - without being tied to the bench, freeing up your research staff for other important tasks. Set up in 30 minutes, walk away, and return to results in as little as 3 hours. Need higher throughput? Process up to 6 plates in a single day - a productivity that would take a full week with a manual workflow.



Performance at-a-glance

The Simoa® Cytokine 4-plex Advantage PLUS Assays detect the target cytokine analytes in serum and EDTA samples down to fg/mL levels, enabling a high rate of detection of endogenous proteins across target analytes in samples spanning the health-disease spectrum. **Visit the [Quanterix website](#) for detailed assay specifications.**

	C4PA Advantage PLUS				C4PB Advantage PLUS				C4PC Advantage PLUS			
	IL-1β	IL-6	IL-10	TNF-α	IL-17A	IL-4	IL-13	IL-5	IL-2	IL-6	IL-8	IFN-γ
LOD (pg/mL)	0.002	0.008	0.033	0.061	0.002	0.003	0.004	0.004	0.005	0.008	0.014	0.012
LLOQ (pg/mL)	0.01	0.028	0.143	0.481	0.008	0.01	0.034	0.01	0.019	0.031	0.07	0.075
Dynamic range* (pg/mL)	0 - 48	0 - 160	0 - 300	0 - 400	0 - 48	0 - 40	0 - 60	0 - 60	0 - 80	0 - 160	0 - 320	0 - 160

LOD: lower limit of detection; LLOQ: lower limit of quantification; * Serum/EDTA plasma at minimal required dilution of 4x

Supercharge your cytokine biomarker analysis with the power of the Simoa® Cytokine Solution.

Simoa® Cytokine 4-Plex Advantage PLUS Assays		
Full Product Name	Target Cytokines	Product Number
Cytokine 4-plex A (C4PA) Advantage PLUS	IL-1β, IL-6, IL-10, TNF-α	104979
Cytokine 4-plex B (C4PB) Advantage PLUS	IL-17A, IL-4, IL-13, IL-5	104997
Cytokine 4-plex C (C4PC) Advantage PLUS	IL-2, IL-6, IL-8, IFN-γ	105066

Quanterix Corporation
900 Middlesex Turnpike, Billerica, MA 01821

©2024 Quanterix Corporation. Simoa® and Quanterix® are registered trademarks of Quanterix Corporation. For research use only. Not for use in diagnostic procedures.

Quanterix®
Discovery Fueled by Ultra-Sensitivity
www.quanterix.com