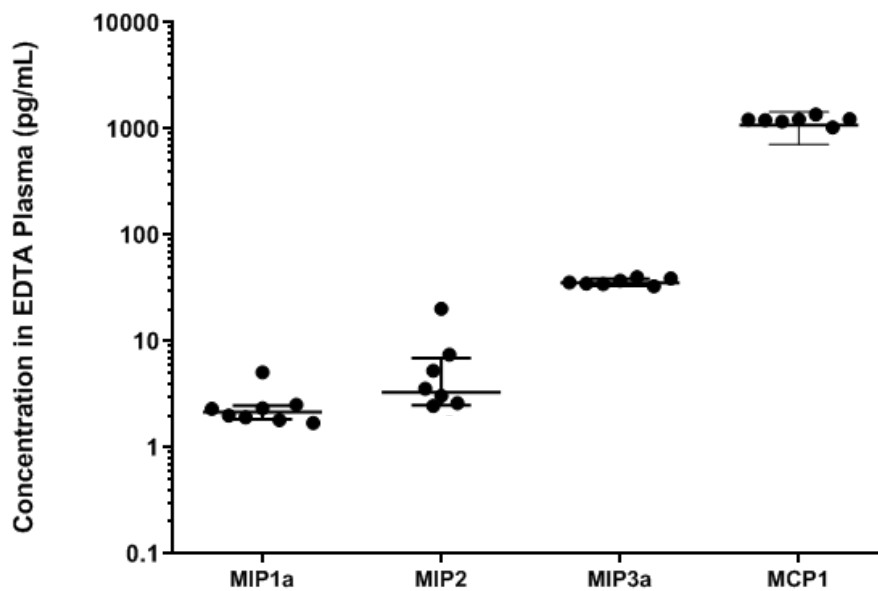


Simoa Planar Array Rat Chemokine Panel 1

Catalog Number 88-0016

MIP-1α, MIP-2, MIP-3α, MCP-1

The Simoa Planar Array Rat Chemokine Panel 1 assay is a multiplex immunoassay designed for the Quanterix SP-X™ Imaging and Analysis system. The assay measures MIP-1α, MIP-2, MIP-3α and MCP-1 in rat serum and plasma samples. Representative normal healthy sample readings are listed in the graph below along with the limit of detection (LoD) and limits of quantitation (LLoQ and ULoQ) for each biomarker in the assay. Simoa assay kits are available for any combination of the biomarkers in the panel from 1-plex to 4-plex.



Rat Chemokine 4-Plex Panel 1				
	MIP-1α	MIP-2	MIP-3α	MCP-1
LoD (pg/ml)	0.088	0.23	1.8	0.21
LLoQ (pg/ml)	0.29	0.39	3.12	1.17
ULoQ (pg/ml)	300	400	3,200	1,200