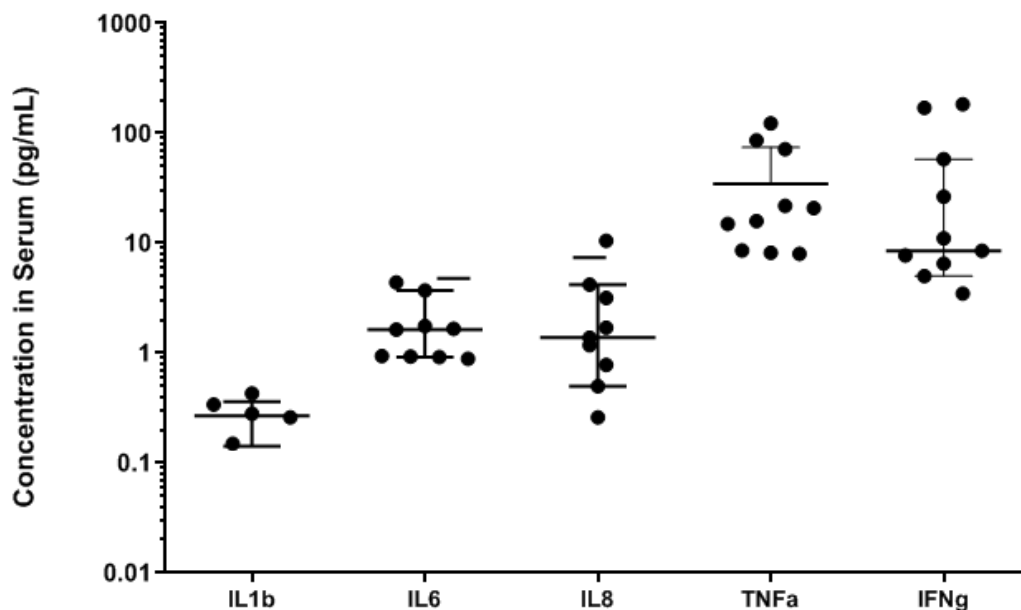


# Simoa Planar Array Non-human Primate Cytokine Panel 1

Catalog Number 85-0013

## IL-1β, IL-6, IL-8, TNF-α, IFN-γ

The Simoa Planar Array NHP Cytokine Panel 1 assay is a multiplex immunoassay designed for the Quanterix SP-X™ Imaging and Analysis system. The assay measures IL-1β, IL-6, IL-8, TNF-α and IFN-γ in non-human primate 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 5-plex.



NHP Cytokine 5-Plex Panel 1					
	IL-1β	IL-6	IL-8	TNF-α	IFN-γ
LoD (pg/ml)	0.09	0.005	0.012	0.47	0.36
LLOQ (pg/ml)	0.12	0.11	0.78	1.95	0.97
ULoQ (pg/ml)	125	114	400	2,000	1,000