

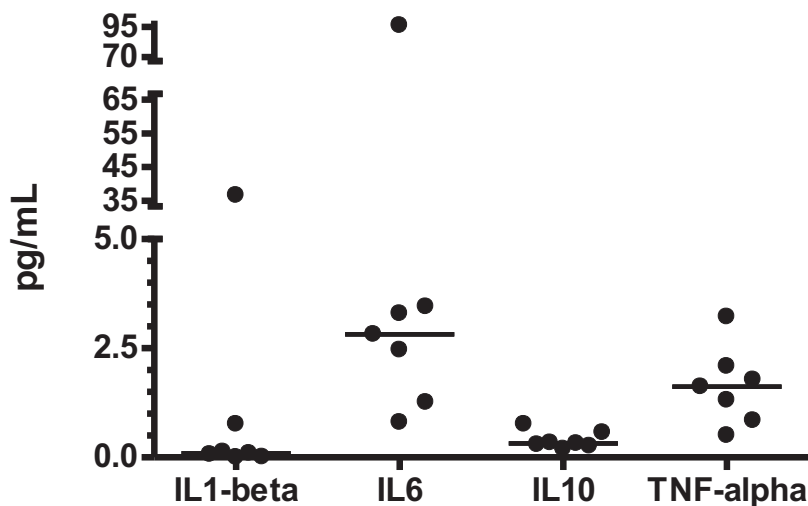


# Ultra-Sensitive Biomarker Quantification Assay Services

powered by Simoa™ technology

- Single Molecule Array (Simoa™) technology enables **femtogram/ml** level measurement of biomarkers in **serum, plasma**, or other matrices
- Selected **multiplex** panels available
- Provides robust and reproducible results with **< 10% CVs** and the capacity for high precision 'digital' or 'traditional' analog measurement
- Ultra-sensitive services covering over 25 analytes including A $\beta$ -40, A $\beta$ -42, GM-CSF, IL-1 $\beta$ , IL-6, IL-8, IL-17A, IL-23, TNF- $\alpha$ , TRAIL and Troponin-I

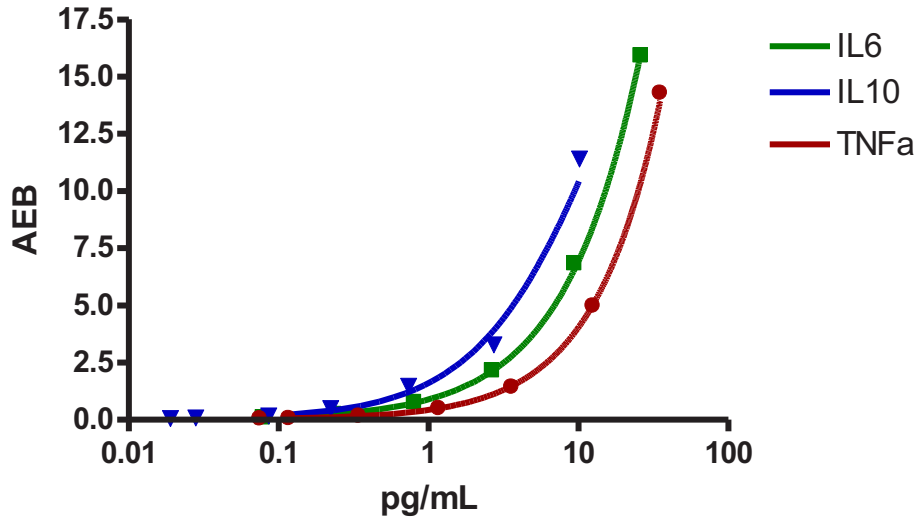
**Figure 1.** Normal human serum endogenous levels of several pro- and anti-inflammatory cytokines



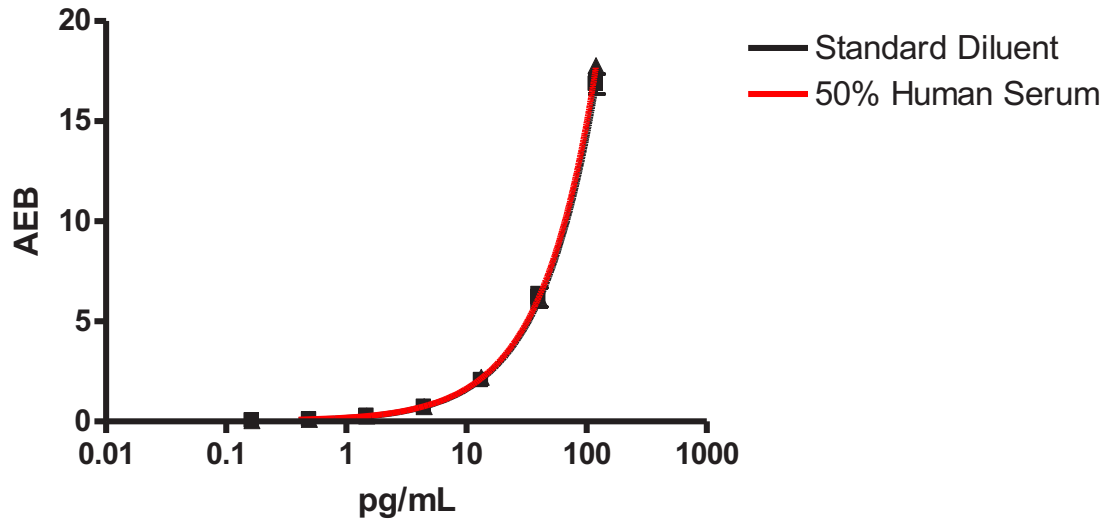
Powered by the Quanterix Simoa™ Technology, Lexington MA, USA



**Figure 2.** Human Cytokine 3-Plex A (IL-10, IL-6, TNF- $\alpha$ ) Immunoassay Standard Curves  
 AEB: Average Enzyme per Bead

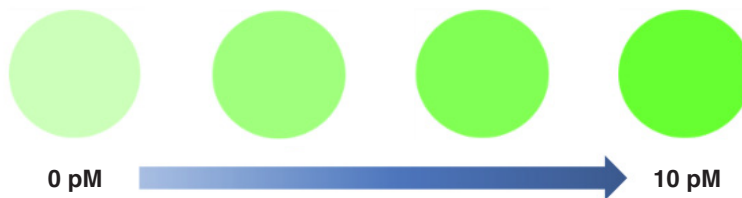


**Figure 3.** IL-1beta Immunoassay Standard Curve



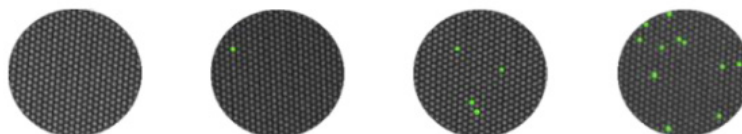
**Figure 4.** Traditional (analog) detection versus high precision Simoa (digital) measurement

**Traditional (analog)**



Analog measurements give increasing intensity as the concentration increases.

**Simoa (digital)**



In contrast, digital measurements are independent of intensity and simply rely on a signal/no signal readout.