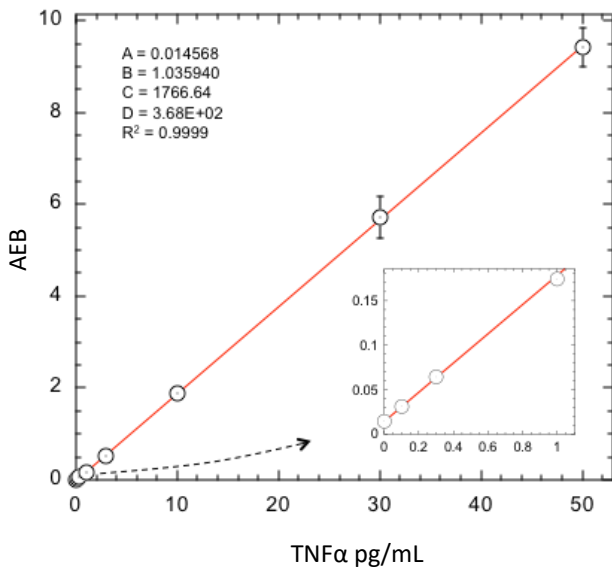


**Description**

Human tumor necrosis factor alpha (TNFα) is a homotrimeric transmembrane protein that functions as a proinflammatory cytokine. It is produced mainly by macrophages but also by a variety of other cell types, including monocytes, neutrophils, and T-cells. The involvement of TNFα in several signal transduction pathways links the protein to such diverse functions as acute inflammation, apoptosis, septic shock, cellular proliferation, and differentiation. Human TNFα is a non-glycosylated protein of 157 amino acids, with a molecular weight of approximately 17,000 daltons. The clinical relevance of TNFα stems from its association with numerous disease states including rheumatoid arthritis, cancer, cachexia, and Crohn’s disease.

**Calibration Curve:** Four-parameter curve fit parameters are depicted.



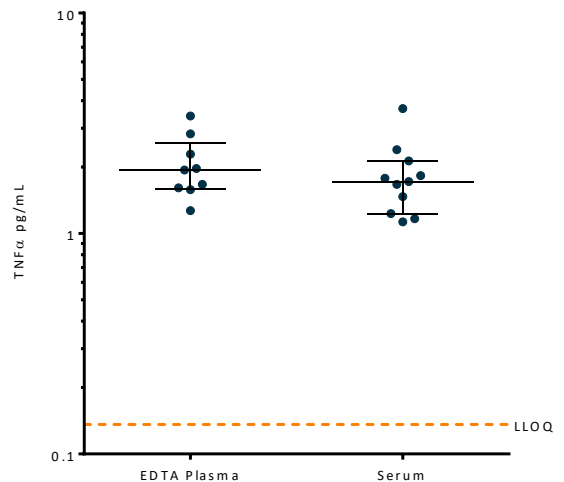
**Lower Limit of Quantification (LLOQ):** Triplicate measurements of serially diluted calibrator were read back on the calibration curve over 2 reagent lots across 3 instruments (12 runs total).

**Limit of Detection (LOD):** Calculated as 2.5 standard deviations from the mean of background signal read back on each calibration curve over 2 reagent lots across 3 instruments (12 runs total).

<b>LLOQ</b>	<b>0.034 pg/mL</b>
<b>LOD</b>	<b>0.016 pg/mL</b> SD 0.0186 pg/mL
<b>Dynamic range (serum and plasma)</b>	0–200 pg/mL
<b>Diluted Sample volume*</b>	100 μL per measurement
<b>Tests per kit</b>	192

\*See Kit Instruction for details

**Endogenous Sample Reading:** Healthy donor matched EDTA plasma (n=9) and serum (n=11) were measured. Error bars depict mean and SEM.



Sample Type	Median TNFα pg/mL	% Above LOD
Serum	1.78	100%
EDTA Plasma	1.94	100%

**Precision:** Five samples consisting of three serum-based panels and two TNFα controls were assayed in replicates of three at two separate times per day for five days using a single lot of reagents and calibrators. Analysis of variance (fully nested ANOVA) results are summarized in the following table.

Sample	Mean (pg/mL)	Within run CV	Between run CV	Between Day
Control 1	2.77	3.4%	4.1%	0.0%
Control 2	29.9	2.8%	4.8%	0.0%
Panel 1	1.71	4.9%	9.1%	0.0%
Panel 2	2.15	5.5%	9.2%	0.0%
Panel 3	43.4	3.5%	0.4%	1.0%

**Spike and Recovery:** TNFα spiked into 4 serum samples at 2 levels.

**Admixture Linearity:** High TNFα serum sample admixed with low TNFα sample, mean of 10 levels.

**Dilution Linearity:** 1 spiked serum sample was diluted 2x serially from MRD (4x) to 64x with Sample Diluent.

**Lot Consistency:** 5 samples tested with 2 reagent lots across 2 runs x 3 instruments.

<b>Spike and Recovery (Serum)</b>	<b>Mean = 77.9%</b> Range: 68.2–87.1%
<b>Admixture Linearity</b>	<b>Mean = 108%</b>
<b>Dilution Linearity (64x)</b>	<b>Mean = 121%</b> Range: 101–139%
<b>Lot Consistency</b>	<b>Mean Difference = -2.7%</b> Range: 1.34–41.1 pg/mL

The Simoa TNFα Discovery assay kit is formulated for use on the SR-X®, HD-1, or HD-X® platform. Data in this document was obtained from runs on the HD-1 platform unless otherwise noted. Some differences in performance claims between SR-X and HD-1/HD-X may be observed when comparing datasheets for these platforms. This may be due to experiments run at different time-points with different reagent lots and different samples, or it may be due to minor differences in antibody and analyte behavior in the different assay formats.