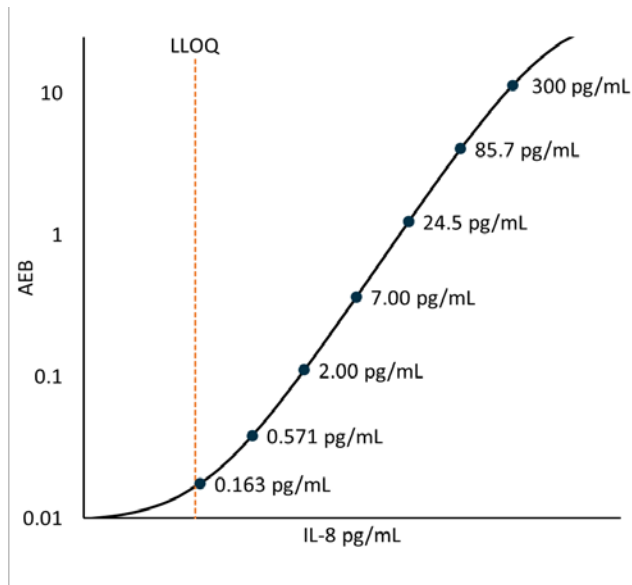


**Description**

Interleukin 8 (IL-8) is a cytokine of 72 amino acids (molecular weight 8 kDa) whose primary role is induction of chemotaxis in neutrophils, basophils, and T-cells, causing them to migrate to the site of infection. IL-8 also induces phagocytosis by the target cells. IL-8 is secreted by cells involved in the immune response to antigens, typically starting with macrophages, which release IL-8 to recruit other cells. Secretion of IL-8 is increased by oxidative stress and results in the recruitment of inflammatory cells. This recruitment further induces oxidative stress mediators, making it a key player in localized inflammation. IL-8 elevation has been associated with a range of clinical conditions, including psoriasis, chronic hepatitis C, and thyroid disease. IL-8 has recently been identified as a potential therapeutic target in inflammatory diseases.

**Calibration Curve:** Calibrator concentrations and Lower Limit of Quantification depicted.



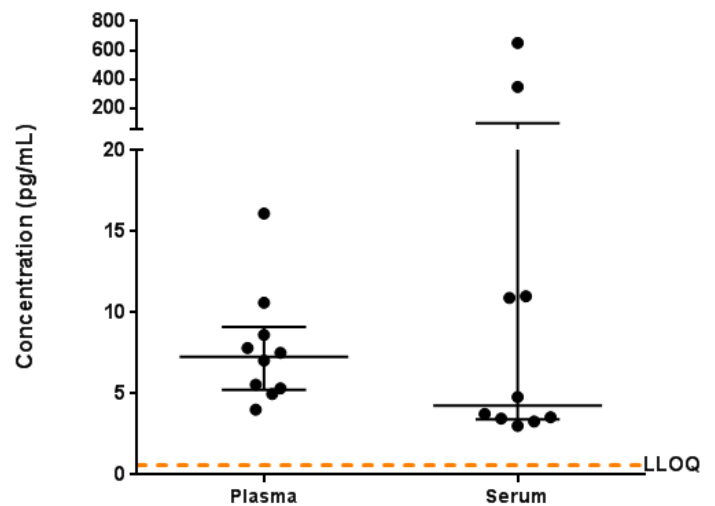
**Lower Limit of Quantification (LLOQ):** Triplicate measurements of serially diluted calibrator were read back on the calibration curve over 6 runs each for 1 reagent lot across 2 instruments (6 runs total).

**Limit of Detection (LOD):** Calculated as 2.5 standard deviations from the mean of background signal read back on each calibration curve over 6 runs each for 1 reagent lot across 2 instruments (6 runs total).

<b>LLOQ</b>	<b>0.146 pg/mL</b> pooled CV 18% mean recovery 101%
<b>LOD</b>	<b>0.0387 pg/mL</b> range 0.0111-0.0568 pg/mL
<b>Dynamic range (serum and plasma)</b>	0-300 pg/mL
<b>Diluted Sample volume*</b>	100 µL per measurement
<b>Tests per kit</b>	96

\*See Kit Instruction for details

**Endogenous Sample Reading:** Healthy donor matched EDTA plasma (n=10), and serum (n=10) were measured. Bars depict median with interquartile range. Orange line represents functional LLOQ.



**Note:** Specificity of highest measured IL-8 concentration was verified with knockdown study.

Sample Type	Mean IL-8 pg/mL	Median IL-8pg/mL	% Above LOD
Serum	104	4.27	100%
Plasma	7.76	7.28	100%

**Precision:** Measurements of 3 serum-based panels and 2 calibrator-based controls. Triplicate measurements were made for 6 runs each for 1 reagent lot across 2 instruments (6 runs total, 18 measurements).

Sample	Mean (pg/mL)	Within run CV	Between run CV	Between inst CV
Control 1	10.8	2.8%	2.3%	0.7%
Control 2	190	4.5%	4.7%	2.2%
Panel 1	5.59	3.0%	3.6%	14%
Panel 2	54.9	2.5%	4.2%	2.1%
Panel 3	335	3.4%	2.4%	5.0%

**Spike and Recovery:** 2 serum and 2 EDTA plasma samples were spiked at high and low concentrations within the range of the assay and analyzed on SR-X.

**Dilution Linearity:** 1 endogenous EDTA plasma and 1 endogenous serum sample were diluted 2X serially from MRD (4x) to 64x with Sample Diluent.

<b>Spike and Recovery (Serum/Plasma)</b>	<b>Mean = 86%</b> Range: 73-98%
<b>Dilution Linearity (64x)</b>	<b>Mean = 107%</b> Range: 99 –119%

The Simoa IL-8 assay kit is formulated for use on either the SR-X or HD-1 platform. Data in this document was obtained from runs on the SR-X platform unless otherwise noted. Some differences in performance claims between the HD-1 and SR-X may be observed when comparing datasheets for the two platforms. This may be due to experiments run at different time-points with different reagent lots and different samples, or may be due to minor differences in antibody and analyte behavior in the different assay formats.