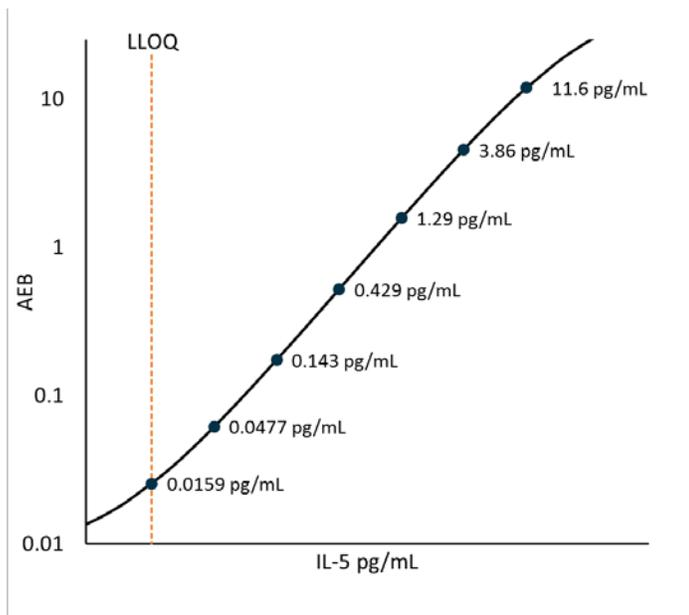


**Description**

Interleukin 5 (IL-5) is a cytokine with a length of 115 amino acids and a molecular weight of 15.2 kDa that is derived from T-cells with hematopoietic functions predominantly associated with antigen-induced eosinophilia. IL-5 induces differentiation of B-cells to immunoglobulin secreting cells and is an important factor in growth, differentiation and activation of eosinophils. IL-5, GM-CSF and IL-3 comprise the  $\beta$ -common ( $\beta$ c) cytokine family, so named because the receptors share a common  $\beta$  chain complexed with cytokine-specific  $\alpha$  chains. IL-5 and IL-5R are the targets of therapeutic antibodies for treatment of eosinophilic asthma and are involved in type 2 inflammation in the mucosal allergic reaction to grass pollen. Activation of the IL-3/IL-5/GM-CSF receptors results in rapid activation of the JAK/STAT pathway.

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



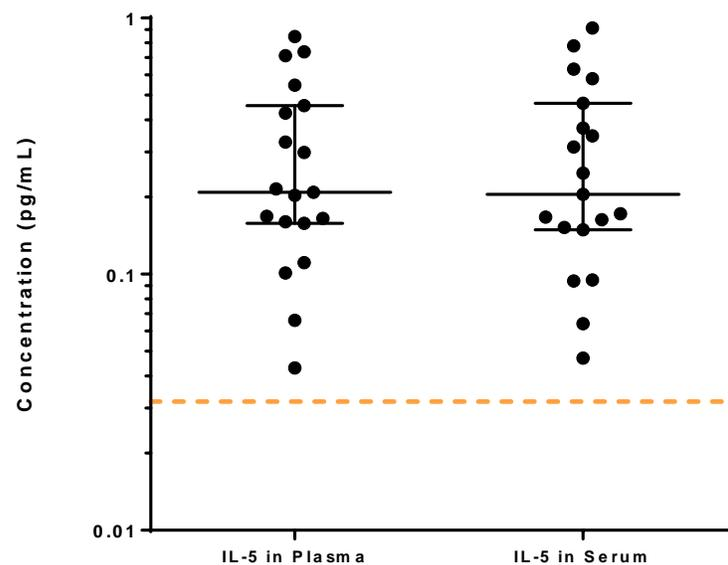
**Lower Limit of Quantification (LLOQ):** Triplicate measurements of serially diluted calibrator were read back on the calibration curve over 13 runs total over 2 reagent lots across 4 instruments.

**Limit of Detection (LOD):** Calculated as 2.5 standard deviations from the mean of background signal read back on each calibration curve over 13 runs total over 2 reagent lots across 4 instruments.

<b>LLOQ</b>	<b>0.0165 pg/mL</b> pooled CV 18% mean recovery 102%
<b>LOD</b>	<b>0.0034 pg/mL</b> range 0.0007-0.0075 pg/mL
<b>Dynamic range (serum and plasma)</b>	0 - ~12 pg/mL
<b>Diluted Sample volume*</b>	152 $\mu$ L per measurement
<b>Tests per kit</b>	96

\*See Kit Instruction for details

**Endogenous Sample Reading:** Healthy donor matched EDTA plasma (n=19), and serum (n=19) were measured. Bars depict median with interquartile range. Orange line represents functional LLOQ.



Sample Type	Mean IL-5 pg/mL	Median IL-5 pg/mL	% Above LOD
Serum	0.31	0.21	100%
Plasma	0.31	0.21	100%

**Precision:** Measurements of 3 serum-based panels and 2 calibrator-based controls. Triplicate measurements were made for 15 runs total over 2 reagent lots across 4 instruments.

Sample	Mean (pg/mL)	Within run CV	Between run CV	Between inst CV	Between Lot CV
Control 1	0.31	8.1%	16.2%	9.9%	17.9%
Control 2	5.68	14.2%	11.7%	8.3%	10.2%
Panel 1	0.53	7.6%	9.6%	8.6%	0.3%
Panel 2	1.47	6.4%	11.9%	7.6%	7.0%
Panel 3	4.97	8.5%	12.3%	8.2%	3.1%

**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:** 2 EDTA plasma (1 endogenous and 1 spiked) and 2 serum samples (1 endogenous and 1 spiked) were diluted 2X serially starting from MRD (2x).

<b>Spike and Recovery (Serum/Plasma)</b>	<b>Mean = 82%</b> Range: 64-92%
<b>Endogenous Dilution Linearity (16x)</b>	<b>Mean = 96%</b> Range: 86-110%
<b>Spiked Dilution Linearity (256x)</b>	<b>Mean = 104%</b> Range: 85-130%

The Simoa IL-5 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.