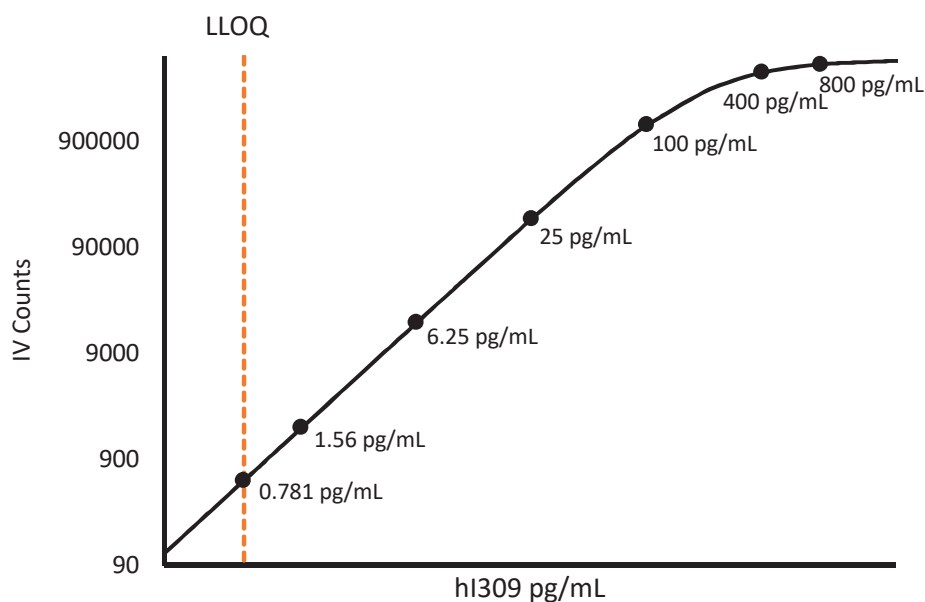


Description – I-309 (CCL-1)

I-309, also known as CCL-1, is a small 15 kDa glycoprotein secreted by activated T cells that belong to a family of inflammatory cytokines known as chemokines. CCL-1 cDNA encodes a 96 amino acid residue precursor protein with a hydrophobic signal peptide that is cleaved to yield a 73 amino acid residue mature protein. I-309 functions to attract monocytes, NK cells, immature B cells, and dendritic cells by interacting with the CCR8 receptor.

Calibration Curve: Calibrator concentrations and Lower Limit of Quantification are depicted in the figure below. This standard curve is for demonstration purposes; end users should prepare a standard curve for each assay run.



Minimum Required Dilution (MRD)

Diluted Sample volume (1:2 Dilution)*	50 µL per measurement
--	--------------------------

*See Kit Instructions for details

Endogenous Serum and Plasma Readings: Healthy EDTA plasma and serum samples (n=8) were measured.

% Above LOD	100%
% Above LLOQ	76%

Assay Range: The upper end of the dynamic range is equal to the top calibrator concentration multiplied by MRD.

Analytical LLOQ	0.781 pg/mL
Functional LLOQ (x MRD)	1.562 pg/mL
LOD	0.644 pg/mL
Assay Range	0 – 1600 pg/mL

Note: Data described were developed during assay development. Under different assay conditions, assay may perform differently than shown. For complex matrices such as serum or plasma, assay diluent optimization (for example by adding blocking agents) may improve performance of these matrices in this assay.