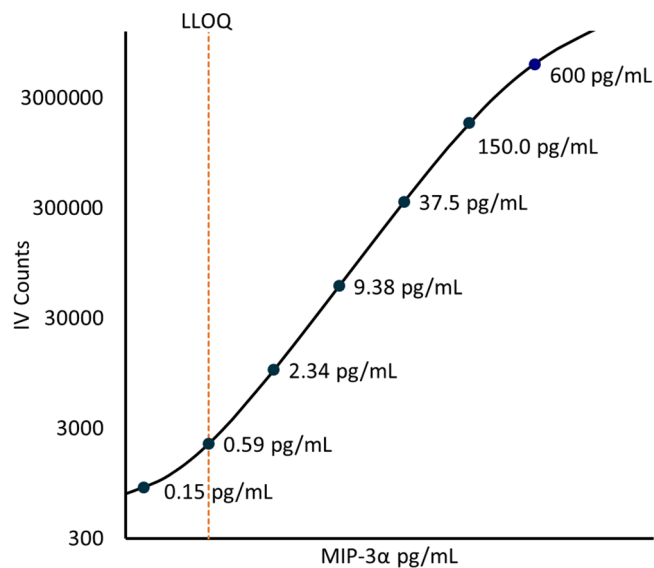


Description – MIP-3α

Macrophage Inflammatory Protein-3 alpha (MIP-3α), also known as C-C Motif Chemokine Ligand 20 (CCL20) is a 70 amino acid chemokine (molecular weight 8.0 kDa). MIP-3α is expressed in many tissues with the highest levels of expression found primarily in the liver, the lymph nodes and on peripheral blood lymphocytes. MIP-3α is upregulated by inflammatory cytokines such as Tumor Necrosis Factor Alpha (TNFα) and Interferon Gamma (IFNγ) while being strongly down regulated by IL-10. MIP-3α binds to target cells via the CCR6 receptor commonly found on Th-17 cells. MIP-3α serves as a strong chemoattractant for both lymphocytes and dendritic cells.

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 (1:4 Dilution) *	12.5 μL per measurement
---	-------------------------

*See Kit Instructions for details

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

Analytical LLOQ	0.5859 pg/mL
Functional LLOQ (x MRD)	2.344 pg/mL
LOD	0.0989 pg/mL

Assay Range 0 - 2400pg/mL

Endogenous Serum and Plasma Readings: Healthy EDTA plasma (n=6) and serum (n=6) samples were measured.

% Above LOD 100%

% Above LLOQ 88%

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.