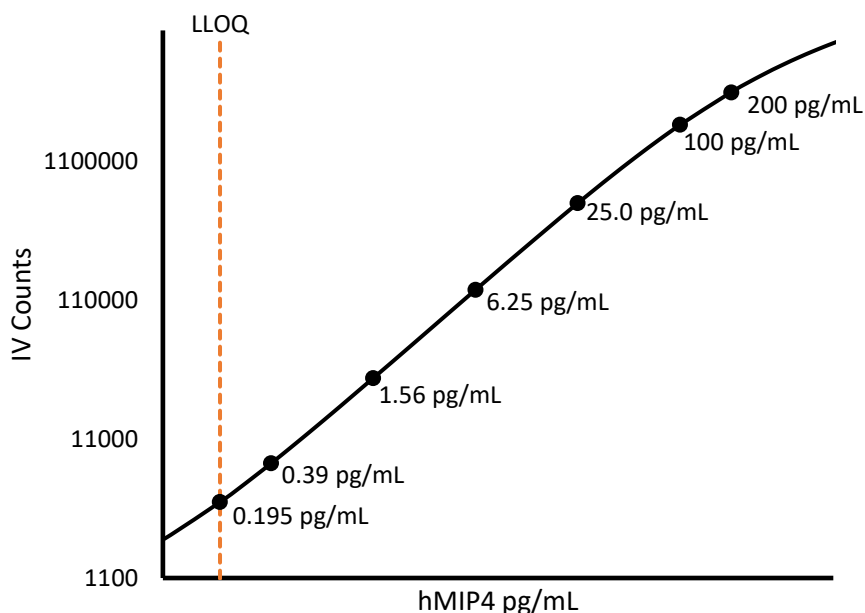


**Description – MIP-4 (PARC)**

MIP-4 (macrophage inflammatory protein-4) also known as CCL18 is a small chemokine of the CC chemokine family that encodes an 89 amino acid residue precursor protein. MIP-4 is produced and secreted mainly by the innate immune system. MIP-4 is highly expressed in lung and placenta but is not expressed in epidermal Langerhans cells. Recombinant MIP-4 has been shown to chemoattract naïve T cells but not monocytes or neutrophils.

**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)**

<b>Diluted Sample volume (1:5000 Dilution)*</b>	50 µL per measurement
---	-----------------------

\*See Kit Instructions for details

**Endogenous Serum and Plasma Readings:** Healthy EDTA plasma and serum samples (n=10) were measured.

<b>% Above LOD</b>	<b>100%</b>
<b>% Above LLOQ</b>	<b>100%</b>

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

<b>Analytical LLOQ</b>	<b>0.195 pg/mL</b>
<b>Functional LLOQ (x MRD)</b>	<b>975 pg/mL</b>
<b>LOD</b>	<b>70.3 fg/mL</b>
<b>Assay Range</b>	<b>0 – 1000 ng/mL</b>

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.