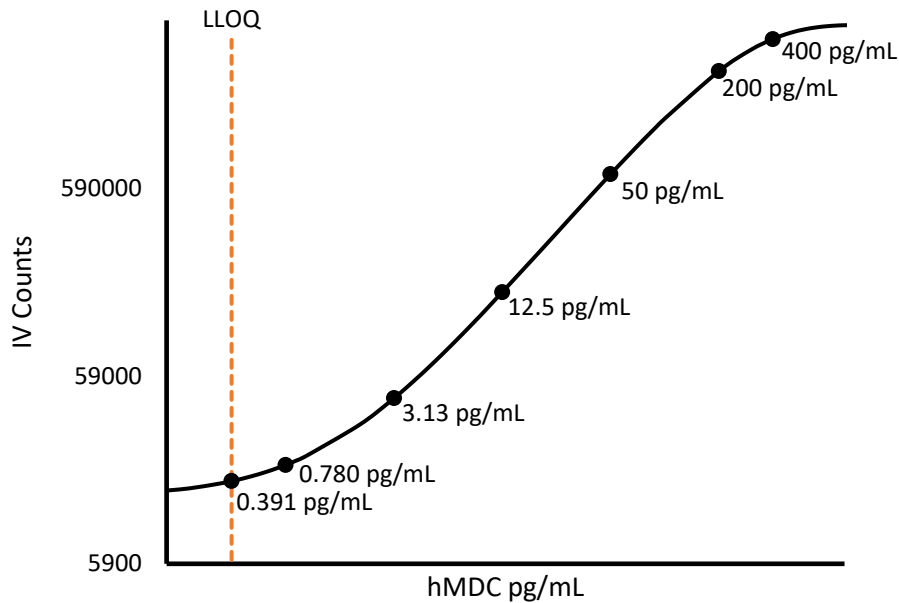


**Description – MDC**

MDC (Macrophage derived chemokine) also known as CCL22, is a CC chemokine that encodes a precursor protein of 93 amino acid residues. These proteins interact with cell surface chemokine receptors like CCR4. Recombinant mature MDC has been shown to induce chemotaxis in dendritic cells, IL-2 activated NK cells, and activated T lymphocytes.

**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 (1:20 Dilution)*</b>	50 µL per measurement
---	-----------------------

\*See Kit Instructions for details

**Endogenous Serum and Plasma Readings:** Healthy EDTA plasma (n=4) and serum (n=2) samples 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>0.391 pg/mL</b>
<b>Functional LLOQ (x MRD)</b>	<b>7.82 pg/mL</b>
<b>LOD</b>	<b>25.6 fg/mL</b>
<b>Assay Range</b>	<b>0 – 8000 pg/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.