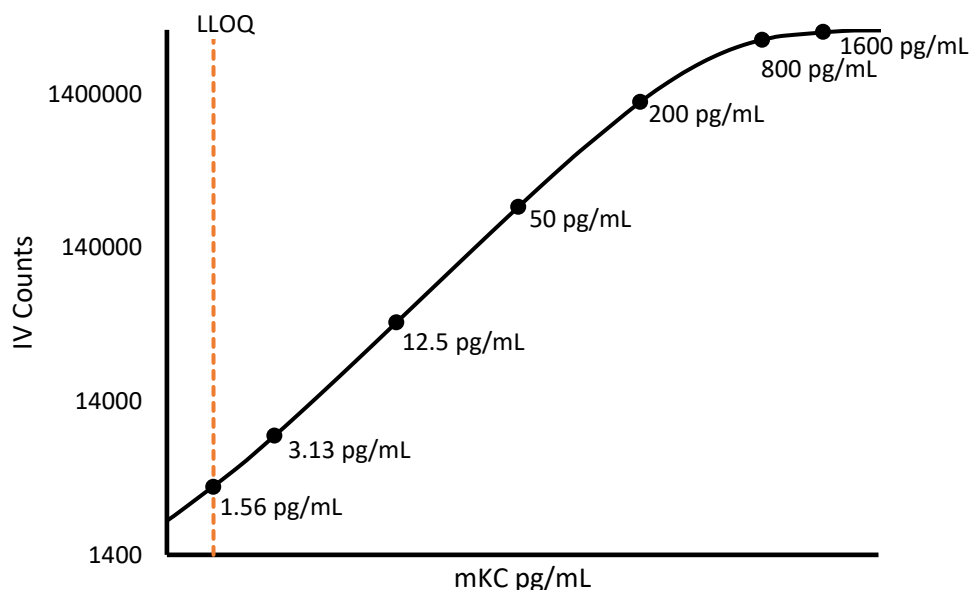


**Description – Mouse KC**

KC, better known as chemokine (C-X-C motif) ligand 1 (CXCL1), is a small cytokine that belongs to the CXC chemokine family. It is expressed by macrophages, neutrophils, and epithelial cells. CXCL1 plays a role in spinal cord development by inhibiting the migration of oligodendrocyte precursors and is involved in the processes of angiogenesis, arteriogenesis, inflammation, wound healing, and tumorigenesis. Dysregulation of CXCL1 has shown to be correlated with tumors, lung infections, pulmonary diseases and more.

**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 (1:2 Dilution)*</b>	50 µL per measurement
--	-----------------------

\*See Kit Instructions for details

**Endogenous Serum and Plasma Readings:** Healthy EDTA plasma and serum samples (n=8) from non-medicated, non-immunized mice 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>1.56 pg/mL</b>
<b>Functional LLOQ (x MRD)</b>	<b>3.12 pg/mL</b>
<b>LOD</b>	<b>25.2 fg/mL</b>
<b>Assay Range</b>	<b>0 – 3200 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.