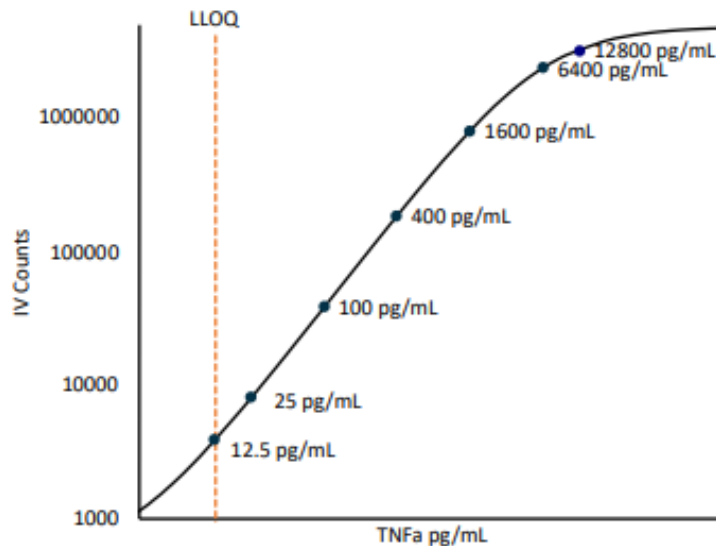


**Description – Mouse TNFα**

Tumor Necrosis Factor alpha (TNFα) is a polypeptide cytokine produced by monocytes and macrophages. The involvement of TNFα in several signal transduction pathways links the protein to such diverse functions as acute inflammation, apoptosis, septic shock, cellular proliferation, and differentiation. The clinical relevance of TNFα is from its association with numerous disease states including rheumatoid arthritis, cancer, cachexia, and Crohn’s disease.

**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 (n=10) and serum (n=10) samples from non-medicated, non-immunized mice were measured.

<b>% Above LOD</b>	<b>100%</b>
<b>% Above LLOQ</b>	<b>20%</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>12.5 pg/mL</b>
<b>Functional LLOQ (x MRD)</b>	<b>25.0 pg/mL</b>
<b>LOD</b>	<b>0.63 pg/mL</b>
<b>Assay Range</b>	<b>0 – 25,600 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.