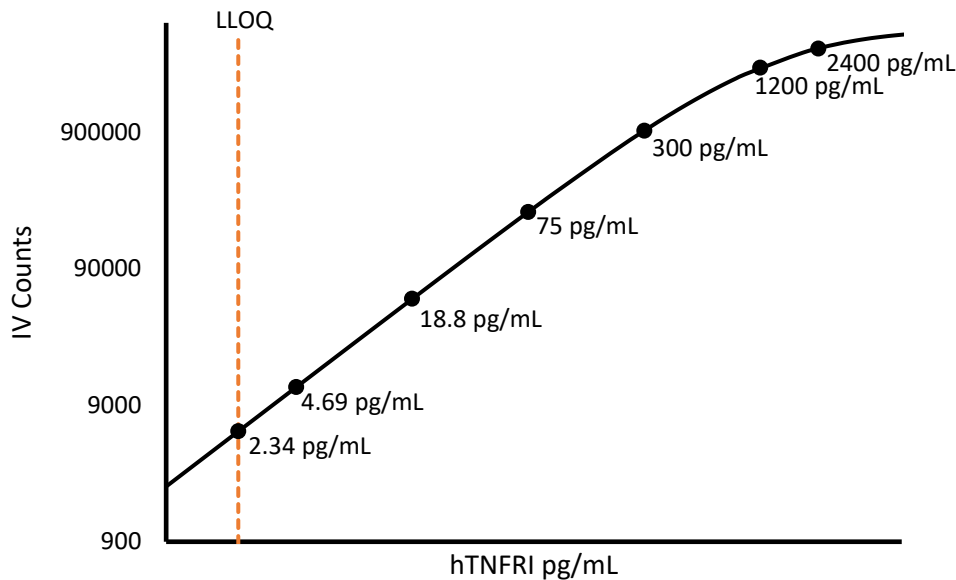


**Description – TNF-RI**

TNF-RI (Tumor necrosis factor receptor 1), also known as CD120a and TNFRSF1A, is a 21 kDa transmembrane protein that functions as a regulator of apoptosis and inflammation. TNF-RI is thought to mediate most of the cellular effects of TNF $\alpha$ . TNF-RI is also involved in neurotoxic processes.

**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 (1:25 Dilution)*</b>	50 $\mu$ L per measurement
---	----------------------------

\*See Kit Instructions for details

**Endogenous Serum and Plasma Readings:** Healthy EDTA plasma (n=4) and serum (n=4) 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>2.34 pg/mL</b>
<b>Functional LLOQ (x MRD)</b>	<b>58.5 pg/mL</b>
<b>LOD</b>	<b>0.208 pg/mL</b>
<b>Assay Range</b>	<b>0 – 60,000 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.