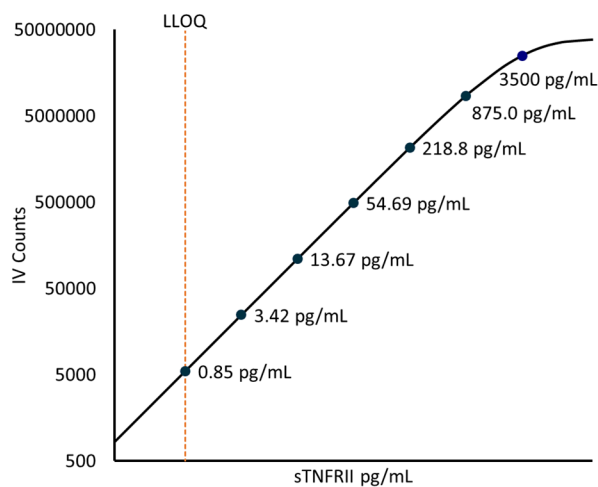


**Description – sTNFRII**

Tumor Necrosis Factor Receptor 2, also known as TNFR2 or CD120, is a type I transmembrane protein belonging to the TNF superfamily. Both TNFR1 and TNFR2 are co-expressed with TNFR1 ubiquitously expressed throughout many cell types. TNFR2 on the other hand is primarily expressed within immune cells and certain endothelial cells. Both receptors can also exist in soluble forms by proteolytic shedding of the extracellular portion of the receptors by the tumor necrosis factor alpha converting enzyme (TACE/ADAM17). The soluble forms of both TNFR1 and TNFR2 have been found to neutralize the activity of TNFα thus inhibiting the TNFα-mediated proinflammatory effects. Circulating sTNFR1 and sTNFR2 are elevated in many disease types including the two autoimmune inflammatory diseases ulcerative colitis 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 (1:8 Dilution)*</b>	6.25 µL per measurement
--	-------------------------

\*See Kit Instructions for details

**Endogenous Serum and Plasma Readings:** Healthy EDTA plasma (n=20) and serum (n=20) 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>0.855 pg/mL</b>
<b>Functional LLOQ (x MRD)</b>	<b>6.84 pg/mL</b>
<b>LOD</b>	<b>0.0965 pg/mL</b>
<b>Assay Range</b>	<b>0 – 28,000 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.