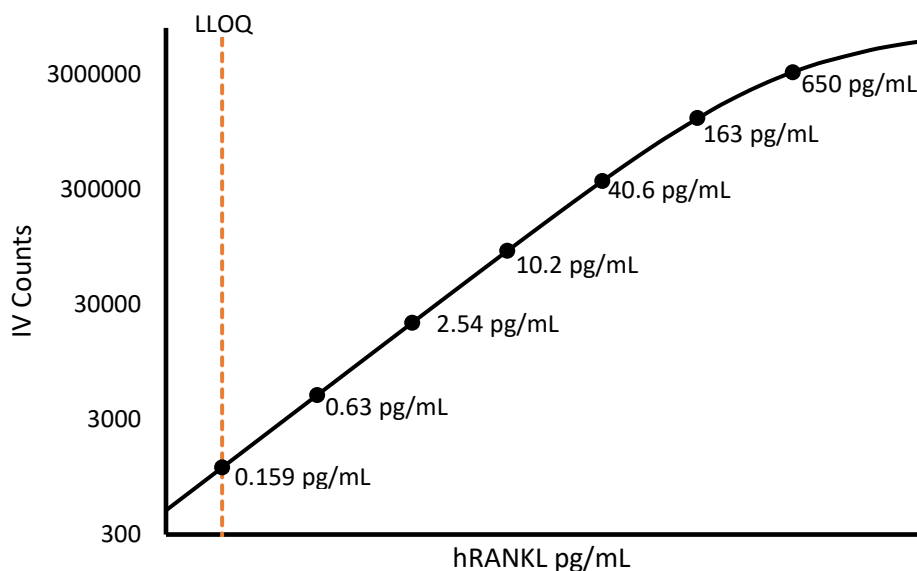


**Description – RANKL**

RANKL also known as TRANCE (TNF-related activation induced cytokine), is a type II membrane protein that is a member of the TNF (tumor necrosis factor) family. RANKL is also referred to as TNSF11 and functions as an inducer of osteoclastogenesis, lymph node organogenesis, enhancement of T cell growth, and dendritic cell function. Osteoprotegerin is a naturally occurring decoy receptor that counterbalances the effects of RANKL.

**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) were measured.

<b>% Above LOD</b>	<b>100%</b>
<b>% Above LLOQ</b>	<b>87.5%</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>0.159 pg/mL</b>
<b>Functional LLOQ (x MRD)</b>	<b>0.318 pg/mL</b>
<b>LOD</b>	<b>55.0 fg/mL</b>
<b>Assay Range</b>	<b>0 – 1300 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.