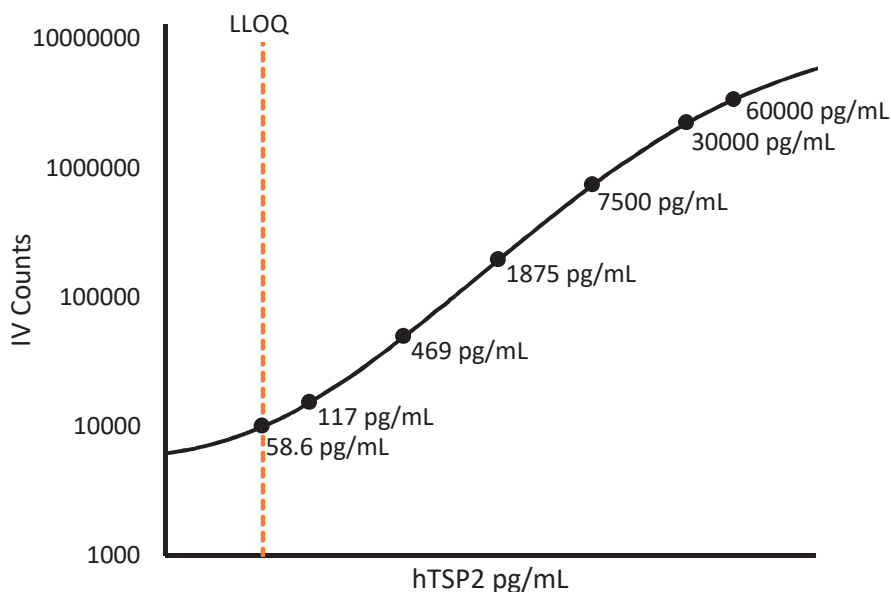


### Description – Thrombospondin-2 (TSP-2)

TSP-2 is a 150 kDa calcium-binding protein that regulates cellular interactions with extracellular matrix. TSP-2 is a potent inhibitor of tumor growth and angiogenesis. TSP-2 regulates collagen matrix formation by altering fibroblast behavior during development and in areas of tissue remodeling in the adult. TSP-2 is involved in cell adhesion and migration as well as modulating the cell surface properties of mesenchymal cells.

**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 (1:50 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>58.6 pg/mL</b>
<b>Functional LLOQ (x MRD)</b>	<b>2930 pg/mL</b>
<b>LOD</b>	<b>12.6 pg/mL</b>
<b>Assay Range</b>	<b>0 – 3000 ng/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.