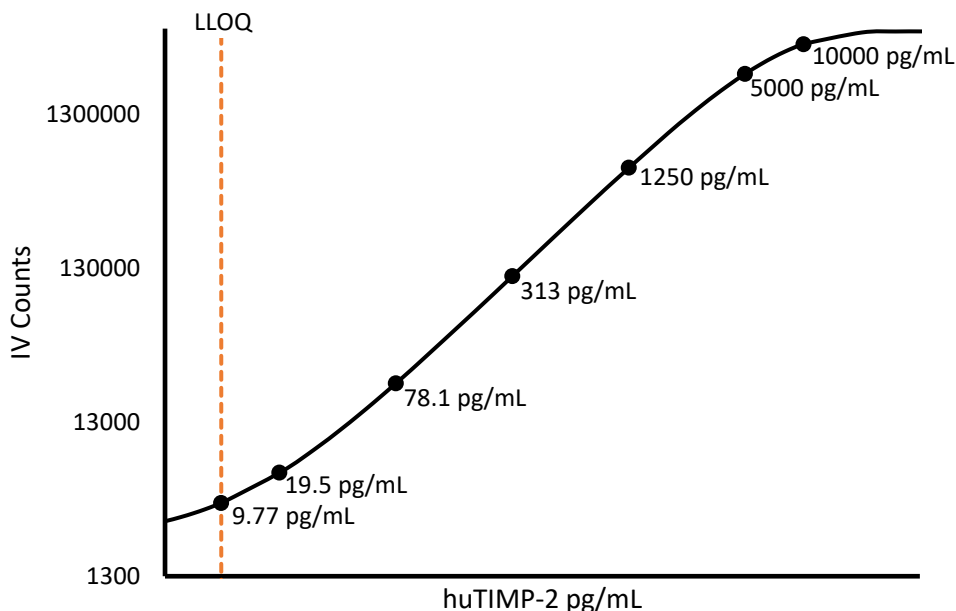


**Description – TIMP-2**

TIMP-2 (Tissue inhibitor of metalloproteinases-2) is a 21 kDa glycoprotein that regulates MMPs (matrix metalloproteinases). In addition to its inhibitory role against most of the known MMPs, TIMP-2 also has erythroid-potentiating and cell growth promoting activities. TIMP-2 may play a significant role in cancer development.

**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 (1:1000 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>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>9.77 pg/mL</b>
<b>Functional LLOQ (x MRD)</b>	<b>9770 pg/mL</b>
<b>LOD</b>	<b>4.63 pg/mL</b>
<b>Assay Range</b>	<b>0 – 10,000 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.