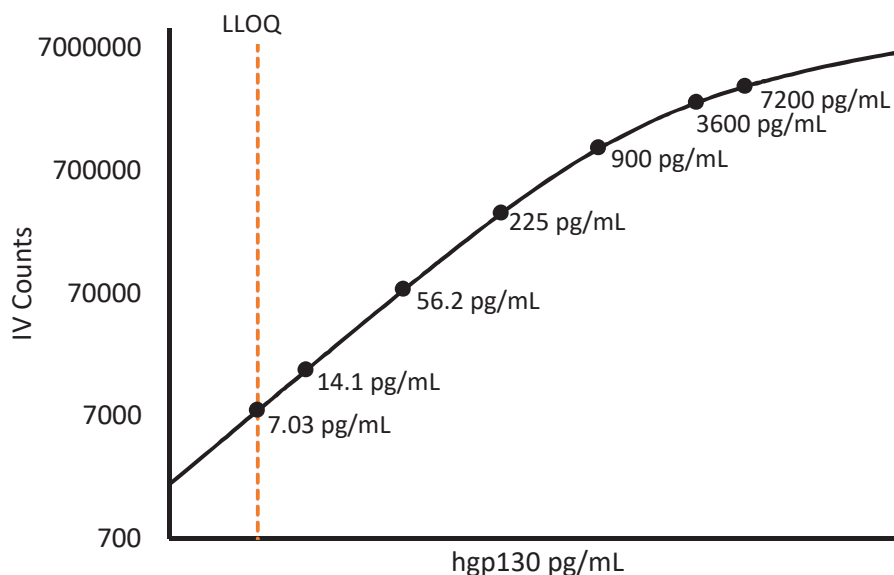


**Description – Glycoprotein 130 (gp130)**

Gp130 is a transmembrane protein within the IL-6 family of receptor cytokines. Gp130 forms signaling complexes to stimulate the production of anti-oxidant and anti-apoptotic proteins. Soluble gp130 has been shown to have inhibitory effects on IL-6 activities.

**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)**

Diluted Sample volume (1:1000 Dilution)*	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.

% Above LOD	100%
% Above LLOQ	100%

**Assay Range:** The upper end of the dynamic range is equal to the top calibrator concentration multiplied by MRD.

Analytical LLOQ	7.03 pg/mL
Functional LLOQ (x MRD)	7030 pg/mL
LOD	0.754 pg/mL
Assay Range	0 – 7200 ng/mL

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.