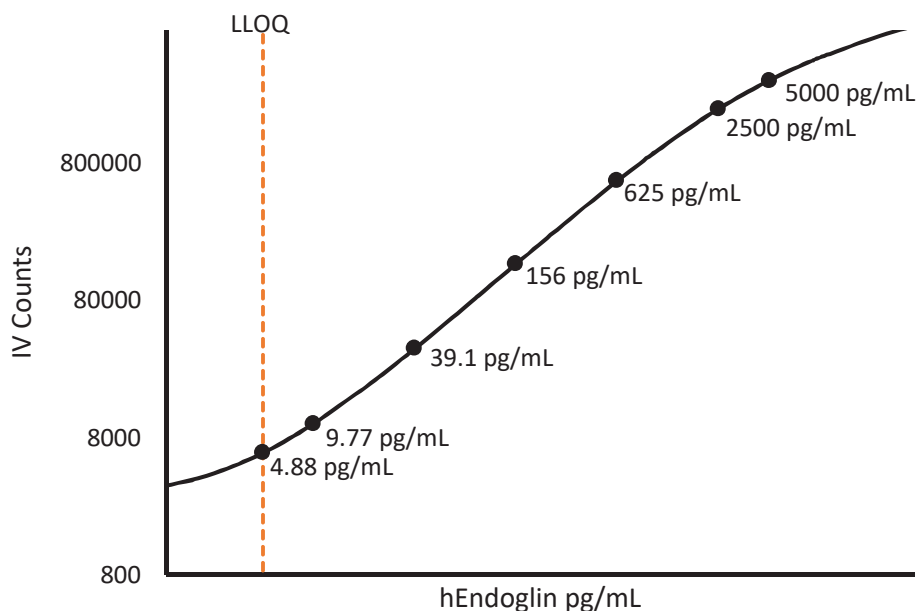


Description – Endoglin

Endoglin is a 180 kDA transmembrane glycoprotein highly expressed on proliferating vascular endothelial cells, chondrocytes, and syncytiotrophoblasts of term placenta. Due to its role in angiogenesis, it is a protein of interest for studying metastasis of cancer cells and tumor growth.

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 (1:100 Dilution)*	50 μ 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	4.88 pg/mL
Functional LLOQ (x MRD)	488 pg/mL
LOD	1.47 pg/mL
Assay Range	0 - 500 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.