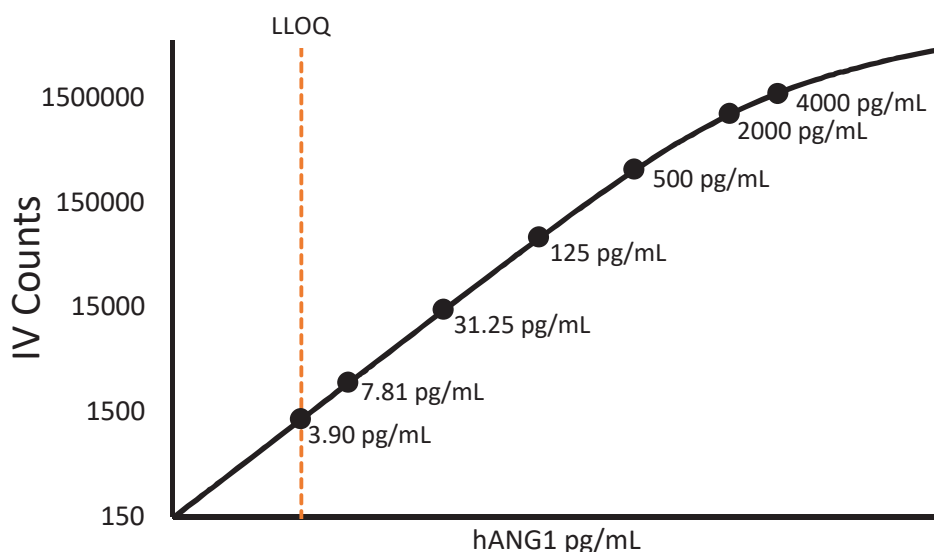


Description – Ang-1

Angiopoietin-1 (Ang-1) is a multimeric ligand involved in the angiopoietin-Tie signaling system that activates Tie2. It is predominantly expressed by endothelial cells and some smooth muscle cells, with increased expression in many inflammatory and angiogenic conditions, such as inflammation-induced vascular remodeling and tumor angiogenesis. Ang-1 expression has been documented in a range of human cancers, including glioblastoma, melanoma, prostate adenocarcinoma, and renal cell carcinoma.

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 (1:10 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	3.90 pg/mL
Functional LLOQ (x MRD)	39.0 pg/mL
LOD	1.99 pg/mL
Assay Range	0 – 40,000 pg/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.