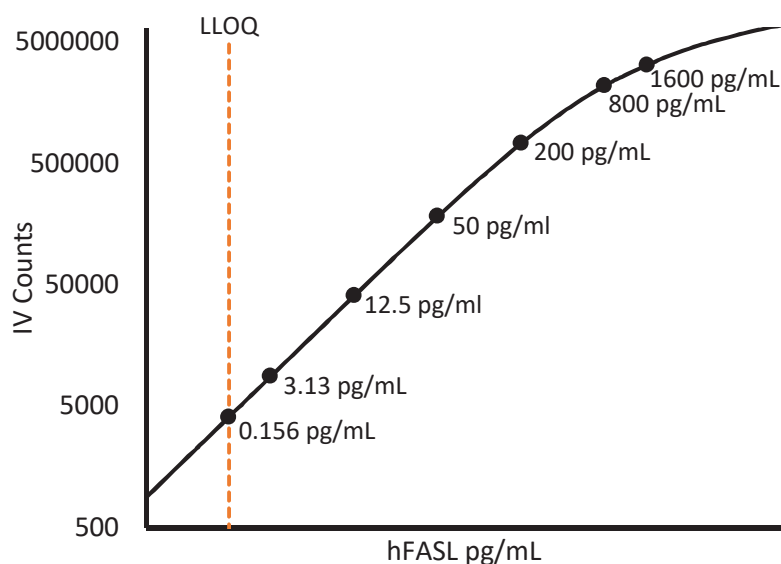


Description – Fas Ligand (FasL)

FasL is a 40 kDa type-II transmembrane protein also known as CD178 or TNFSF6. FasL plays a crucial role in the development, homeostasis, and function of the immune system due to its ability to induce apoptosis in target cells by binding to its receptor Fas/CD95 on adjacent cells. Because of this apoptotic property, FasL may be a protein of interest for cancer research.

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 (1:2 Dilution)*	50 μ L per measurement
--	----------------------------

*See Kit Instructions for details

Endogenous Serum and Plasma Readings: Healthy EDTA plasma (n=6) and serum (n=8) 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	1.56 pg/mL
Functional LLOQ (x MRD)	3.12 pg/mL
LOD	0.158 pg/mL
Assay Range	0 – 3200 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.