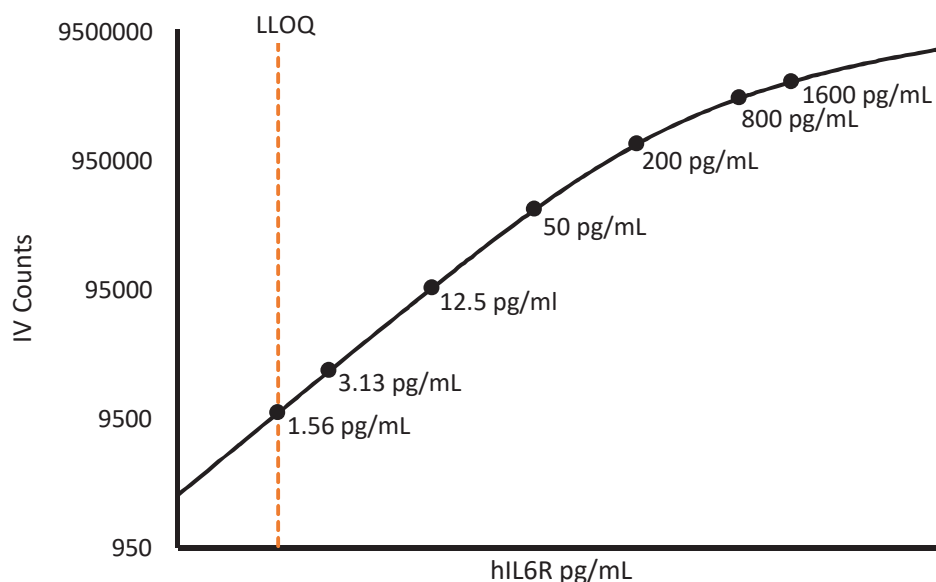


### Description – Interleukin-6 Receptor (IL-6R)

Interleukin-6 Receptor (IL-6R) is a multifunctional factor interleukin which functions by binding to a high-affinity receptor complex consisting of two membrane glycoproteins. The two components are a signal-transducing gp130 and a component receptor that binds IL-6 with low affinity, as known as IL-6R alpha. IL-6R is a subunit of the receptor for IL-6. Low concentration of a soluble form of IL-6R acts as an agonist of IL-6 activity. It interacts with Interleukin 6 and ciliary neurotrophic factor. Soluble interleukin-6 receptor plays a role in immunoregulation and cellular processes.

**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	1.56 pg/mL
Functional LLOQ (x MRD)	1.56 ng/mL
LOD	0.0771 pg/mL
Assay Range	0 – 1600 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.