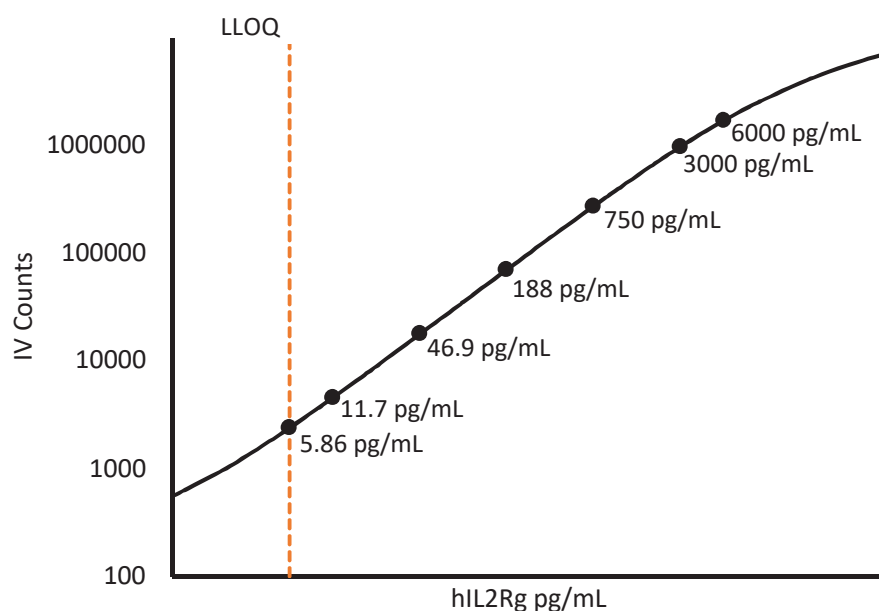


### Description – Interleukin-2 Receptor Gamma (IL-2RG)

Interleukin-2 receptor gamma (IL-2RG) is a glycoprotein of the type I cytokine receptor family expressed on most lymphocyte populations. The gamma chain is the signaling subunit of the receptor complex. Lymphocytes that express this chain can form functional receptors for these cytokine proteins, which transmit signals from one cell to another and direct programs of cellular differentiation. Mutations in IL-2RG have been associated with immunodeficiency disorders.

**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)

<b>Diluted Sample volume (1:2 Dilution)*</b>	50 µL per measurement
--	--------------------------

\*See Kit Instructions for details

**Endogenous Serum and Plasma Readings:** Healthy EDTA plasma (n=4) and serum (n=4) samples were measured.

<b>% Above LOD</b>	<b>100%</b>
<b>% Above LLOQ</b>	<b>12.5%</b>

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

<b>Analytical LLOQ</b>	<b>5.86 pg/mL</b>
<b>Functional LLOQ (x MRD)</b>	<b>11.7 pg/mL</b>
<b>LOD</b>	<b>1.69 pg/mL</b>
<b>Assay Range</b>	<b>0 – 12,000 pg/mL</b>

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.