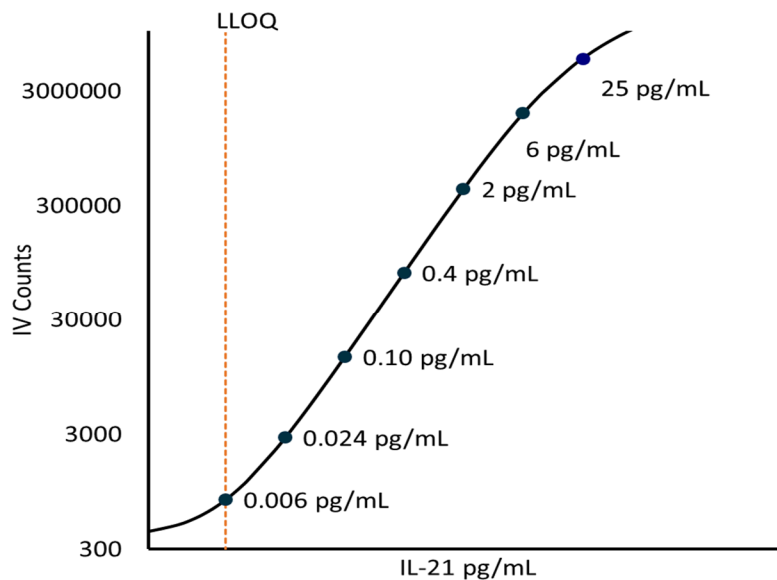


**Description – IL-21**

Interleukin 21 (IL-21) is a 133 amino acid (molecular weight 15.4 kDa) cytokine that plays a key role in enhancing antigen specific immune responses. It is primarily secreted by CD4+ T cells after antigenic stimulation. IL-21 binds to a heterodimeric receptor composed of IL-21R and a common gamma chain subunit used by other cytokines such as IL-2 and IL-4. IL-21 has been shown to promote the activation and proliferation of CD8+ T cells, B cell maturation and differentiation, as well as driving TH-17 cell based immune responses.

**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 (1:4 Dilution)*</b> | 12.5 µL per measurement |
|--|-------------------------|

\*See Kit Instructions for details

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

|                                |              |
|--------------------------------|--------------|
| <b>Analytical LLOQ</b>         | 0.0061 pg/mL |
| <b>Functional LLOQ (x MRD)</b> | 0.0244 pg/mL |
| <b>LOD</b>                     | 0.0008 pg/mL |

|                    |               |
|--------------------|---------------|
| <b>Assay Range</b> | 0 - 100 pg/mL |
|--------------------|---------------|

**Endogenous Serum and Plasma Readings:** Healthy EDTA plasma (n=8) and serum (n=8) samples were measured.

|                     |       |
|---------------------|-------|
| <b>% Above LOD</b>  | 93.8% |
| <b>% Above LLOQ</b> | 31.3% |

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.