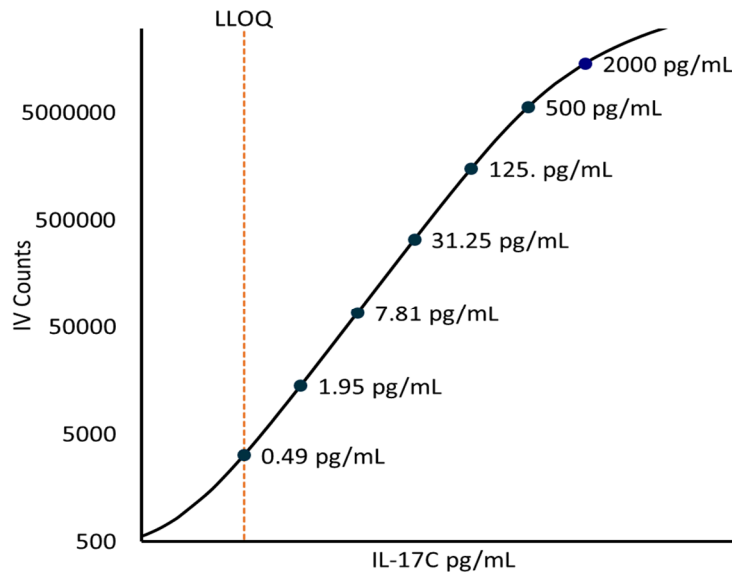


**Description – IL-17C**

Interleukin 17C (IL-17C) is a glycosylated 179 amino acid cytokine (molecular weight 20 kDa) which is expressed as a disulfide linked dimer. IL-17C is member of a family of structurally related proteins (IL-17A – IL-17F) which play pivotal roles in inflammatory and immunoregulatory responses. In contrast to other members of the family IL-17C is expressed primarily in non-lymphoid cells such as keratinocytes and epithelial cells. It binds to the receptor IL-17RE which is expressed on TH-17 cells. This cytokine is thought to play important role in TH-17 mediated autoimmunity by stimulating the production of IL-17A, IL-22 and MIP-3α.

**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>	<b>0.4883 pg/mL</b>
<b>Functional LLOQ (x MRD)</b>	<b>1.953 pg/mL</b>
<b>LOD</b>	<b>0.0429 pg/mL</b>

<b>Assay Range</b>	<b>0 - 8000 pg/mL</b>
--------------------	-----------------------

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

<b>% Above LOD</b>	<b>100%</b>
<b>% Above LLOQ</b>	<b>88%</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.