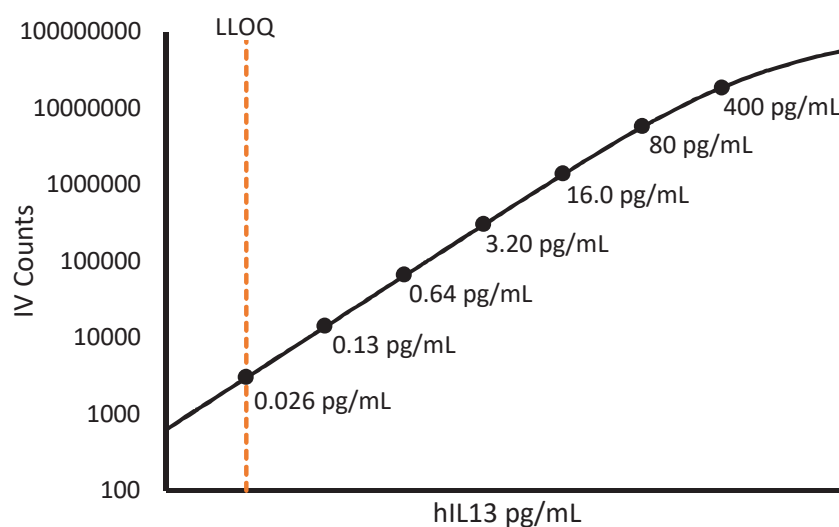


### Description – Interleukin-13 (IL-13)

Interleukin 13 (IL-13) is a cytokine of 111 amino acids (molecular weight 15.8 kDa) whose major roles include down-modulation of macrophage activity (lowering the production of pro-inflammatory cytokines) and mediation of allergic responses. It is secreted by many cell types, but primarily by activated T-cells, in particular T-helper type 2 cells. IL-13 affects immune cells in a manner similar to IL-4 but is more associated with physiological changes induced by allergic inflammation. The effects of IL-13 are induced through a receptor that includes the alpha chain of the IL-4 receptor and at least one or two known IL-13 specific binding chains. IL-13 is the central mediator of allergic asthma, where it regulates eosinophilic inflammation, mucus secretion, and airway hyperresponsiveness.

**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 $\mu$ L per measurement
--	-------------------------------

\*See Kit Instructions for details

**Endogenous Serum and Plasma Readings:** Healthy EDTA plasma and serum samples (n=6) 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	0.0256 pg/mL
Functional LLOQ (x MRD)	0.0512 pg/mL
LOD	0.00876 pg/mL
Assay Range	0 – 800 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.