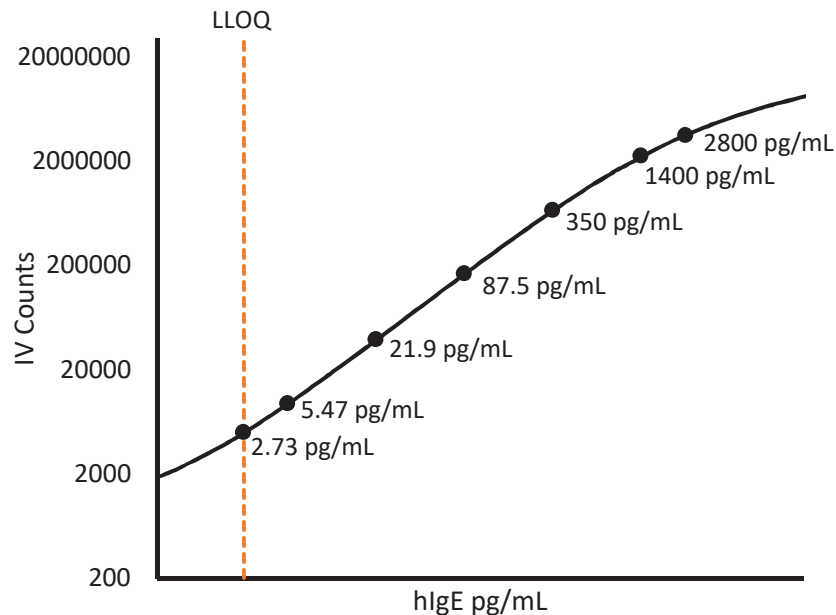


Description – Immunoglobulin E (IgE)

IgE is an antibody synthesized by plasma cells and made up of two heavy chains and two light chains. IgE primarily functions to defend against parasitic invasion and is responsible for allergic reactions. IgE mediates inflammatory reactions as a response to allergens, releasing chemicals like histamine, leukotrienes, and certain interleukins.

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 (1:2000 Dilution)*	50 µL per measurement
---	--------------------------

*See Kit Instructions for details

Endogenous Serum and Plasma Readings: Healthy EDTA plasma and serum samples (n=8) 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	2.734 pg/mL
Functional LLOQ (x MRD)	5468 pg/mL
LOD	0.216 pg/mL
Assay Range	0 – 5600 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.