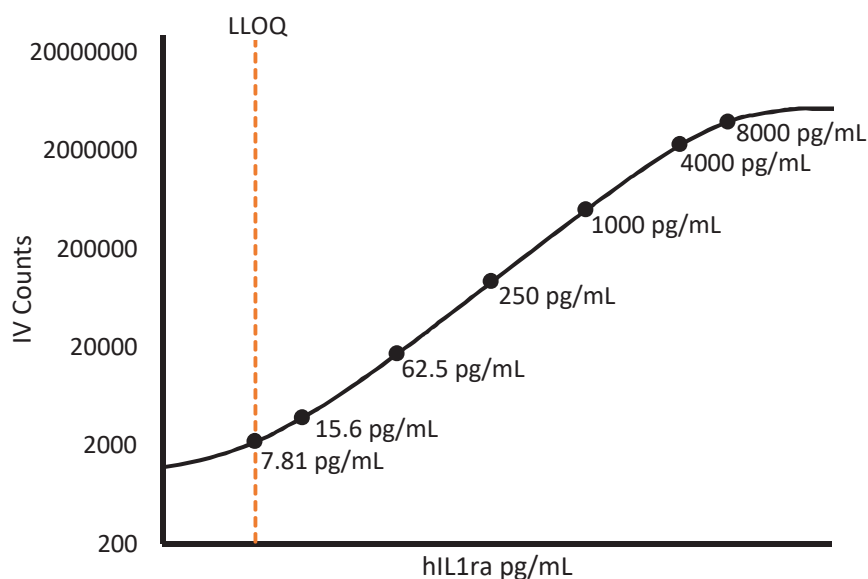


Description – Interleukin-1 receptor antagonist (IL-1RA)

Interleukin-1 receptor antagonist (IL-1RA) is an anti-inflammatory cytokine that inhibits the inflammatory responses of IL-1 through competitive binding of IL-1 type I receptors. IL-1RA is critical to the development and progression of diseases such as sepsis and rheumatoid arthritis. A recombinant form of IL-1RA has been investigated as a treatment for rheumatoid arthritis and a potential treatment for brain injury.

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 µL per measurement
--	--------------------------

*See Kit Instructions for details

Endogenous Serum and Plasma Readings: Healthy EDTA plasma (n=5) and serum (n=5) samples 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	7.81 pg/mL
Functional LLOQ (x MRD)	15.6 pg/mL
LOD	4.04 pg/mL
Assay Range	0 – 16,000 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.