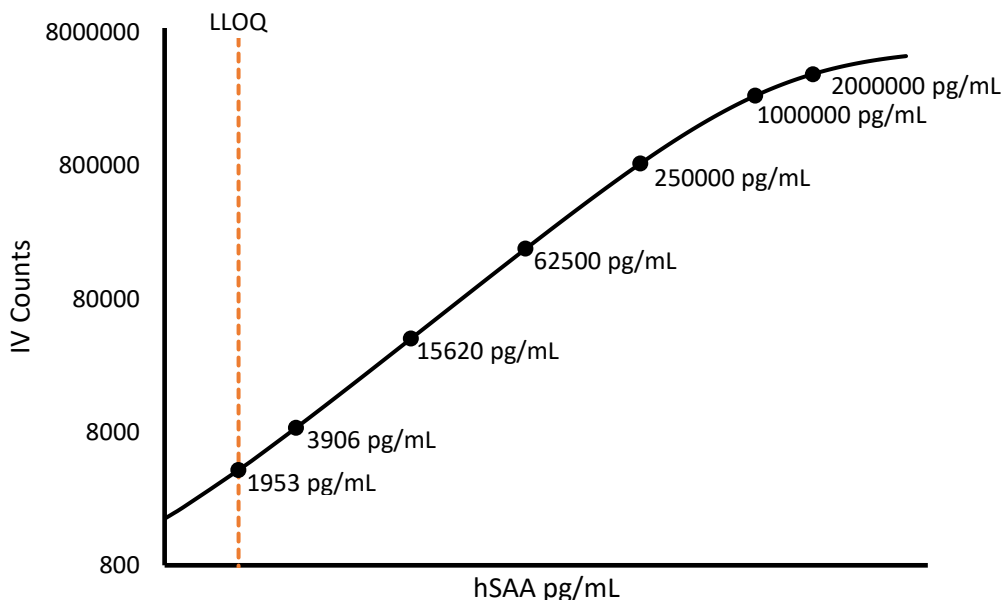


Description – SAA

SAA (serum amyloid A) are multifunctional apolipoproteins produced by hepatocytes in response to proinflammatory cytokines. SAA has been shown to have four different isoforms in humans which are expressed constitutively at different levels in response to inflammatory stimuli.

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 (1:200 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	1953 pg/mL
Functional LLOQ (x MRD)	391 ng/mL
LOD	65.4 pg/mL
Assay Range	0 – 400 µg/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.