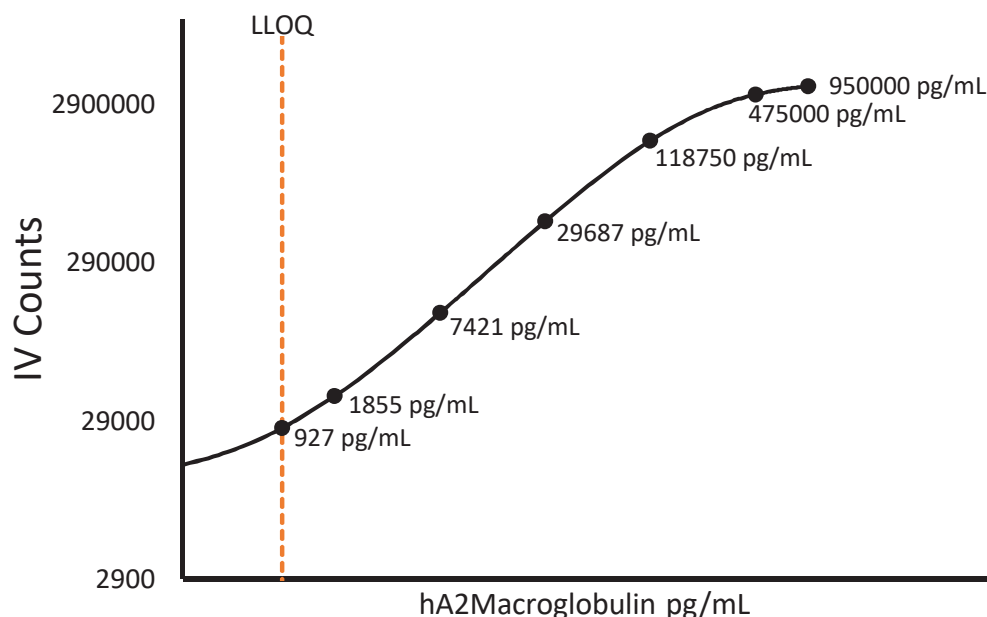


Description – Alpha-2 Macroglobulin

Alpha-2 macroglobulin is a large protein found in serum and plasma and synthesized primarily by the liver. It functions as a protein inhibitor in a variety of biological functions including coagulation and fibrinolysis. Levels of alpha-2 macroglobulin in the blood have been associated with nephrotic syndrome and Alzheimer's disease.

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 (1:100,000 Dilution)*	50 µL per measurement
--	--------------------------

*See Kit Instructions for details

Endogenous Serum and Plasma Readings: Healthy EDTA plasma (n=4) and serum (n=4) 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	927 pg/mL
Functional LLOQ (x MRD)	92.7 µg/mL
LOD	10.1 pg/mL
Assay Range	0 – 95,000 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.