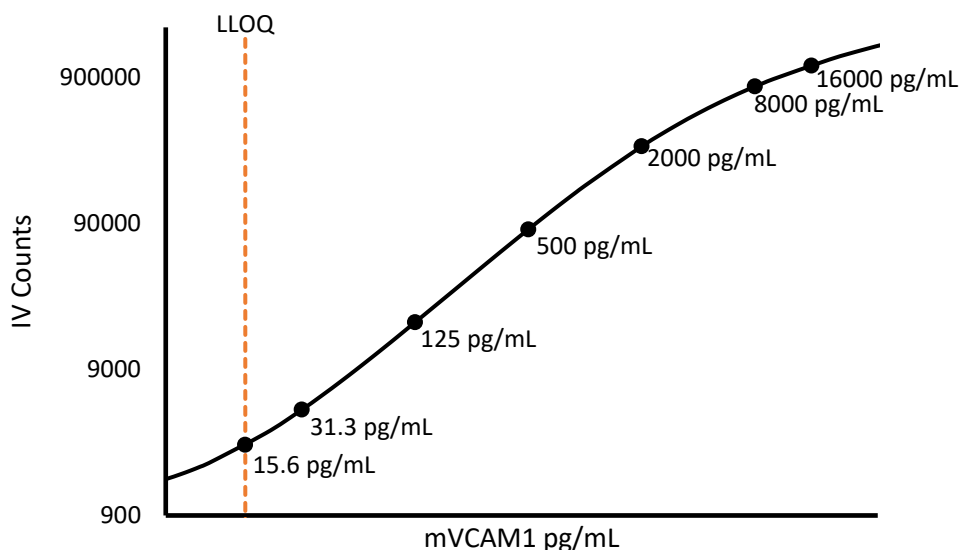


Description – Mouse VCAM-1

Vascular Cellular Adhesion Molecule 1 (VCAM-1) also known as CD106 is a 110 kDa transmembrane glycoprotein expressed exclusively on cytokine-activated vascular endothelium. VCAM-1 protein mediates the adhesion of lymphocytes, monocytes, eosinophils, and basophils to vascular endothelium, and also functions in leukocyte-endothelial cell signal transduction. VCAM-1 has been shown to play an important role in the development of atherosclerosis and rheumatoid arthritis, and soluble VCAM-1 is a useful biomarker for many inflammatory diseases.

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 and serum samples (n=8) from non-medicated, non-immunized mice 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	15.6 pg/mL
Functional LLOQ (x MRD)	31.2 pg/mL
LOD	2.29 pg/mL
Assay Range	0 – 32,000 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.