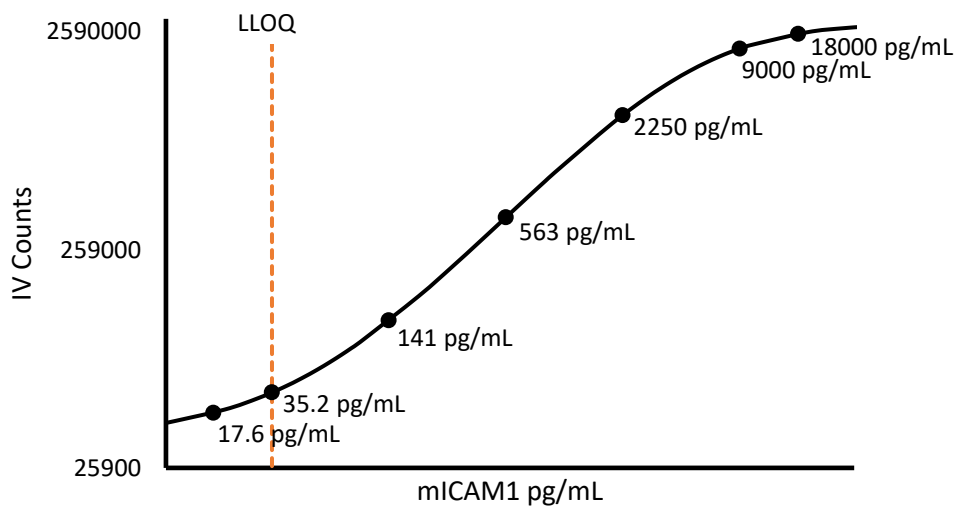


Description – Mouse ICAM-1

Intercellular adhesion molecule-1 (ICAM-1; also known as CD54) is a 90kDa member of the immunoglobulin (Ig) superfamily and plays a central role in leukocyte trafficking, activation of lymphocytes, and numerous additional immune functions. ICAM-1 is constitutively expressed on many immune, endothelial and epithelial cell types, and its expression is upregulated in inflammatory conditions, specifically by TNF- α , IL-1 β , IFN- γ and other cytokines. By interacting with lymphocyte function associated molecule (LFA)-1 and others, ICAM-1 acts as a leukocyte adhesion molecule. ICAM-1 is highly expressed in inflammatory conditions, chronic diseases and a number of malignancies.

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 μ L 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 | 35.2 pg/mL |
| Functional LLOQ (x MRD) | 70.4 pg/mL |
| LOD | 8.73 pg/mL |
| Assay Range | 0 – 36,000 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.