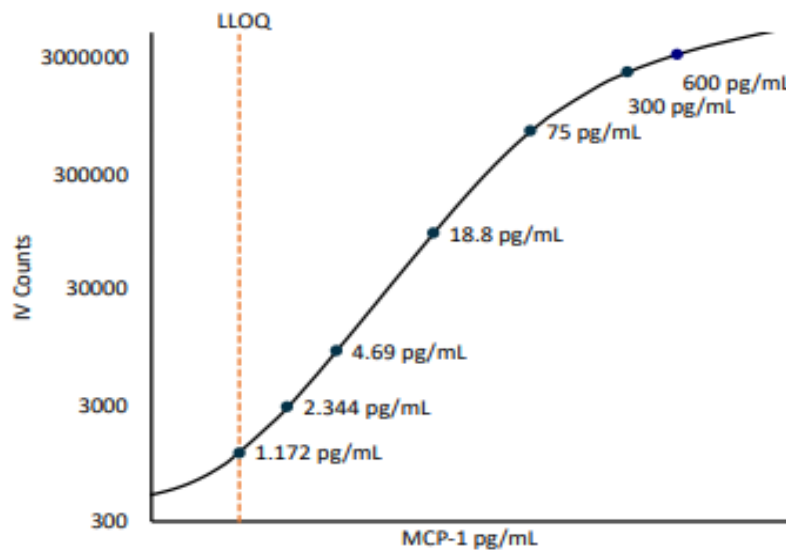


Description – MCP-1

Monocyte chemoattractant protein-1 (MCP-1) is a 13 kDa chemokine that regulates migration and infiltration of monocytes/macrophages. The idiopathic inflammatory myopathies, including dermatomyositis (DM), polymyositis and inclusion body myositis, are a group of autoimmune diseases characterized by chronic lymphocytic and macrophagic infiltration in muscle. MCP-1 is believed to play a major role in the recruitment of these cells, especially in DM. Plasma levels of MCP-1 have also been shown to be increased in patients with high viral load compared with HIV-seropositive subjects with undetectable plasma viral RNA and healthy controls.

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 (1:10 Dilution)*	50 µL per measurement
---	-----------------------

*See Kit Instructions for details

Endogenous Serum and Plasma Readings: Healthy EDTA plasma (n=20) and serum (n=20) 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	1.172 pg/mL
Functional LLOQ (x MRD)	11.719 pg/mL
LOD	0.511 pg/mL
Assay Range	0 – 6000 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.