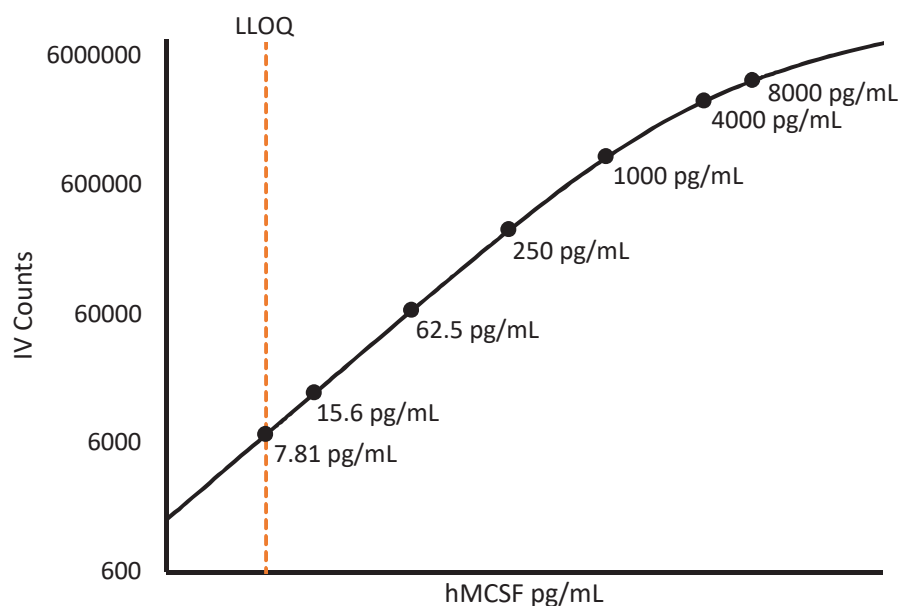


Description – Macrophage Colony-Stimulating Factor (M-CSF)

M-CSF is a secreted cytokine that is a primary regulator of macrophage survival, proliferation, and differentiation. M-CSF is responsible for the differentiation of HSCs to macrophages and other similar cell types, and is essential for the survival and proliferation of osteoclast progenitors. M-CSF is clinically used in the treatment of infection, malignancies, and atherosclerosis and in facilitating hematopoietic recovery after bone marrow transplantation.

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 (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	7.81 pg/mL
Functional LLOQ (x MRD)	15.6 pg/mL
LOD	0.896 pg/mL
Assay Range	0 – 16,000 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.