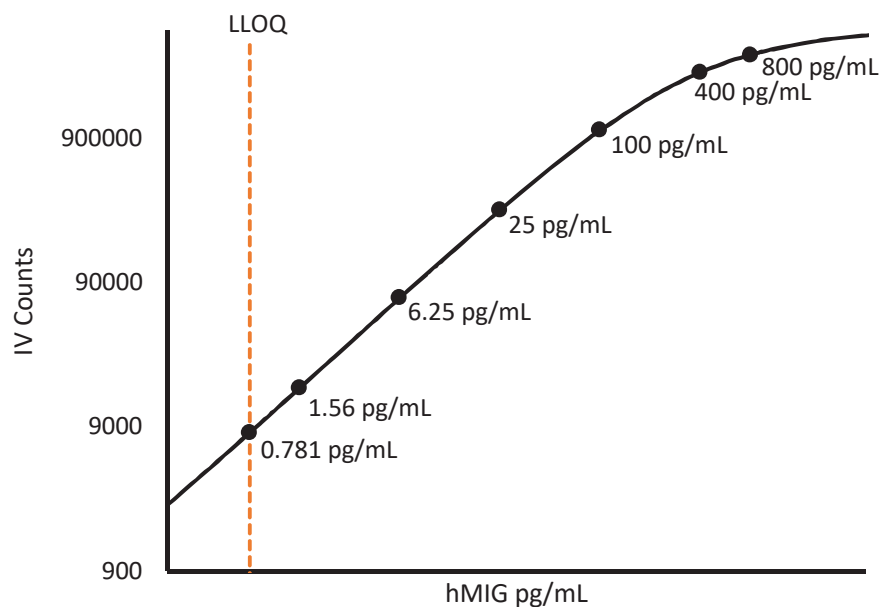


Description – Monokine induced by gamma interferon (MIG)

Monokine induced by interferon gamma (MIG), also known as CXCL9, is a ~12 kDa cytokine belonging to the subfamily of CXC chemokines. MIG is produced by IFN- γ stimulated monocytes, macrophages, and endothelial cells. MIG is involved in inhibiting tumor growth and colony formation of hematopoietic progenitors.

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 (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	0.781 pg/mL
Functional LLOQ (x MRD)	1.56 pg/mL
LOD	89.0 fg/mL
Assay Range	0 – 1600 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.