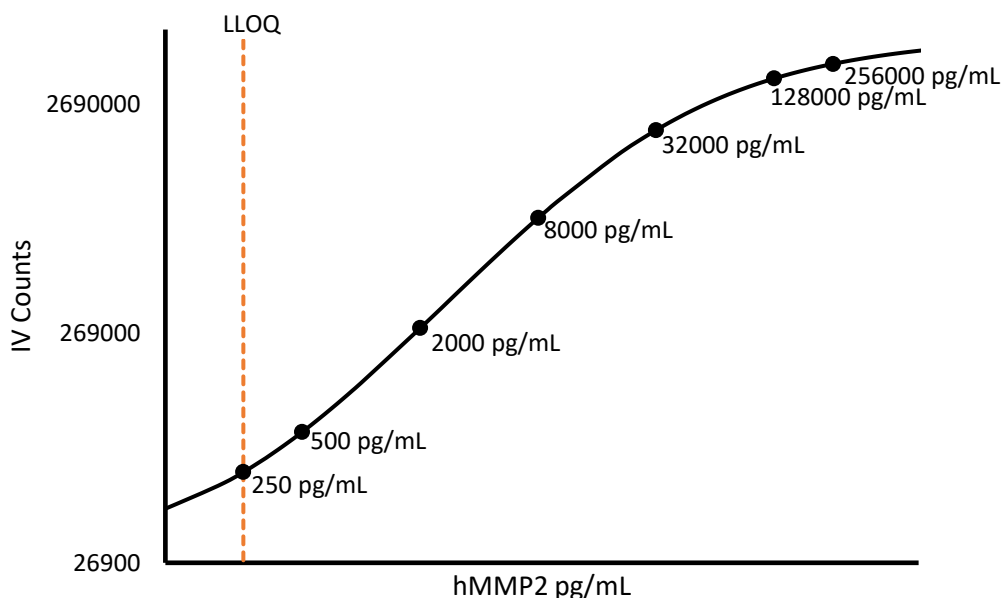


Description – MMP-2

MMP-2 (matrix metalloproteinase-2), also known as gelatinase A, is a 72 kDa enzyme which is encoded by the MMP2 gene. Its main function is to degrade collagens to switch on different patterning mechanisms for VEGF-C. It can also degrade matrix proteins decorin, elastin, and fibronectin. MMP-2 modulates the functions of many vasoactive and proinflammatory molecules such as Adrenomedullin, MCP-3, and SDF-1.

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

*See Kit Instructions for details

Endogenous Serum and Plasma Readings: Healthy EDTA plasma and serum samples (n=8) 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	250 pg/mL
Functional LLOQ (x MRD)	12,500 pg/mL
LOD	68.0 pg/mL
Assay Range	0 – 12,800 ng/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.