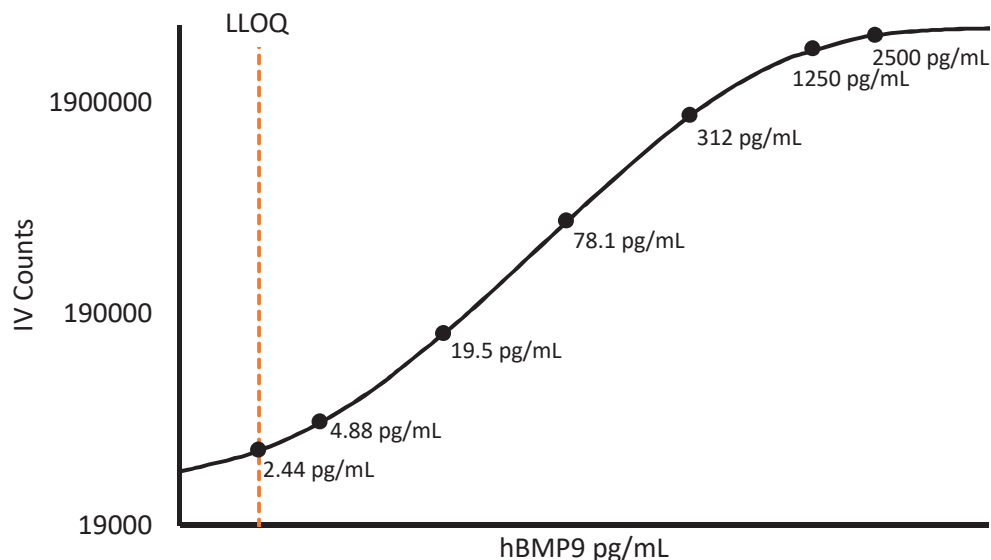


### Description – BMP-9

Bone morphogenetic protein-9 (BMP-9), also known as growth and differentiation factor 2 (GDF-2), is a 12kDa protein that regulates the development and function of numerous embryonal and adult tissues. BMP-9 is a member of the TGF- $\beta$  superfamily which initiates signals through heterodimeric complexes.

**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 $\mu$ 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	2.44 pg/mL
Functional LLOQ (x MRD)	4.88 pg/mL
LOD	0.634 pg/mL
Assay Range	0 – 5000 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.