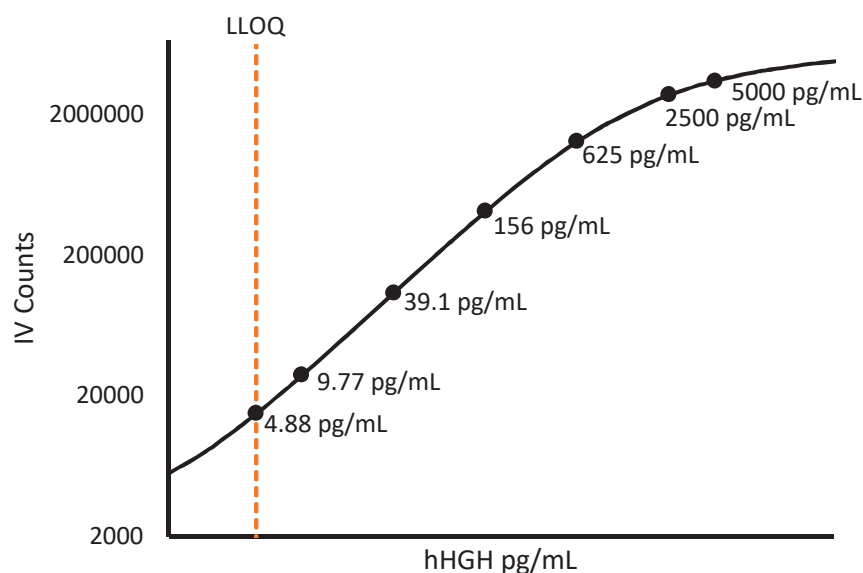


Description – Human Growth Hormone (HGH)

Human growth hormone (HGH) is a peptide produced by the pituitary gland and stored in secretory granules. HGH stimulates the liver and other tissues to produce IGF-1, which regulates growth and metabolism. HGH has also been shown to have direct effects on growth that is independent of IGF-1. HGH can act directly on various cell types to induce lipolysis, lactation, amino acid uptake, and protein synthesis.

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 serum (n=10) samples were measured.

% Above LOD	100%
% Above LLOQ	90%

Assay Range: The upper end of the dynamic range is equal to the top calibrator concentration multiplied by MRD.

Analytical LLOQ	4.88 pg/mL
Functional LLOQ (x MRD)	9.76 pg/mL
LOD	0.545 pg/mL
Assay Range	0 – 10 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.