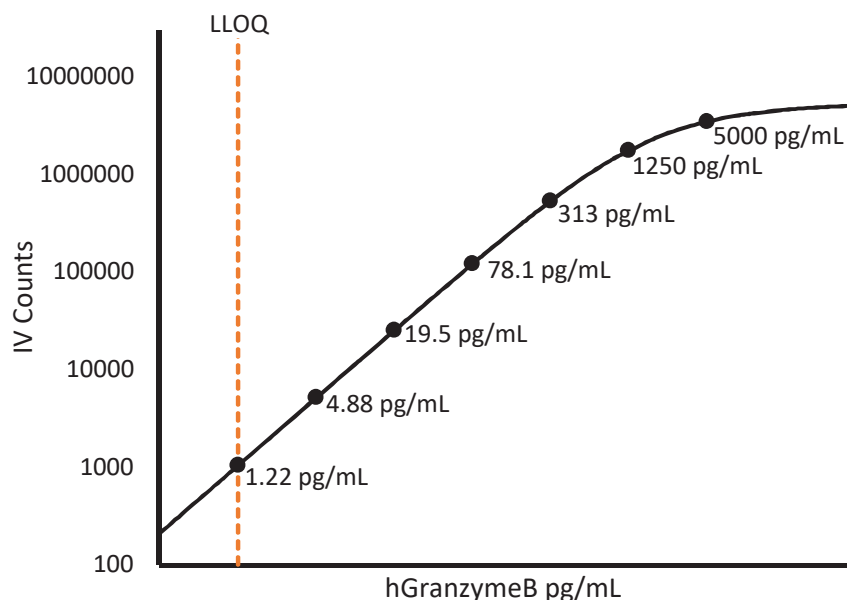


Description – Granzyme B

Granzyme B is a member of the granzyme family of serine proteases found specifically in the cytotoxic granules of cytotoxic T lymphocytes and natural killer cells. Granzyme B plays an important role in mediating apoptosis in target cells, inducing inflammation by stimulating cytokine release, and remodeling extracellular matrix.

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=5) and serum (n=5) samples were measured.

% Above LOD	100%
% Above LLOQ	30%

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

Analytical LLOQ	1.22 pg/mL
Functional LLOQ (x MRD)	2.44 pg/mL
LOD	0.547 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.