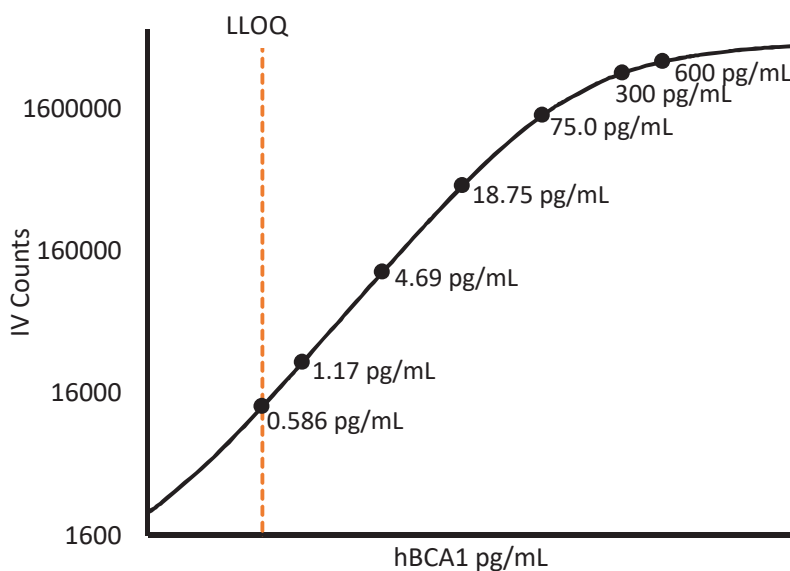


### Description – BCA-1 (CXCL13)

The CXCL13 (CXC Ligand 13) protein, also known as B cell attracting chemokine 1 (BCA-1) or B lymphocyte chemo-attractant (BLC) is a chemokine ligand. It is expressed in secondary lymphoid tissue by stromal cells within B-cell follicles and by follicular dendritic cells, macrophages, and germinal center T cells. CXCL13 modulates homeostatic immune cell trafficking and may become up-regulated under pathological conditions.

**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=5) and serum (n=5) 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	0.586 pg/mL
Functional LLOQ (x MRD)	1.17 pg/mL
LOD	0.0085 pg/mL
Assay Range	0 – 1,200 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.