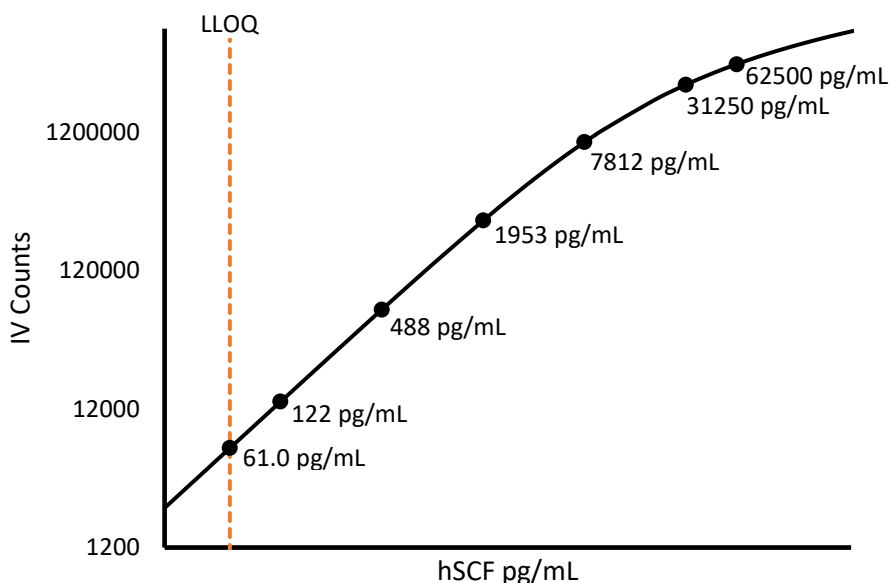


Description – SCF (KIT-Ligand)

SCF (Stem cell factor), also known as MGF (mast cell growth factor), SF (steel factor) and KL (kit ligand), is a 18.5 kDa transmembrane glycoprotein. SCF promotes survival, differentiation, and mobilization of multiple cell types including lymphoid, germ cell, and melanocyte progenitors. SCF is a primary growth and activation factor for mast cells and eosinophils.

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 µL 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	61.0 pg/mL
Functional LLOQ (x MRD)	122 pg/mL
LOD	5.84 pg/mL
Assay Range	0 – 125 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.