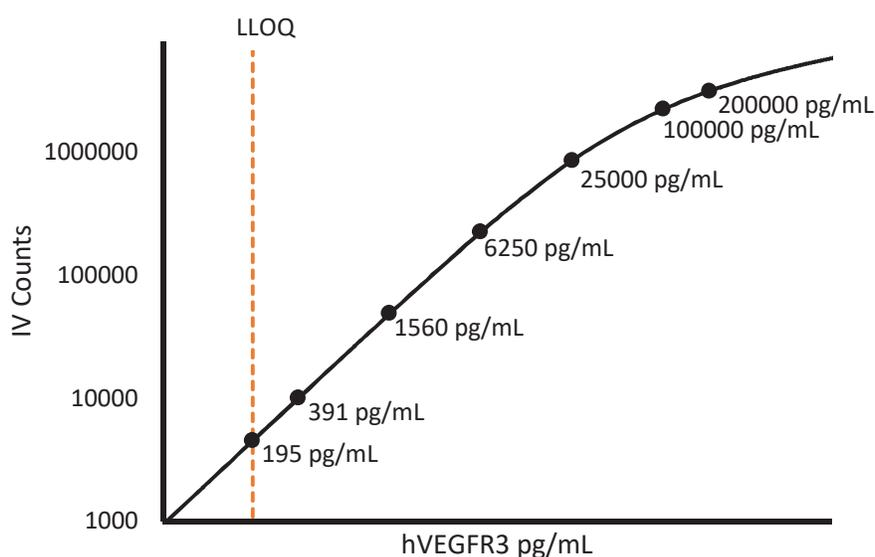


Description – Vascular Endothelial Growth Factor Receptor 3 (VEGF-R3)

VEGF-R3 is one of the three main subtypes of VEGFR. The ligands for VEGF-R3 are VEGF-C, and VEGF-D. Vascular endothelial growth factor (VEGF) is a 27KDa signaling protein produced by cells that stimulates vasculogenesis and angiogenesis. VEGF is a heparin binding protein and exists as a disulfide linked homo-dimer. It has at least 6 isoforms produced by alternative splicing. VEGF induces endothelial cell proliferation, promotes cell migration, inhibits apoptosis, and induces permeabilization of blood vessels.

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 (1:50 Dilution)*	50 μ 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	195 pg/mL
Functional LLOQ (x MRD)	9750 pg/mL
LOD	48.3 pg/mL
Assay Range	0 – 10 μg/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.