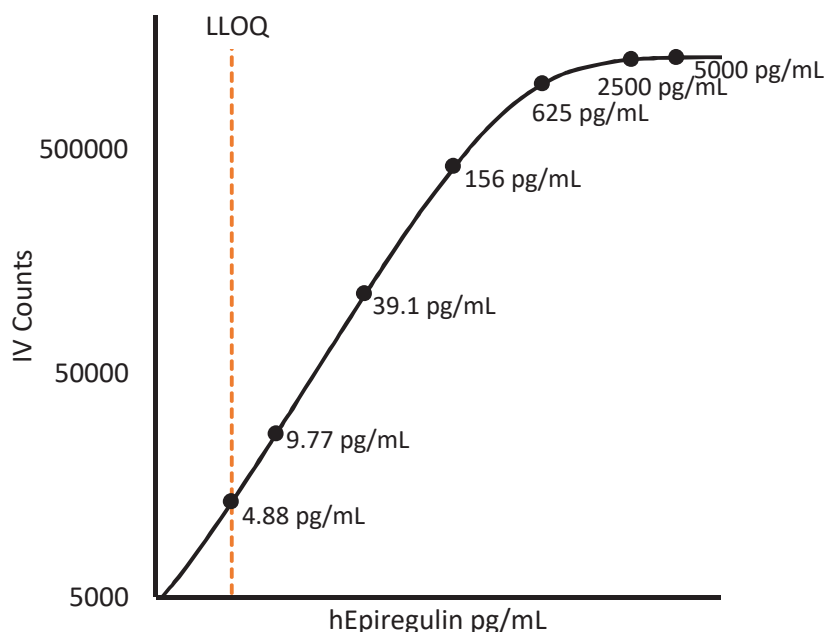


### Description – Epiregulin (ER)

Epiregulin is a 5.4 kDa protein and member of the epidermal growth factor (EGF) family. Epiregulin binds to ErbB proteins to regulate signaling in development, including the development of tumors. Epiregulin regulates inflammation, wound healing, tissue repair, and oocyte maturation through stimulus of angiogenesis, vascular remodeling, and cell proliferation.

**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=4) and serum (n=4) samples were measured.

% Above LOD	100%
% Above LLOQ	37.5%

**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	82.1 fg/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.