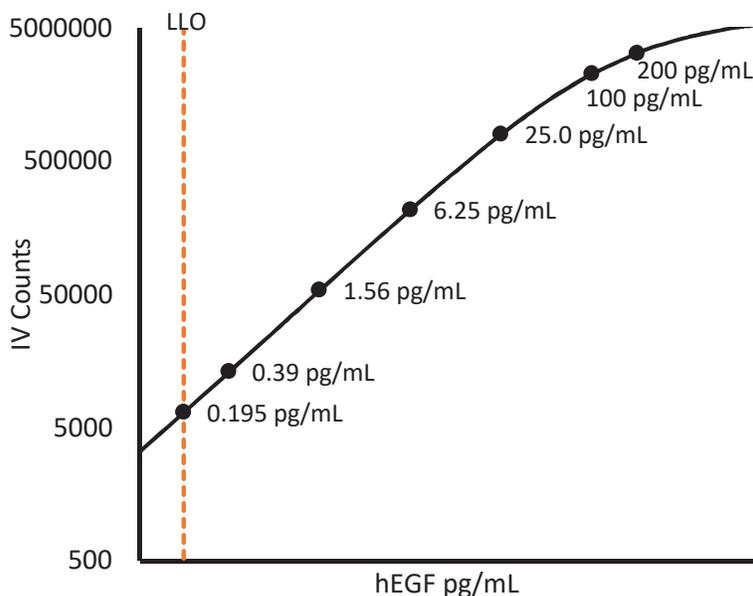


### Description – EGF

Epidermal Growth Factor (EGF) is a 6 kDa protein secreted primarily by the submaxillary gland. EGF consists of three intramolecular disulfide bonds that are essential in binding to other protein members of the EGF family. EGF binds to and activates the epidermal growth factor receptor to regulate cell growth and survival.

**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)

<b>Diluted Sample volume (1:2 Dilution)*</b>	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.

<b>% Above LOD</b>	<b>100%</b>
<b>% Above LLOQ</b>	<b>100%</b>

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

<b>Analytical LLOQ</b>	<b>0.195 pg/mL</b>
<b>Functional LLOQ (x MRD)</b>	<b>0.390 pg/mL</b>
<b>LOD</b>	<b>45.8 fg/mL</b>
<b>Assay Range</b>	<b>0 - 400 pg/mL</b>

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.