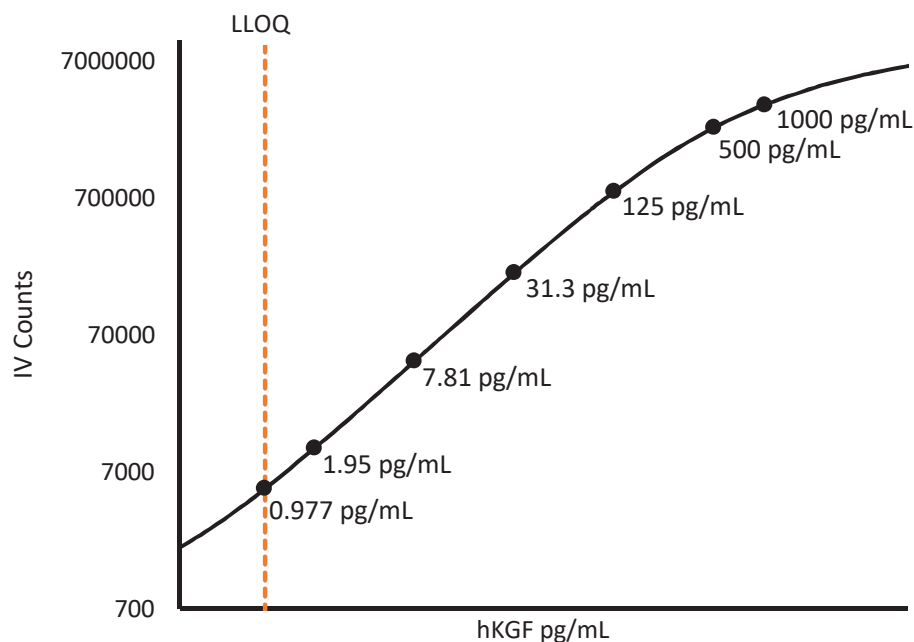


Description – Keratinocyte Growth Factor (KGF)

Keratinocyte growth factor (KGF) is a 19 kDA protein encoded by the fibroblast growth factor (FGF-7) gene. When secreted, KGF acts as a paracrine growth factor for nearby epithelial cells. KGF plays a key role in development, morphogenesis, angiogenesis, wound healing, and tumorigenesis.

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 | 50% |

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

| | |
|-------------------------|----------------|
| Analytical LLOQ | 0.977 pg/mL |
| Functional LLOQ (x MRD) | 1.95 pg/mL |
| LOD | 0.00287 pg/mL |
| Assay Range | 0 – 2000 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.