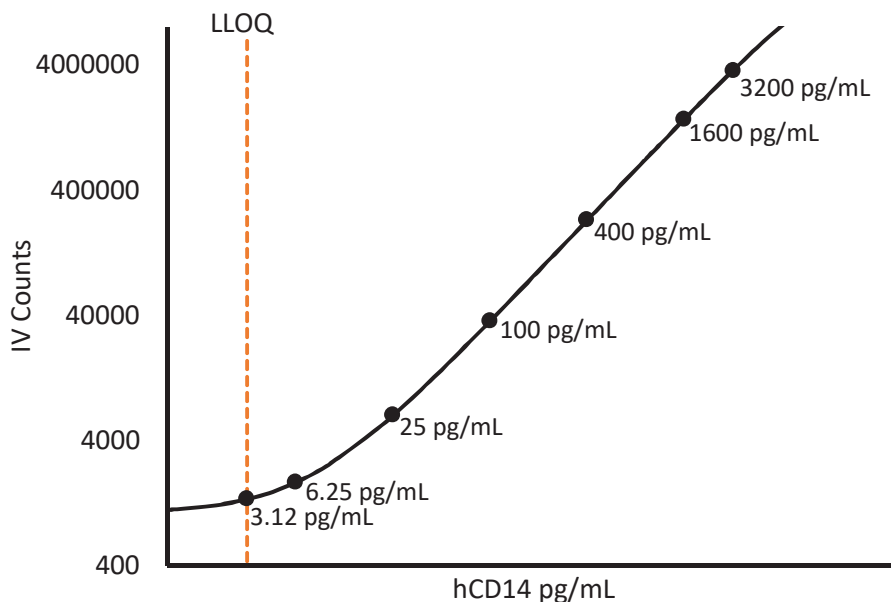


### Description – CD-14

CD-14 is a 55 kDa cell surface glycoprotein expressed on monocytes and most tissue macrophages. CD-14 is a GPI-anchored protein that functions as a pattern recognition receptor, binding lipopolysaccharides and many different ligands from a variety of microbial sources. When soluble, CD-14 functions as a monocyte activator.

**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 (1:5000 Dilution)*	50 $\mu$ L per measurement
---	-------------------------------

\*See Kit Instructions for details

**Endogenous Serum and Plasma Readings:** Healthy EDTA plasma (n=5) and serum (n=5) 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	3.13 pg/mL
Functional LLOQ (x MRD)	15.7 ng/mL
LOD	0.258 pg/mL
Assay Range	0 – 16 $\mu$ g/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.