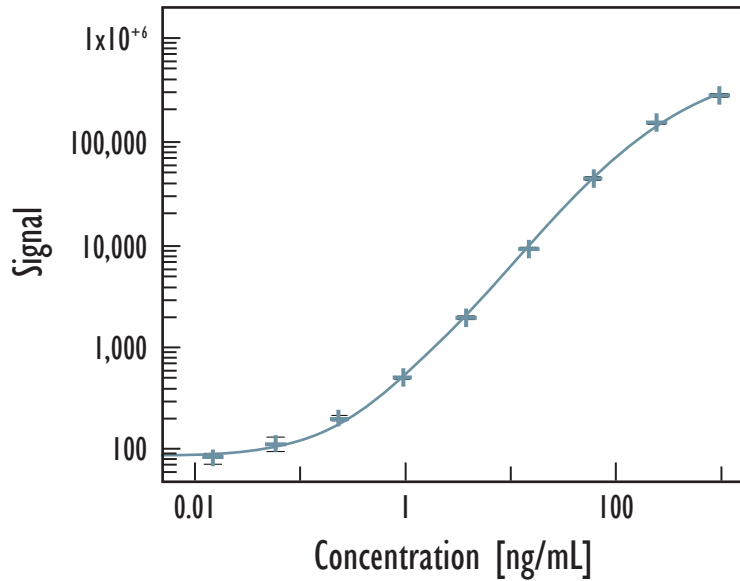


# MULTI-ARRAY<sup>®</sup> VCAM-1 Assay

## Detection of VCAM-1 in Serum and Plasma Samples



Standard curve data is from a representative experiment

Detection Limits (ng/mL)	VCAM-1
Average LLOD	0.07
Average LLOQ	0.7

LLOD is defined as 2.5x stdev above the background

LLOQ is designated by the lowest concentration on the standard curve where % CV is less than 20% and recovery of predicted concentration is within 20% of 100%

Kit Size	Catalog Number
1 plate	K151EQC-1
5 plates	K151EQC-2
20 plates	K151EQC-3
20 plates (Base)	K151EQA-3

### Endogenous Levels in Human Samples

- 16 normal human donors; matched sera and plasmas
- Detected level was above LLOQ in all samples
- Average CVs for measured samples was less than 10%
- Samples diluted 1:200 for use in assay

		ng/mL
Serum	Mean	198
	Median	160
	Range	118 - 444
EDTA Plasma	Mean	173
	Median	147
	Range	72 - 326
Heparin Plasma	Mean	159
	Median	141
	Range	83 - 349

# MULTI-ARRAY<sup>®</sup> VCAM-1 Assay

## Detection of VCAM-1 in Serum and Plasma Samples

### Dilutional Linearity

- Measured endogenous analyte levels in samples diluted 1:200 in assay diluent followed by subsequent dilution

$$\% \text{ recovery} = \frac{(\text{measured value} * \text{dilution factor} * 100)}{\text{predicted value}}$$

	Dilution Factor	% Recovery
Serum	1/2	112
	1/4	96
	1/8	93
EDTA Plasma	1/2	115
	1/4	102
	1/8	103
Heparin Plasma	1/2	103
	1/4	97
	1/8	94

Values presented are averages across two pooled samples

### Spike Recovery

- Measured analyte spiked into 1:200 diluted human samples

$$\% \text{ recovery} = \frac{(\text{measured value} * 100)}{\text{expected value}}$$

	% Recovery
Serum	101
EDTA Plasma	101
Heparin Plasma	100

Values presented are averages across two pooled samples and including spike levels of 12.5, 25, and 50 ng/mL