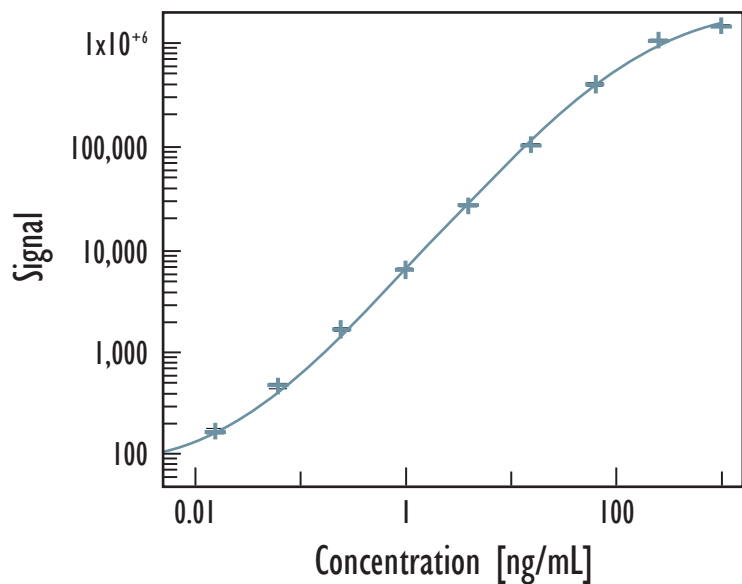


# MULTI-ARRAY<sup>®</sup> ICAM-1 Assay

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



Standard curve data is from a representative experiment

Detection Limits (ng/mL)	ICAM-1
Average LLOD	0.01
Average LLOQ	0.09

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	K151ERC-1
5 plates	K151ERC-2
20 plates	K151ERC-3
20 plates (Base)	K151ERA-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	152
	Median	107
	Range	71 - 294
EDTA Plasma	Mean	140
	Median	102
	Range	51 - 301
Heparin Plasma	Mean	122
	Median	91
	Range	43 - 256

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

## Detection of ICAM-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	100
	1/4	96
	1/8	99
EDTA Plasma	1/2	101
	1/4	97
	1/8	96
Heparin Plasma	1/2	98
	1/4	100
	1/8	98

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	93
EDTA Plasma	94
Heparin Plasma	91

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