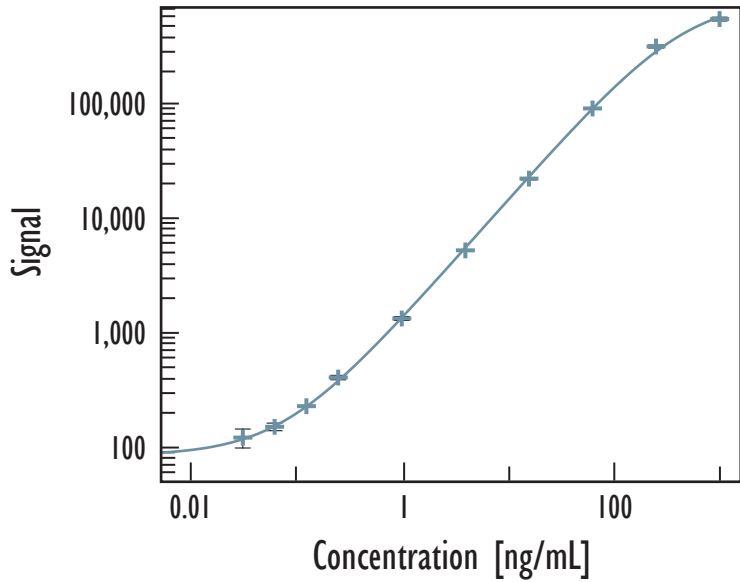


MULTI-ARRAY[®] ICAM-3 Assay

Detection of ICAM-3 in Serum and Plasma Samples



Standard curve data is from a representative experiment

Detection Limits (ng/mL)	ICAM-3
Average LLOD	0.05
Average LLOQ	0.6

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	K151ELC-1
5 plates	K151ELC-2
20 plates	K151ELC-3
20 plates (Base)	K151ELA-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% for all assays

		ng/mL
Serum	Mean	8.8
	Median	7.8
	Range	4.3 - 17.8
EDTA Plasma	Mean	6.6
	Median	5.6
	Range	2.8 - 14.8
Heparin Plasma	Mean	6.4
	Median	6.0
	Range	2.9 - 9.4

MULTI-ARRAY[®] ICAM-3 Assay

Detection of ICAM-3 in Serum and Plasma Samples

Dilutional Linearity

- Measured endogenous analyte levels in samples diluted into assay diluent

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

	Dilution Factor	% Recovery
Serum	1/2	98
	1/4	85
	1/8	101
EDTA Plasma	1/2	116
	1/4	95
	1/8	96
Heparin Plasma	1/2	108
	1/4	95
	1/8	105

Values presented are averages across two pooled samples

Spike Recovery

- Measured analyte spiked into undiluted human samples

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

	% Recovery
Serum	77
EDTA Plasma	71
Heparin Plasma	76

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