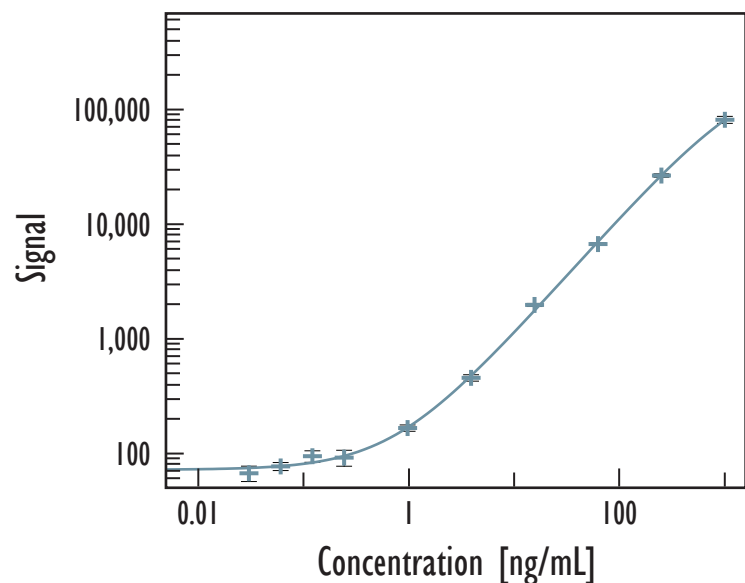


MULTI-ARRAY[®] E-Selectin Assay

Detection of E-Selectin in Serum and Plasma Samples



Standard curve data is from a representative experiment

Detection Limits (ng/mL)	E-Selectin
Average LLOD	0.2
Average LLOQ	0.8

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	K151EMC-1
5 plates	K151EMC-2
20 plates	K151EMC-3
20 plates (Base)	K151EMA-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	15.8
	Median	17.6
	Range	6.5 - 23.4
EDTA Plasma	Mean	8.3
	Median	9.1
	Range	2.9 - 13.1
Heparin Plasma	Mean	11.9
	Median	12.1
	Range	3.4 - 20.3

MULTI-ARRAY[®] E-Selectin Assay

Detection of E-Selectin 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	114
	1/4	95
	1/8	106
EDTA Plasma	1/2	143
	1/4	118
	1/8	107
Heparin Plasma	1/2	127
	1/4	123
	1/8	101

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	97
EDTA Plasma	69
Heparin Plasma	91

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