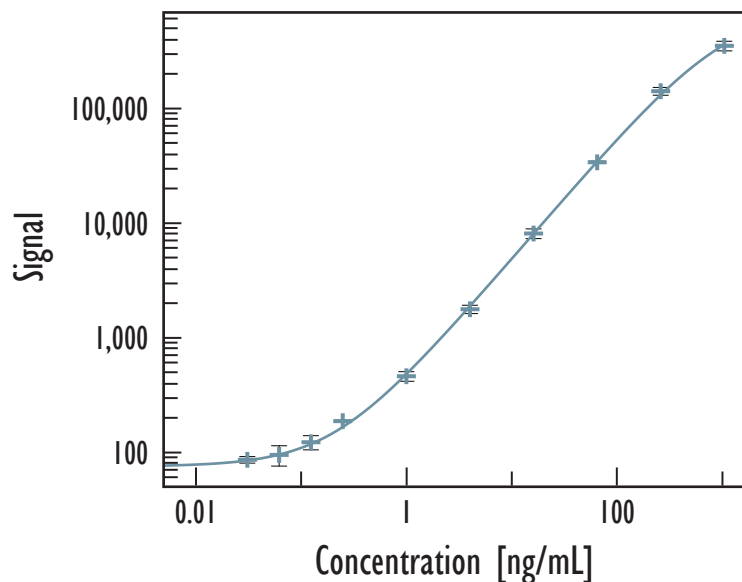


# MULTI-ARRAY<sup>®</sup> P-Selectin Assay

## Detection of P-Selectin in Serum and Plasma Samples



Standard curve data is from a representative experiment

Detection Limits (ng/mL)	P-Selectin
Average LLOD	0.1
Average LLOQ	1.0

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	K151ENC-1
5 plates	K151ENC-2
20 plates	K151ENC-3
20 plates (Base)	K151ENA-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	40.2
	Median	39.7
	Range	21.7 - 58.3
EDTA Plasma	Mean	50.2
	Median	44.2
	Range	28.5 - 102.8
Heparin Plasma	Mean	62.3
	Median	55.7
	Range	31.1 - 124

# MULTI-ARRAY<sup>®</sup> P-Selectin Assay

## Detection of P-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	95
	1/4	89
	1/8	111
EDTA Plasma	1/2	115
	1/4	110
	1/8	116
Heparin Plasma	1/2	113
	1/4	108
	1/8	115

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	89
EDTA Plasma	85
Heparin Plasma	91

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