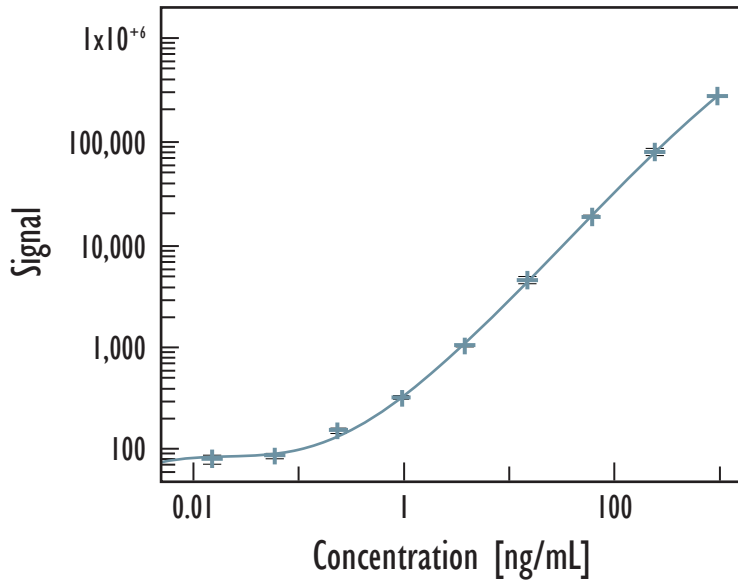


MULTI-ARRAY[®] SAA Assay

Detection of SAA in Serum and Plasma Samples



Standard curve data is from a representative experiment

Detection Limits (ng/mL)	SAA
Average LLOD	0.09
Average LLOQ	0.5

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	K151EOC-1
5 plates	K151EOC-2
20 plates	K151EOC-3
20 plates (Base)	K151EOA-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	33,086
	Median	11,001
	Range	3,899 - 147,438
EDTA Plasma	Mean	30,105
	Median	9,286
	Range	3,085 - 132,933
Heparin Plasma	Mean	26,114
	Median	9,057
	Range	2,586 - 106,487

MULTI-ARRAY[®] SAA Assay

Detection of SAA 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	97
	1/4	97
	1/8	97
EDTA Plasma	1/2	107
	1/4	99
	1/8	101
Heparin Plasma	1/2	95
	1/4	100
	1/8	92

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	96
Heparin Plasma	85

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