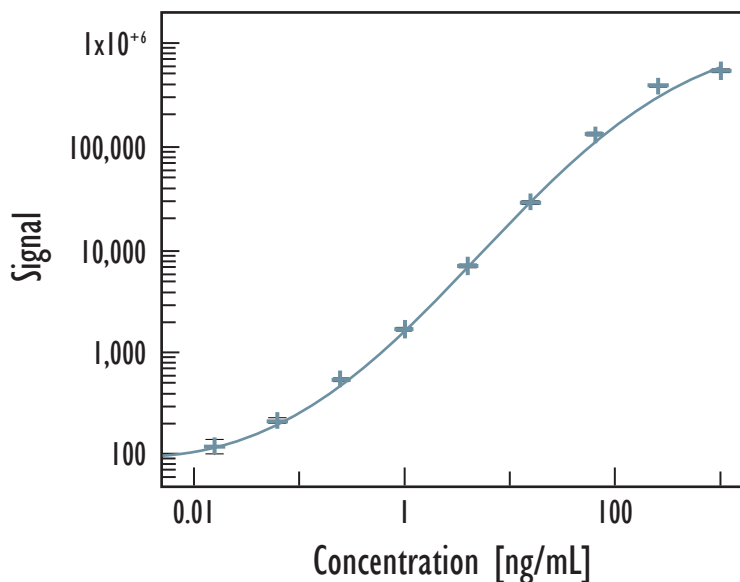


MULTI-ARRAY[®] CRP Assay

Detection of CRP in Serum and Plasma Samples



Standard curve data is from a representative experiment

Detection Limits (ng/mL)	CRP
Average LLOD	0.1
Average LLOQ	0.7

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	SI2400	SI6000
1 plate	K151EPC-1	K111EPC-1
5 plates	K151EPC-2	K111EPC-2
20 plates	K151EPC-3	K111EPC-3
20 plates (Base)	K151EPA-3	K111EPA-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
- Samples diluted 1:200 for use in assay

		ng/mL
Serum	Mean	11,036
	Median	7,007
	Range	756 - 56,143
EDTA Plasma	Mean	7,065
	Median	2,394
	Range	623 - 58,543
Heparin Plasma	Mean	6,553
	Median	4,028
	Range	482 - 45,493

MULTI-ARRAY[®] CRP Assay

Detection of CRP 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	102
	1/4	99
	1/8	96
EDTA Plasma	1/2	99
	1/4	96
	1/8	101
Heparin Plasma	1/2	100
	1/4	105
	1/8	96

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	103
EDTA Plasma	102
Heparin Plasma	93

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