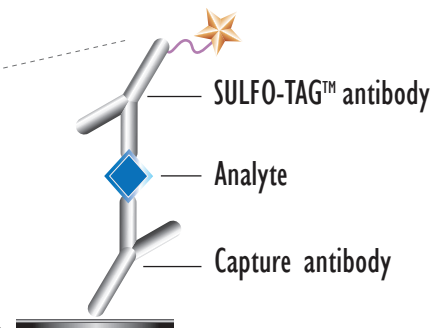
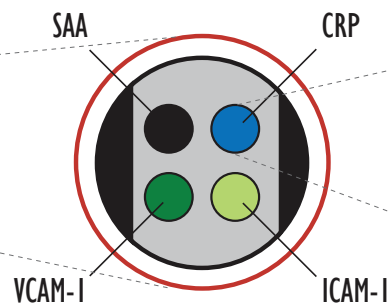
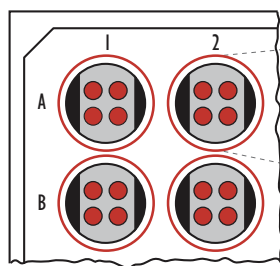


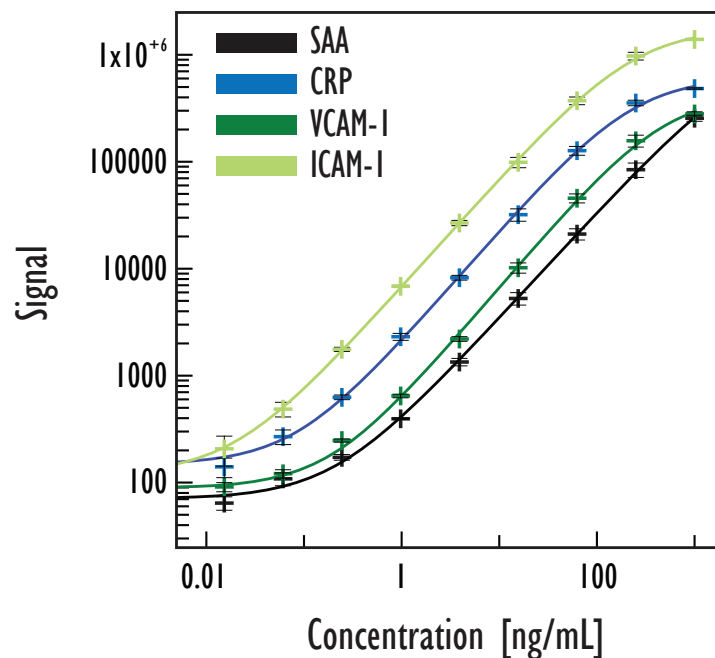
# MULTI-SPOT® Vascular Injury II Assay

## Detection of Vascular Biomarkers in Human Serum Samples

MSD MULTI-SPOT®  
96-Well 4-Spot Plate



## Assay Performance



Detection Limits (ng/mL)	SAA	CRP	VCAM-I	ICAM-I
Average LLOD	0.09	0.1	0.07	0.01
Average LLOQ	0.5	0.7	0.7	0.09

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%

Standard curve data is from a representative experiment

Kit Size	Catalog Number
1 plate	K15136C-1
5 plates	K15136C-2
20 plates	K15136C-3
20 plates (Base)	K15136A-3

# MULTI-SPOT® Vascular Injury II Assay

## Endogenous Levels in Human Samples

- 16 normal human donors; matched sera and plasmas
- Detected level was above LLOQ for all analytes 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	SAA	CRP	VCAM-I	ICAM-I
Serum	Mean	33086	11036	198	152
	Median	11001	7007	160	107
	Range	3899 - 147438	756 - 56143	118 - 444	71 - 294
EDTA Plasma	Mean	30105	7065	173	140
	Median	9286	2394	147	102
	Range	3085 - 132933	623 - 58543	72 - 326	51 - 301
Heparin Plasma	Mean	26114	6553	159	122
	Median	9057	4028	141	91
	Range	2586 - 106487	482 - 45493	83 - 349	43 - 256

## Dilution 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	SAA	CRP	VCAM-I	ICAM-I
Serum	1/2	97	102	112	100
	1/4	97	99	96	96
	1/8	97	96	93	99
EDTA Plasma	1/2	107	99	115	101
	1/4	99	96	102	97
	1/8	101	101	103	96
Heparin Plasma	1/2	95	100	103	98
	1/4	100	105	97	100
	1/8	92	96	94	98

Values presented are averages across two pooled samples

## Spike Recovery

- Measured analyte spiked into 1:200 diluted human samples

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

	SAA	CRP	VCAM-I	ICAM-I
Serum	93	103	101	93
EDTA Plasma	96	102	101	94
Heparin Plasma	85	93	100	91

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