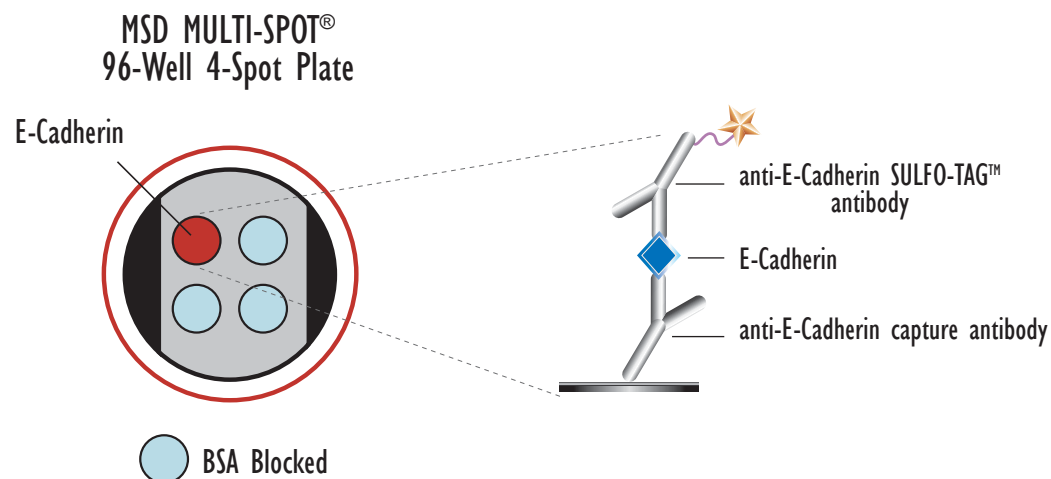
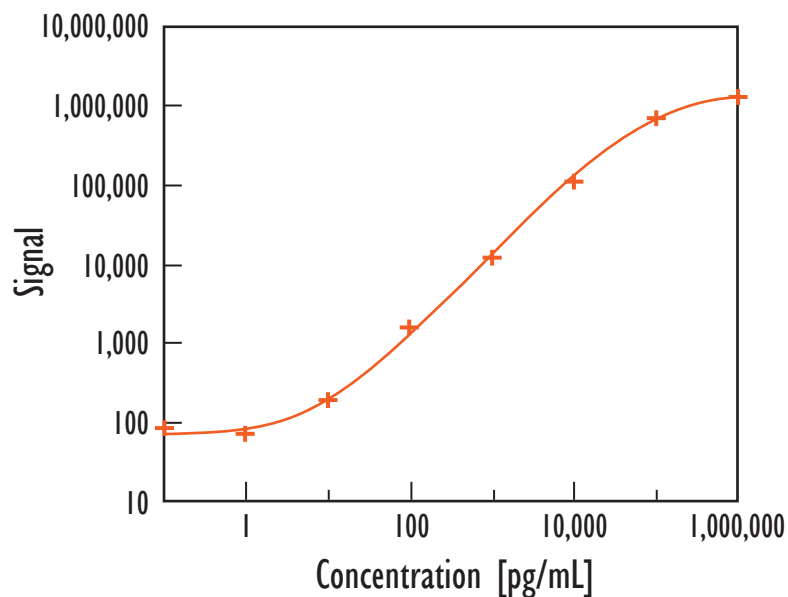


# MULTI-ARRAY<sup>®</sup> Human E-Cadherin Assay

Detection of E-Cadherin in Human Serum and Plasma Samples



Concentration (pg/mL)	Average	%CV
0	85	19
1	71	31
10	193	13
100	1,616	13
1,000	12,272	10
10,000	112,665	9
100,000	713,947	8
1,000,000	1,351,953	9

Standard curve data is from a representative experiment

1:10 dilution of serum and plasma samples is recommended for this assay

E-Cadherin LLOD	4 (pg/mL)
-----------------	-----------

LLOD (Lower Limit of Detection) is defined as 2.5x stdev above the background

Kit Size	SI2400	SI6000
1 plate	K1511ZC-1	K1111ZC-1
5 plates	K1511ZC-2	K1111ZC-2
20 plates	K1511ZC-3	K1111ZC-3
20 plates (Base)	K1511ZA-3	K1111ZA-3

# MULTI-ARRAY<sup>®</sup> Human E-Cadherin Assay

## Detection of E-Cadherin in Human Serum and Plasma Samples

### Dilutional Linearity

- Samples from 7 apparently healthy donors were diluted in Calibrator Diluent

$$\% \text{ recovery} = \frac{(\text{measured value} * \text{dilution factor} * 100)}{\text{predicted value}}$$

- 1X dilution refers to the dilution recommended for serum, i.e. a 10-fold dilution

Dilution Factor	Percent Recovery (%)
2X	119
0.5X	82
0.25X	67

### Endogenous Levels in Human Samples

- 95 normal human donors, Serum
- Average CVs for measured samples was less than 10%

N (ng/mL)	Mean (ng/mL)	Median (ng/mL)	Range (ng/mL)
95	53	50	20 - 106

### Spike Recovery

- Measured analyte spiked into apparently normal human samples

$$\% \text{ recovery} = \frac{(\text{measured spiked value} - \text{measured native})}{\text{spike}}$$

Sample	Neat (ng/mL)	Spiked (ng/mL)	Percent Recovery (%)
S1	14	83	71
S2	8	69	61
S3	5	103	98
S4	6	69	63
S5	17	91	75
S6	14	93	80
S7	14	92	80
Average Percent Recovery (%)			76