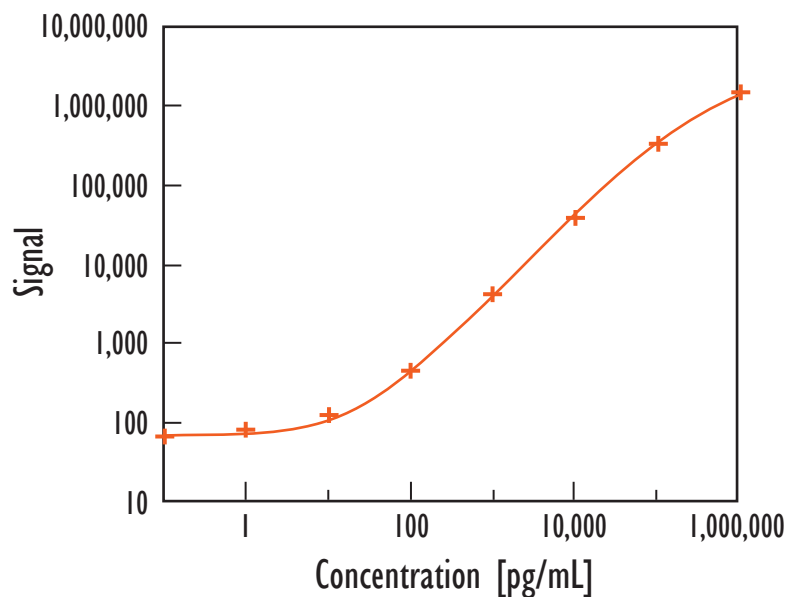


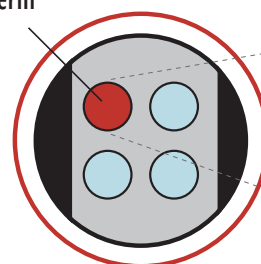
# MULTI-ARRAY<sup>®</sup> Human P-Cadherin Assay

Detection of P-Cadherin in Human Serum and Plasma Samples

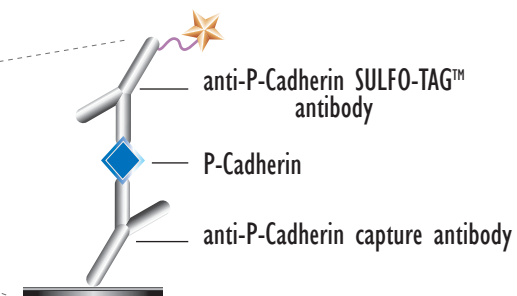


MSD MULTI-SPOT<sup>®</sup>  
96-Well 4-Spot Plate

P-Cadherin



BSA Blocked



Concentration (pg/mL)	Average	%CV
0	63	68
1	80	56
10	122	7
100	436	21
1,000	4,218	3
10,000	39,393	10
100,000	329,464	9
1,000,000	1,528,596	11

Standard curve data is from a representative experiment

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

P-Cadherin LLOD	13 (pg/mL)
-----------------	------------

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

Kit Size	SI2400	SI6000
1 plate	K151JAC-1	K111JAC-1
5 plates	K151JAC-2	K111JAC-2
20 plates	K151JAC-3	K111JAC-3
20 plates (Base)	K151JAA-3	K111JAA-3

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

## Detection of P-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	116
0.5X	102
0.25X	91

### 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	38	37	17 - 75

### 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	11	90	80
S2	43	120	81
S3	33	129	99
S4	21	89	70
S5	38	132	98
S6	40	127	91
S7	33	119	89
Average Percent Recovery (%)			87