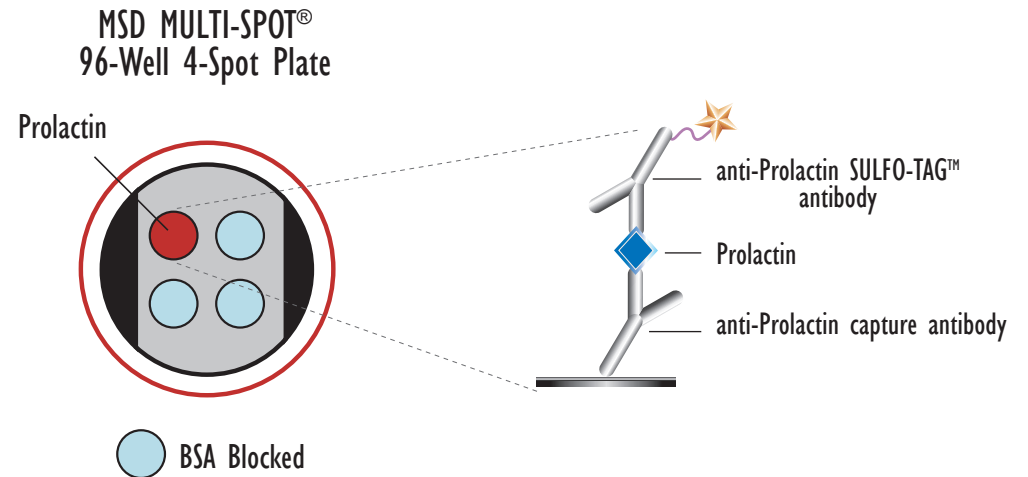
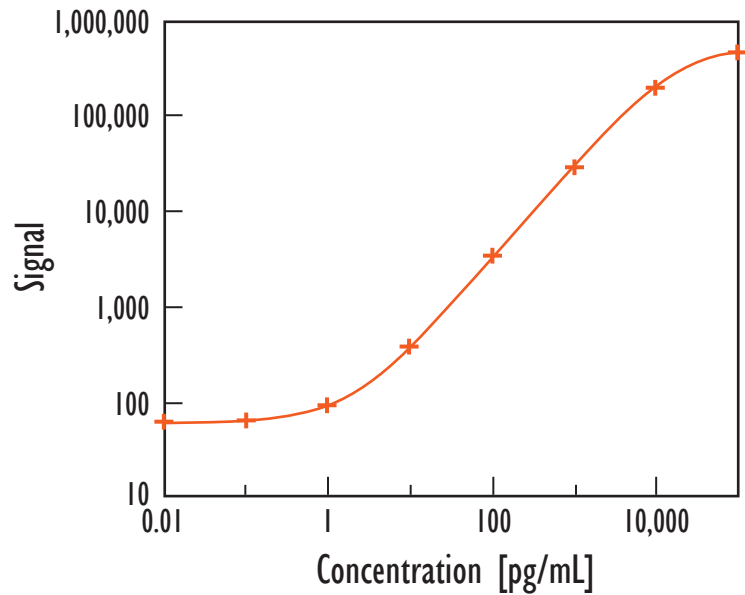


MULTI-ARRAY[®] Human Prolactin Assay

Detection of Prolactin in Human Serum and Plasma Samples



Concentration (pg/mL)	Average	%CV
0	61	15
0.1	62	30
1	89	22
10	364	11
100	3,442	6
1,000	28,418	8
10,000	199,784	8
100,000	460,288	11

Standard curve data is from a representative experiment

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

Prolactin LLOD	2 (pg/mL)
----------------	-----------

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

Kit Size	Catalog Number
1 plate	K151JBC-1
5 plates	K151JBC-2
20 plates	K151JBC-3
20 plates (Base)	K151JBA-3

MULTI-ARRAY[®] Human Prolactin Assay

Detection of Prolactin 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	102
0.5X	98
0.25X	96

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	3.7	2.9	0.7 - 21

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	2.8	12	91
S2	8.5	16	83
S3	5.8	14	87
S4	6.4	16	98
S5	4.2	14	106
S6	4.3	16	123
S7	19	31	135
Average Percent Recovery (%)			103