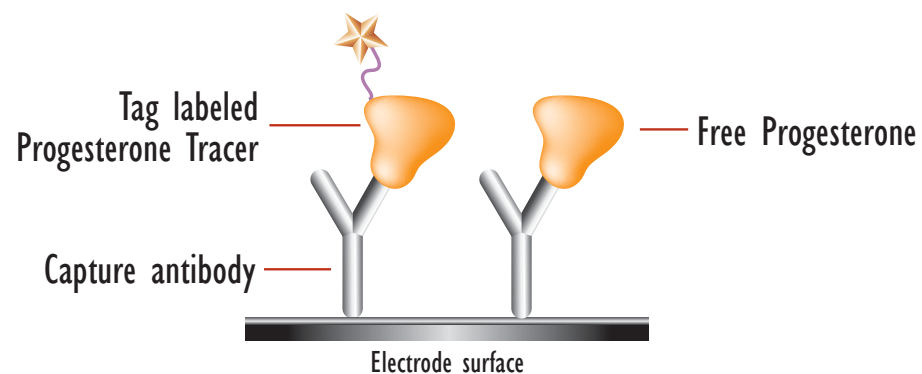
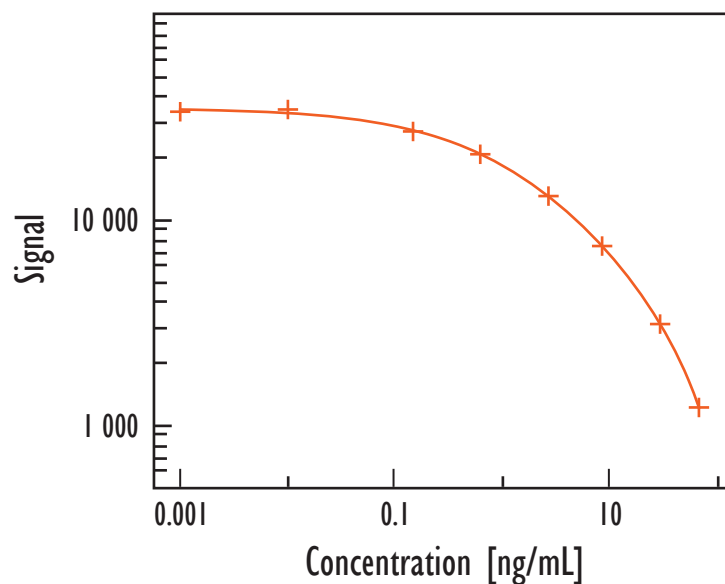


# MULTI-ARRAY<sup>®</sup> Human Progesterone Assay

*Detection of Progesterone in Human Serum and Plasma Samples*



Concentration (ng/mL)	Average	%CV
0	37925	4.7
0.01	37473	3.9
0.1	29310	2.6
1	22705	12.6
3	14212	1.3
8	8122	6.7
29	3389	3.9
66	1332	7.3

Standard curve data is from a representative experiment

**Progesterone LLOD 0.16 (ng/mL)**

LLOD (Lower Limit of Detection) is defined as 2.5X stdev below the zero calibrator signal

Kit Size	Catalog Number
1 plate	K151EUC-1
5 plates	K151EUC-2
20 plates	K151EUC-3
20 plates (Base)	K151EUA-3