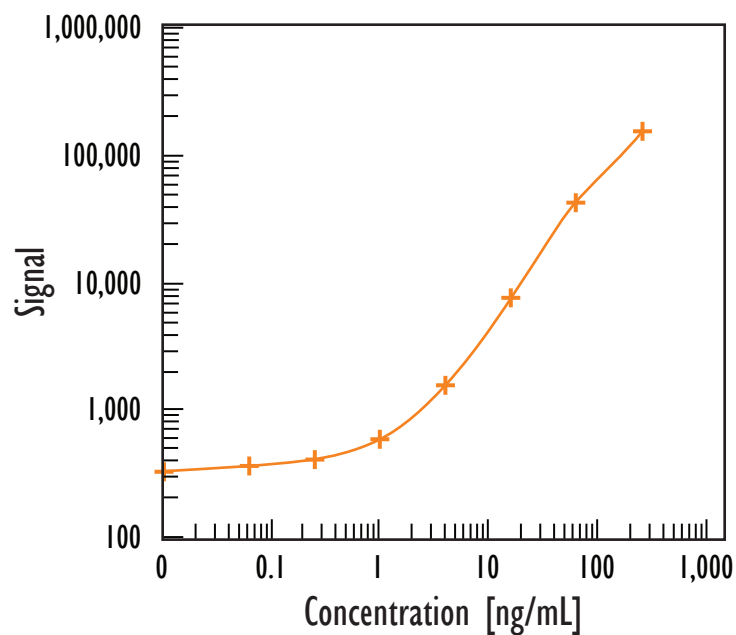
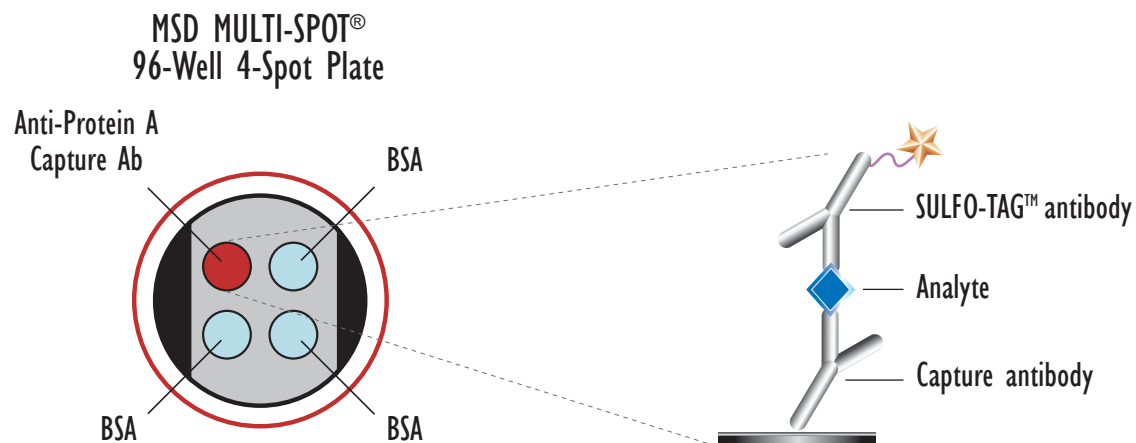


# MULTI-ARRAY<sup>®</sup> Protein A Contamination Assay

## Detection of Protein A



Concentration (ng/mL)	Protein A
0	335
0.061	375
0.24	424
1.0	614
3.9	1,616
16	7,775
63	44,595
250	159,081



## Detection Limits

	Protein A
LLOD (ng/mL)	0.4

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

Standard curve data is from a representative experiment

Kit Size	Catalog Number
1 plate	K150HMF-1
5 plates	K150HMF-2
20 plates	K150HMF-3
20 plates (Base)	K150HMA-3