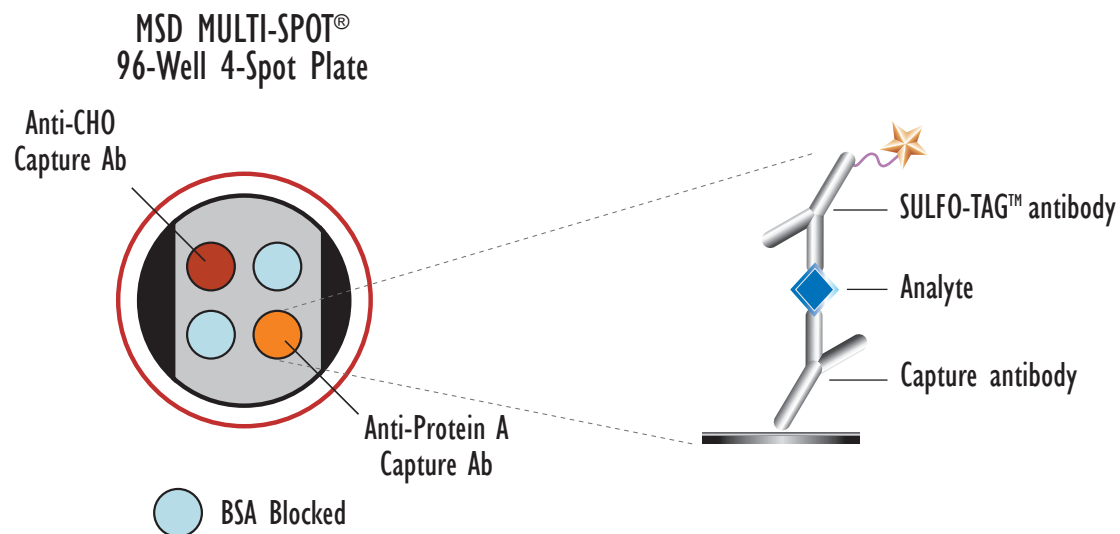
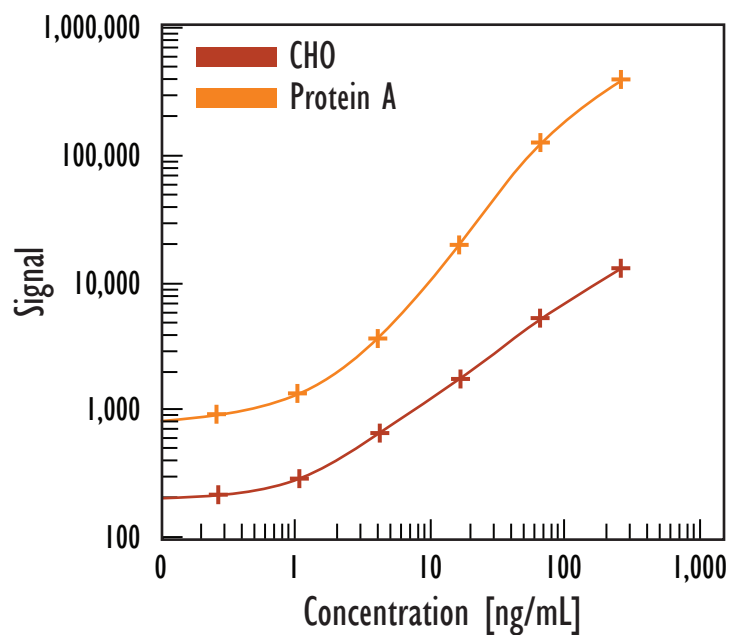


MULTI-SPOT[®] Bioprocess Panel I Assay

Detection of CHO Host Cell Protein and Protein A



Concentration (ng/mL)	CHO	Protein A
0	184	726
0.061	205	775
0.24	220	899
1.0	298	1,323
3.9	675	3,574
16	1,754	19,710
63	5,301	124,592
250	12,896	389,813
Avg % CV	6.8	3.6

Detection Limits

	CHO	Protein A
LLOD (ng/mL)	0.2	0.2

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	K15149F-1
5 plates	K15149F-2
20 plates	K15149F-3
20 plates (Base)	K15149A-3