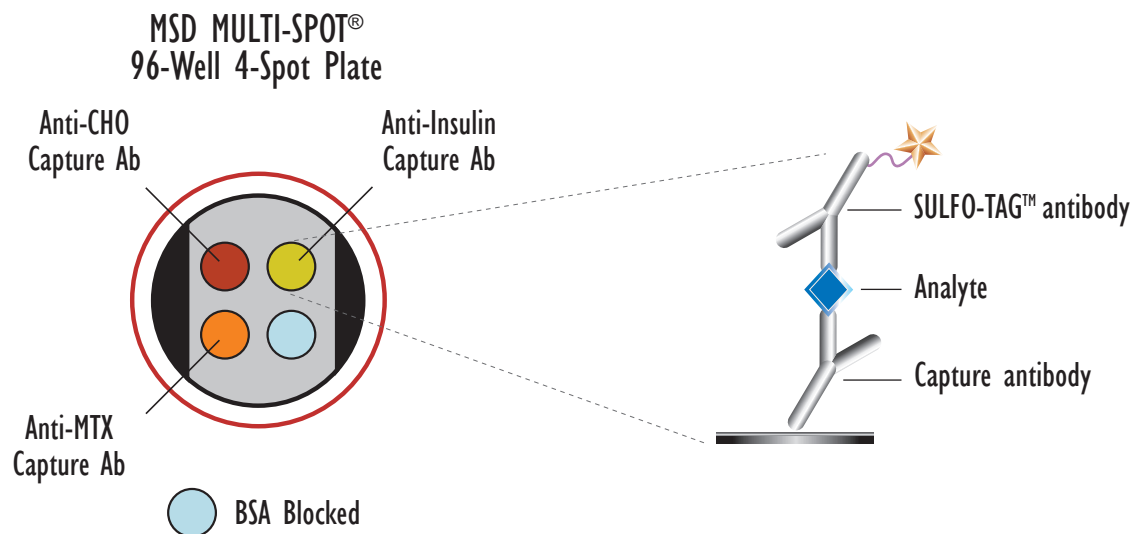
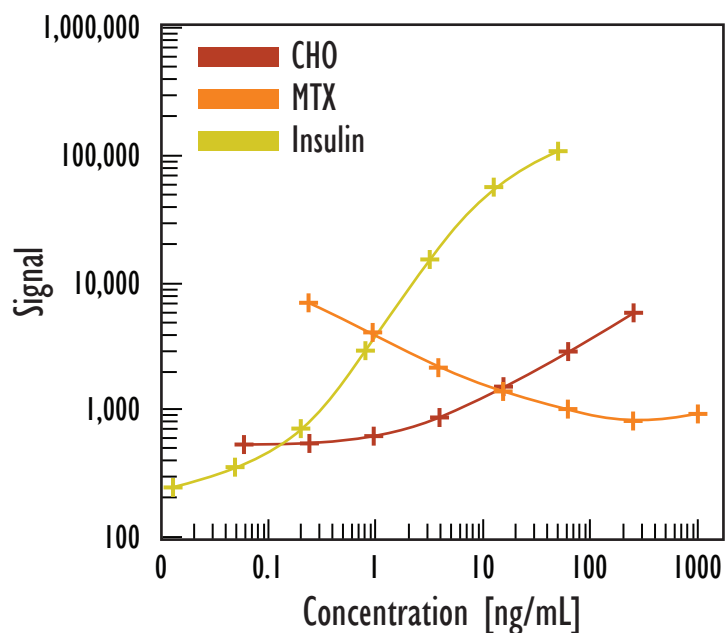


# MULTI-SPOT<sup>®</sup> Bioprocess Panel II Assay

Detection of CHO Host Cell Protein, Insulin, and Methotrexate



Concentration (ng/mL)	CHO
0	486
0.061	546
0.24	551
1.0	633
3.9	877
16	1,538
63	2,877
250	5,680
LLOD (ng/mL)	0.6

Concentration (ng/mL)	MTX
0	11,708
0.24	6,844
1.0	4,044
3.9	2,134
16	1,386
63	1,007
250	815
1000	924
LLOD (ng/mL)	0.06

Concentration (ng/mL)	Insulin
0	215
0.012	251
0.049	362
0.20	716
0.78	2,929
3.1	15,184
13	56,060
50	106,864
LLOD (ng/mL)	0.03

Kit Size	Catalog Number
1 plate	K15150F-1
5 plates	K15150F-2
20 plates	K15150F-3
20 plates (Base)	K15150A-3

Standard curve data is from a representative experiment

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