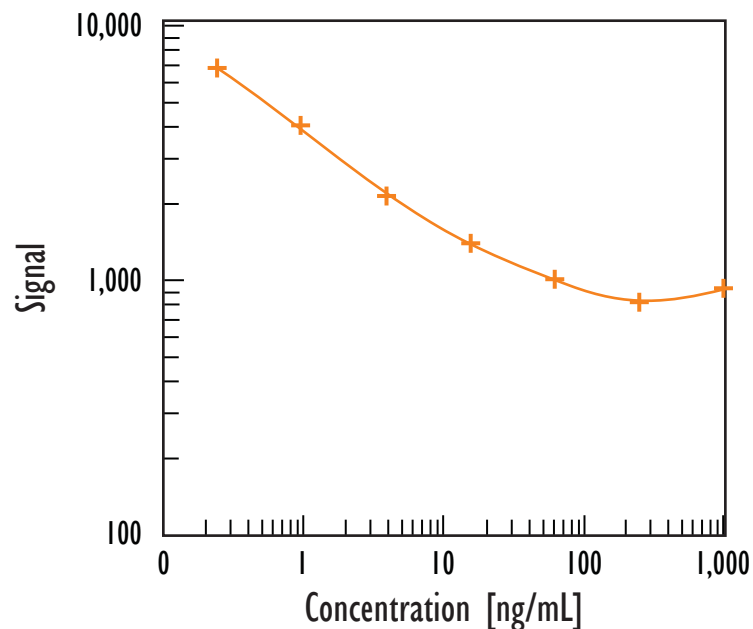


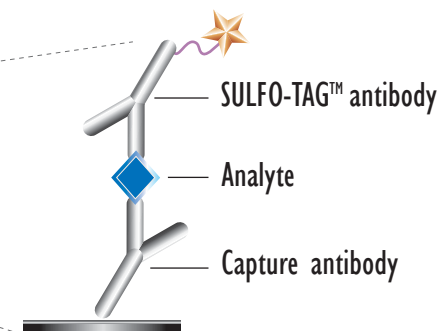
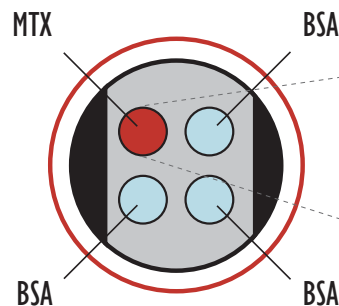
MULTI-ARRAY[®] Methotrexate Assay

Detection of Methotrexate (MTX) in Bioproduction Samples



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
1,000	924

MSD MULTI-SPOT[®]
96-Well 4-Spot Plate



Detection Limits

	MTX
LLOD (ng/mL)	0.06

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

Standard curve data is from a representative experiment

Kit Size	SI2400	SI6000
1 plate	K150HOF-1	K110HOF-1
5 plates	K150HOF-2	K110HOF-2
20 plates	K150HOF-3	K110HOF-3
20 plates (Base)	K150HOA-3	K110HOA-3