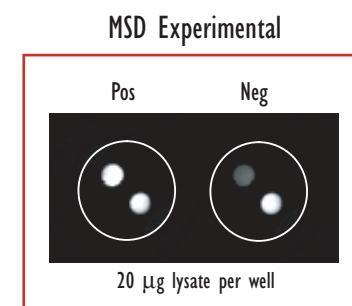
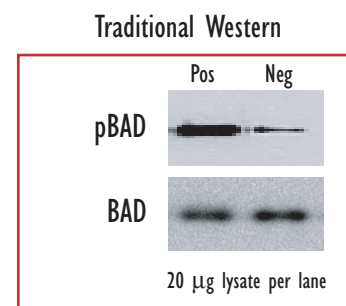
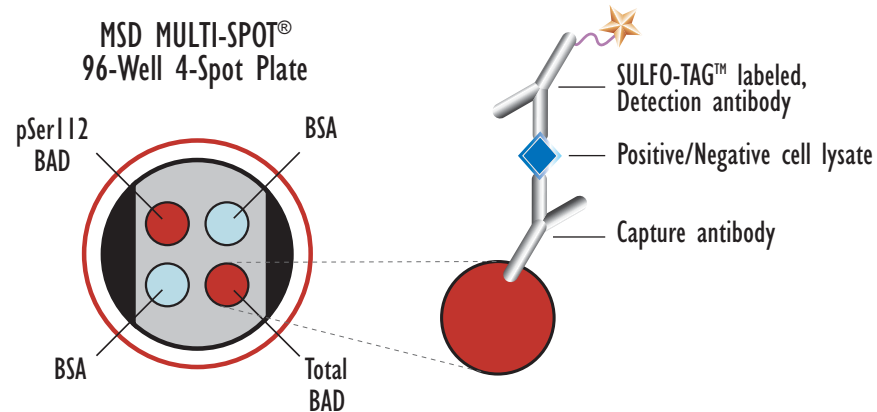
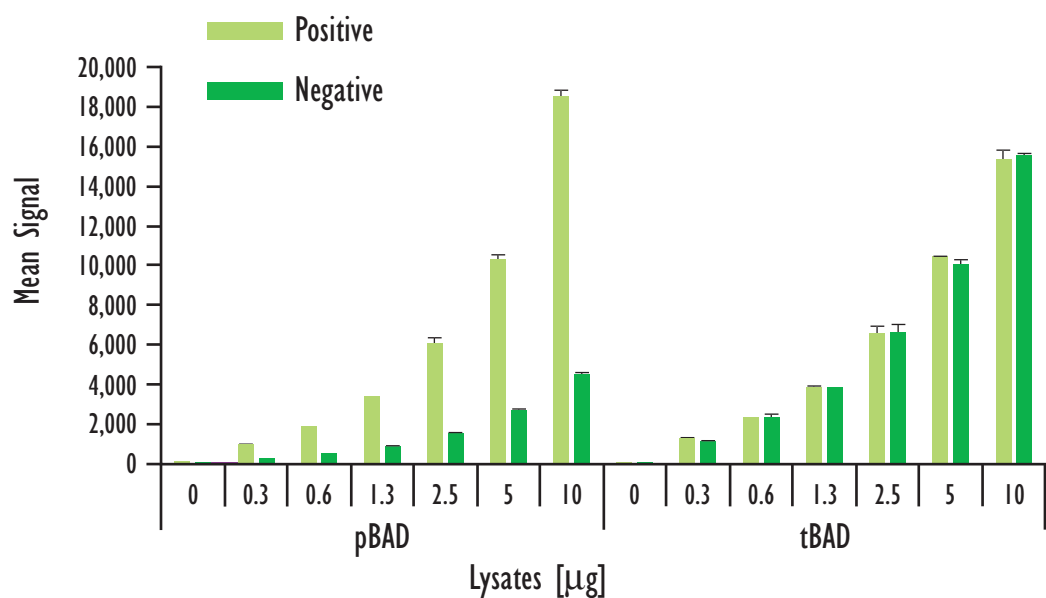


Multiplex BAD Assay: Detection of Phosphorylated and Total BAD in the Same Well



	Lysates (µg)	pBAD Positive			pBAD Negative			P/N
		Average	StdDev	%CV	Average	StdDev	%CV	
pBAD	0	44	13	30	41	9	22	
	0.3	937	45	5	265	5	2	3.5
	0.6	1,823	28	2	489	15	3	3.7
	1.3	3,314	57	2	870	9	1	3.8
	2.5	6,015	350	6	1,522	67	4	4
	5	10,311	254	2	2,658	68	3	3.9
	10	18,483	336	2	4,491	113	3	4.1
tBAD	0	36	13	35	37	13	35	
	0.3	1,222	93	8	1,112	26	2	1.1
	0.6	2,323	28	1	2,316	164	7	1
	1.3	3,840	83	2	3,820	48	1	1
	2.5	6,551	369	6	6,626	365	6	1
	5	10,375	99	1	10,015	273	3	1
	10	15,311	497	3	15,522	120	1	1

Serum deprived COS-7 cells (negative) were treated with PMA (200 nM; 1 hr)(positive). Whole cell lysates were added to MSD MULTI-SPOT 4-Spot plates coated with anti-phospho-BAD antibody and anti-total-BAD antibody on two of the four spatially distinct electrodes per well. Phosphorylated and total BAD were detected with anti-total-BAD antibody labeled with MSD SULFO-TAG reagent.