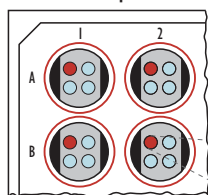
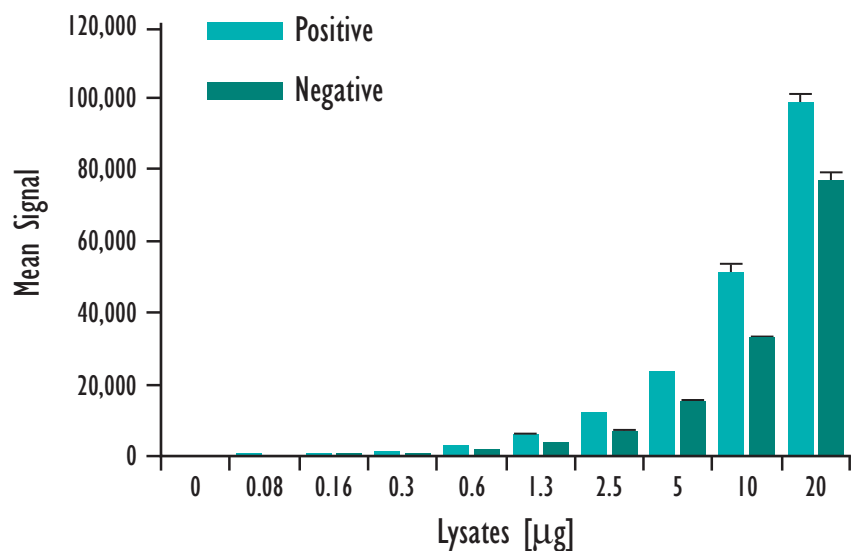
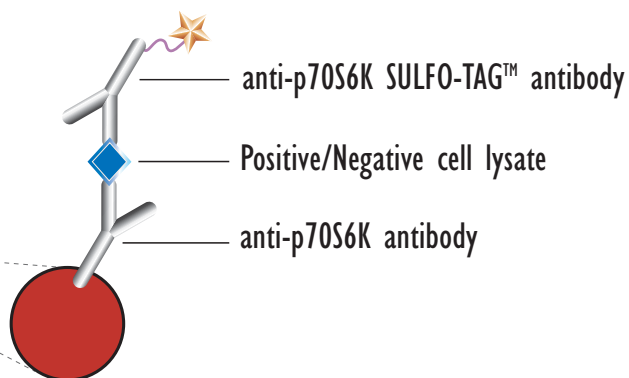


Detection of Total p70S6K

MSD MULTI-SPOT®
96-Well 4-Spot Plate

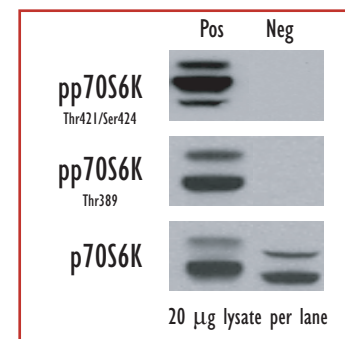


BSA

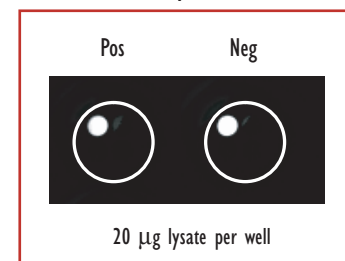


Lysates (µg)	pp70S6K Positive			pp70S6K Negative			P/N
	Average	StdDev	%CV	Average	StdDev	%CV	
0				134	9	7	
0.08	304	11	4	255	4	2	1.2
0.16	532	10	2	382	12	3	1.4
0.3	1,042	6	1	682	38	6	1.5
0.6	2,527	33	1	1,459	25	2	1.7
1.3	5,779	205	4	3,386	58	2	1.7
2.5	11,725	284	2	6,990	146	2	1.7
5	23,393	96	0	15,017	745	5	1.6
10	50,678	2,826	6	32,738	728	2	1.5
20	98,081	2,865	3	76,592	2,439	3	1.3

Traditional Western



MSD Experimental



Growing HEK293 cells were treated with rapamycin (1 µM; 3 hr) (negative) or Calyculin A (50 nM, 30 min)(positive). Whole cell lysates were added to MSD MULTI-SPOT 4-Spot plates coated with anti-total-p70S6K antibody on one of the four spatially distinct electrodes per well. Total p70S6K was detected with anti-total-p70S6K antibody labeled with MSD SULFO-TAG reagent.



Meso Scale Discovery

A division of Meso Scale Diagnostics, LLC.
www.mesoscale.com