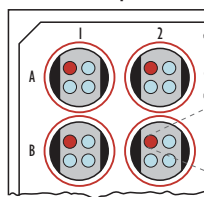


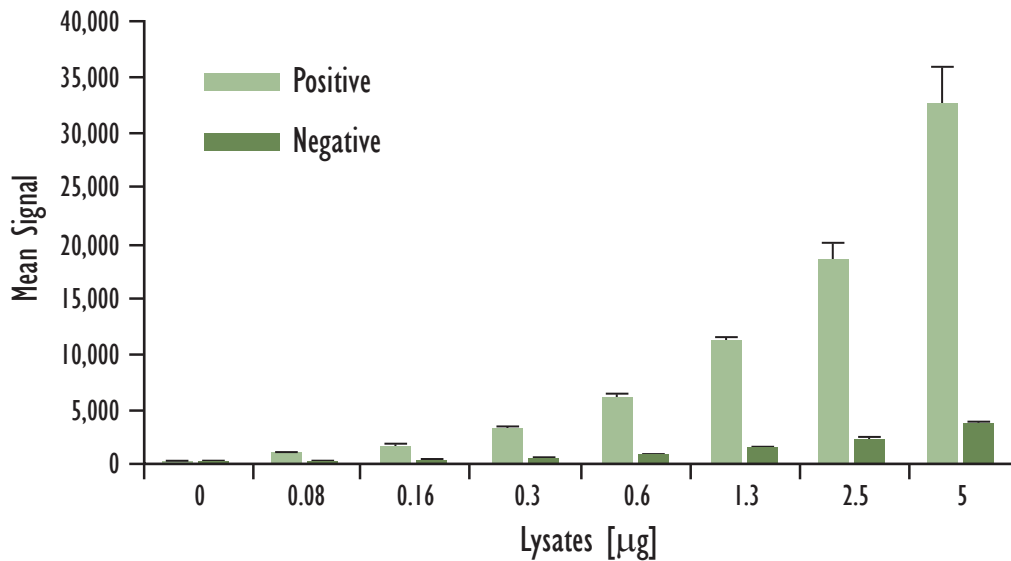
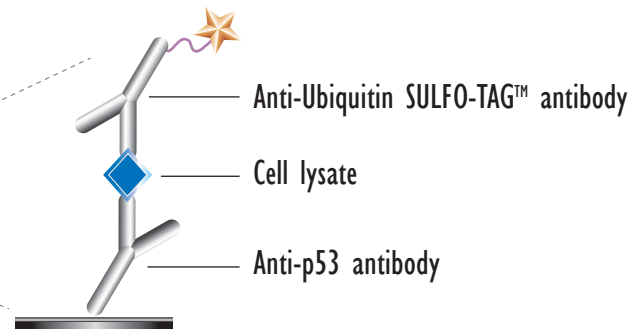
# MULTI-ARRAY<sup>®</sup> Ubiquitinated p53 Assay

Detection of Directly Ubiquitinated p53 & p53 in Complexes with Other Ubiquitinated Proteins in Whole Cell Lysate Samples

MSD MULTI-SPOT<sup>®</sup>  
96-Well 4-Spot Plate

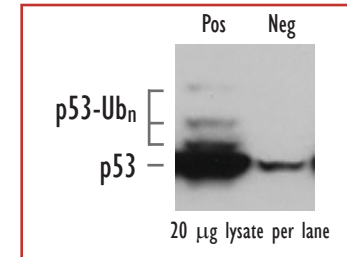


● BSA

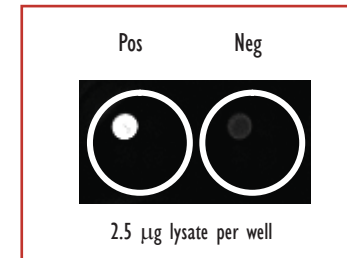


Lysates (µg)	Positive			Negative			P/N
	Average	StdDev	%CV	Average	StdDev	%CV	
0	245	8	3	243	14	6	1.0
0.08	1,018	10	1	321	11	3	3.2
0.16	1,754	50	3	420	11	3	4.2
0.3	3,327	138	4	554	53	10	6.0
0.6	6,006	312	5	858	51	6	7.0
1.3	11,281	277	2	1,482	57	4	7.6
2.5	18,596	1522	8	2,362	173	7	7.9
5	32,744	3009	9	3,714	141	4	8.8

Anti-p53 Western Blot



MSD Experimental



Growing HCT116 cells (negative) were treated with doxorubicin (1 µM, 21 hours) and epoxomicin (1 µM, 6 hours)(positive). Whole cell lysates were added to MSD MULTI-SPOT 4-Spot plates coated with anti-p53 antibody on one of the four spatially distinct electrodes per well. Ubiquitinated p53 was detected with antibody against ubiquitinated proteins labeled with MSD SULFO-TAG reagent.

Kit Size	Catalog Number
1 plate	K151FKD-1
5 plates	K151FKD-2
20 plates	K151FKD-3
20 plates (Base)	K151FKA-3