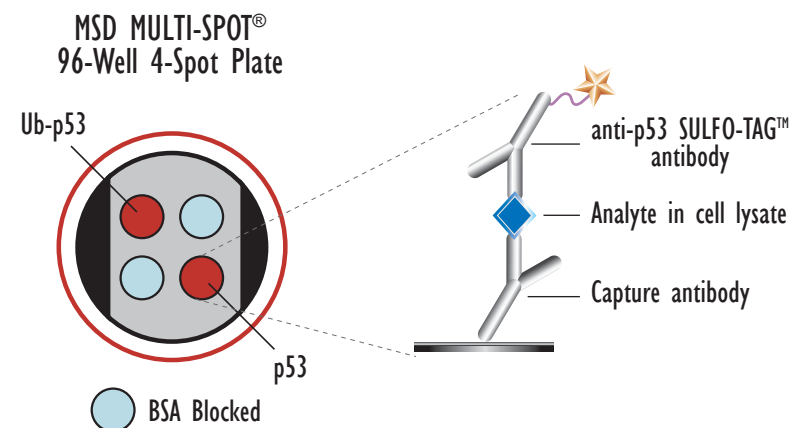
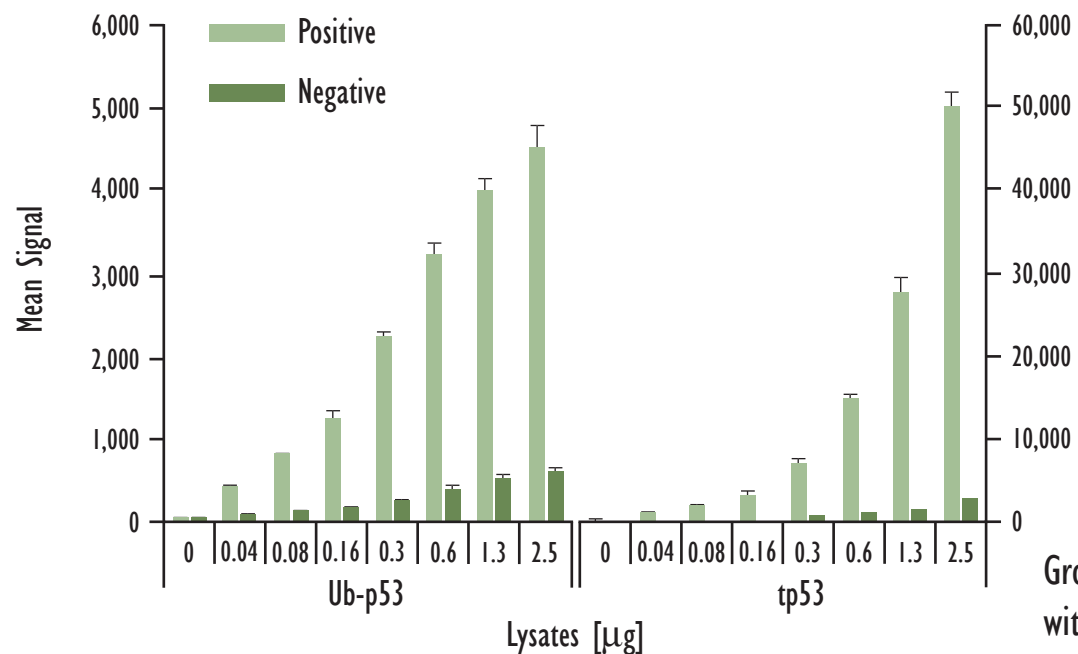


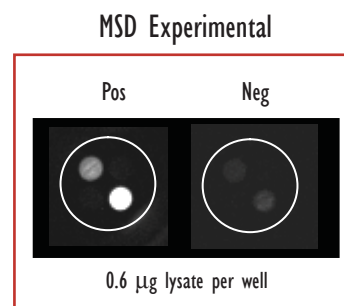
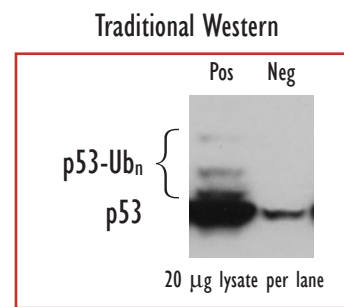
# MULTI-SPOT<sup>®</sup> Ubiquitinated and Total p53 Assay

Detection of Ubiquitinated and Total p53 in Whole Cell Lysate Samples



	Lysates (µg)	p53 Positive			p53 Negative			P/N
		Average	StdDev	%CV	Average	StdDev	%CV	
Ub-p53	0	41	10	26	47	11	23	
	0.04	411	15	4	78	19	24	5.3
	0.08	808	14	2	115	3	2	7.0
	0.16	1,235	86	7	152	2	1	8.1
	0.3	2,260	33	1	244	12	5	9.3
	0.6	3,252	108	3	404	20	5	8.0
	1.3	3,994	154	4	516	62	12	7.7
	2.5	4,550	238	5	587	39	7	7.8
tp53	0	38	16	41	37	14	38	
	0.04	897	21	2	90	1	1	10.0
	0.08	1,861	35	2	132	8	6	14.1
	0.16	3,100	315	10	200	27	13	15.5
	0.3	6,984	469	7	362	21	6	19.3
	0.6	14,481	638	4	713	12	2	20.3
	1.3	27,489	1,782	6	1,263	155	12	21.8
	2.5	50,122	1,751	3	2,383	226	9	21.0

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 and antibody against ubiquitinated proteins on two of the four spatially distinct electrodes per well. Total and ubiquitinated p53 were detected with an anti-total-p53 antibody labeled with MSD SULFO-TAG reagent.



Kit Size	SI2400	SI6000
1 plate	K15169D-1	K11169D-1
5 plates	K15169D-2	K11169D-2
20 plates	K15169D-3	K11169D-3
20 plates (Base)	K15169A-3	K11169A-3



Meso Scale Discovery  
A Division of Meso Scale Diagnostics, LLC.  
www.meso-scale.com