

MAPK Cascade as a Signaling Convergence Point in Cancer Therapy: Simultaneous Detection of Activated p38, ERK and JNK

Nisar Pampori, Paula Denney Eason,
Laura K. Schaefer, Jenny T. Ly,
Bruk G. Leta, Staci M. Leffel,
Paul J. Goodwin, Erin K. Teagan,
Shayla R. Workman, Robert M. Umek
and Jacob N. Wohlstadter



Meso Scale Discovery

A division of Meso Scale Diagnostics, LLC.

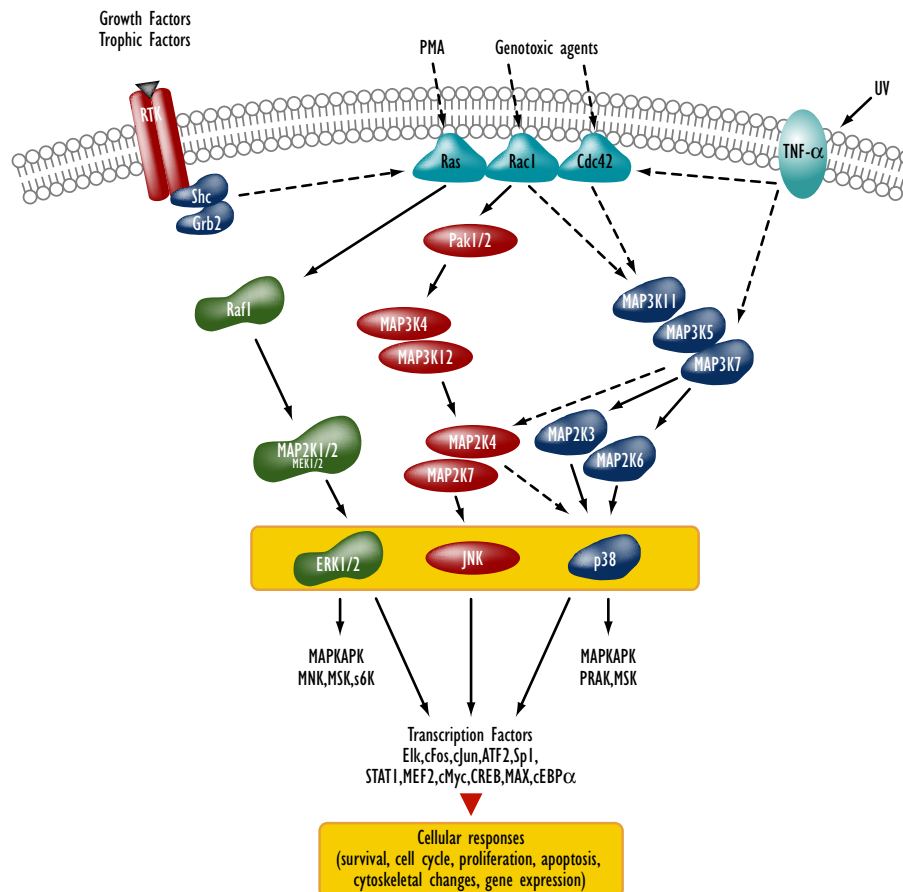
9238 Gaither Road, Gaithersburg, MD 20877
Phone: 240.631.2522 Fax: 240.632.2219
www.meso-scale.com

MAPK Cascade as a Signaling Convergence Point in Cancer Therapy: Simultaneous Detection of Activated p38, ERK and JNK

1 Abstract

The mitogen-activated protein kinase (MAPK) pathways, through major groupings like p38, ERK and JNK, constitute interrelated signal transduction cascades that regulate several diseases. In meeting the challenges of system biology and linear drug discovery, we present MSD's MULTI-SPOT® platform for simultaneous detection of phosphorylated forms of p38(T180/Y182), ERK1/2(T202/Y204)/(T185/Y187) and JNK(T183/Y185). The assays were performed in multiplex format and showed high sensitivity; phospho-p38 was quantifiable at attomoles, or from about 5,000 stimulated cells. High signal/background ratios were registered for phospho-p38 and phospho-JNK targets in UV treated HEK293 or NIH3T3 cells, and from the phospho-ERK1/2 target in Phorbol-ester treated Jurkat cells. Total pools of ERK1/2 and p38 were also detected quantitatively from cell lysates in the same well. These tools enable the rapid identification of drug candidates and promote the comprehensive understanding of the role of MAPK signaling pathway proteins.

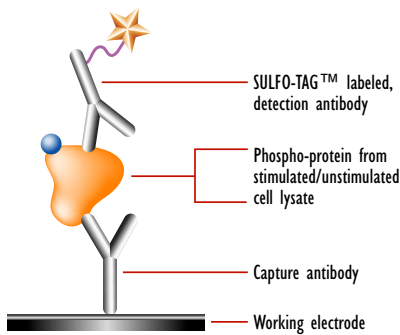
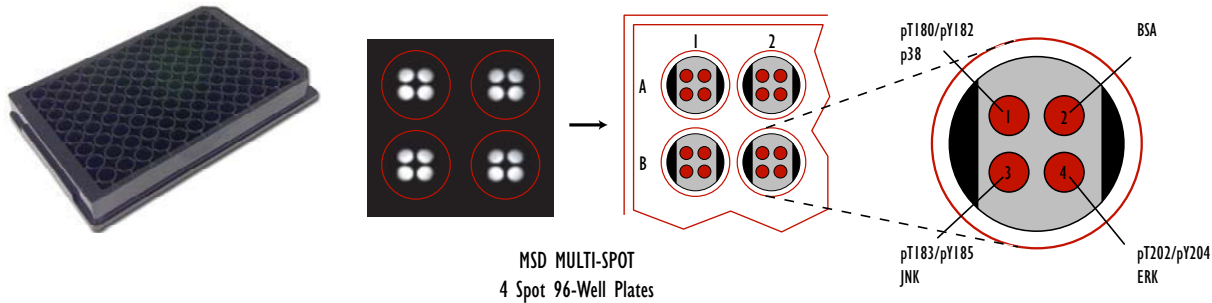
2 MAPK Pathway Convergence: ERK1/2, JNK and p38



Meso Scale Discovery
A division of Meso Scale Diagnostics, LLC.

MAPK Cascade as a Signaling Convergence Point in Cancer Therapy: Simultaneous Detection of Activated p38, ERK and JNK

3 MSD MULTI-ARRAY™ and MULTI-SPOT® Plates



1. MULTI-SPOT 4 Spot 96-Well Plates precoated with capture antibodies are blocked with 200 μ L of MSD Blocker-A for 1h and washed with TBS.
2. 25 μ L of cell lysates are added to the wells and incubated for 3h at RT or overnight at 4 $^{\circ}$ C with shaking, followed by washing with TBS.
3. 25 μ L MSD SULFO-TAG labeled antibodies (in Antibody dilution buffer) are added to the wells and incubated for 1h with shaking, followed by washing with TBS.
4. 150 μ L MSD Read Buffer T (with surfactant) are added to the wells and analyzed on the SECTOR™ 6000 instrument.

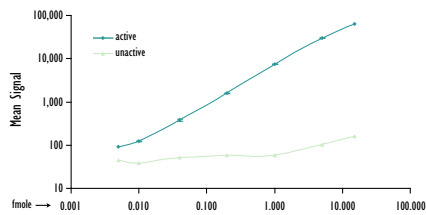


Meso Scale Discovery
A division of Meso Scale Diagnostics, LLC.

MAPK Cascade as a Signaling Convergence Point in Cancer Therapy: Simultaneous Detection of Activated p38, ERK and JNK

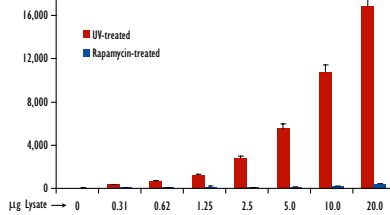
4 MSD Singleplex Phospho-p38 (T180/Y182) assay

in purified recombinant preps



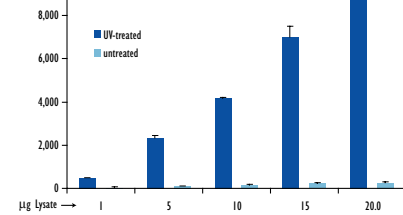
| Pure fmole | Active | | Inactive | | Active/ Inactive |
|---------------|--------|-----|----------|-----|---------------------|
| | Ave. | %CV | Ave. | %CV | |
| 0.0 | | | 48 | 22 | |
| 0.005 | 92 | 2 | 45 | 9 | 2.0 |
| 0.01 | 123 | 14 | 39 | 24 | 3.2 |
| 0.04 | 379 | 5 | 53 | 9 | 7.2 |
| 0.2 | 1,575 | 2 | 58 | 7 | 27.1 |
| 1 | 7,416 | 3 | 60 | 44 | 124.6 |
| 5 | 30,269 | | 104 | 16 | 292.5 |
| 15 | 62,617 | 0 | 164 | 13 | 383.0 |

in HEK293 Whole Cell Lysates

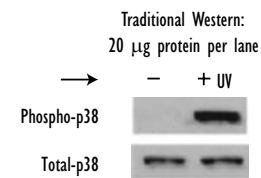
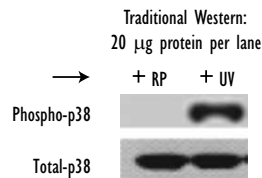


| HEK293 µg | UV-Treated | | RP-Treated | | Active/ Inactive |
|--------------|------------|-----|------------|-----|---------------------|
| | Ave. | %CV | Ave. | %CV | |
| 0.0 | | | 36 | 8 | |
| 0.31 | 334 | 7 | 45 | 27 | 7.5 |
| 0.62 | 623 | 15 | 47 | 30 | 13.3 |
| 1.25 | 1,271 | 1 | 123 | 69 | 10.3 |
| 2.5 | 2,852 | 5 | 85 | 7 | 33.6 |
| 5.0 | 5,531 | 8 | 124 | 13 | 44.8 |
| 10.0 | 10,759 | 6 | 199 | 5 | 54.1 |
| 20.0 | 16,929 | 5 | 402 | 5 | 42.1 |

in NIH3T3 Whole Cell Lysates



| NIH3T3 µg | UV-Treated | | Untreated | | Active/ Inactive |
|--------------|------------|-----|-----------|-----|---------------------|
| | Ave. | %CV | Ave. | %CV | |
| 1.0 | 488 | 3 | 53 | 17 | 9.3 |
| 5.0 | 2,329 | 5 | 114 | 8 | 20.4 |
| 10.0 | 4,159 | 2 | 155 | 15 | 26.9 |
| 15.0 | 7,007 | 7 | 235 | 13 | 29.8 |
| 20.0 | 9,467 | 0 | 279 | 4 | 34.0 |



Growing HEK293 cells were treated with Rapamycin (1 µM; 3 hr)(+RP), while serum deprived HEK293 cells were harvested 30 minutes after UV irradiation (40mJ/cm²)(+UV). Serum deprived NIH3T3 cells (-) were harvested 30 minutes after UV irradiation (40mJ/cm²)(+UV). Whole cell lysates were added to MSD MULTI-SPOT 4 Spot 96-well plates coated with phospho-p38 antibody on one of the four spatially distinct electrodes per well. Phosphorylated p38 was detected with 10 nM anti-p38 antibody labeled with MSD SULFO-TAG reagents.

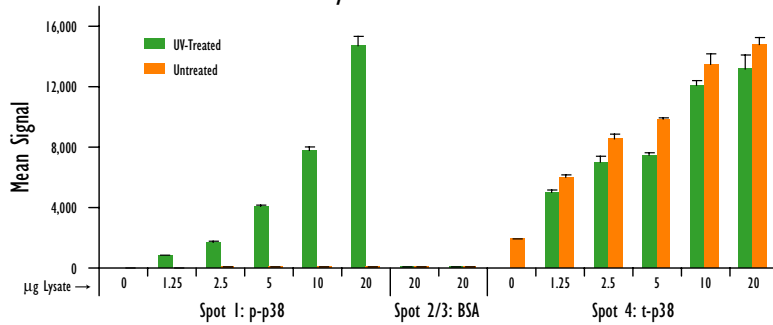


Meso Scale Discovery
A division of Meso Scale Diagnostics, LLC.

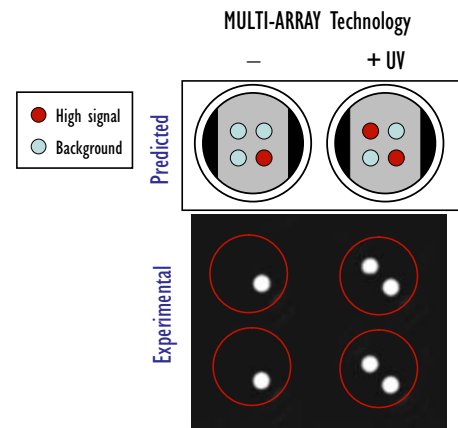
MAPK Cascade as a Signaling Convergence Point in Cancer Therapy: Simultaneous Detection of Activated p38, ERK and JNK

5 MSD Multiplexed Phospho-p38 / Total-p38 assay

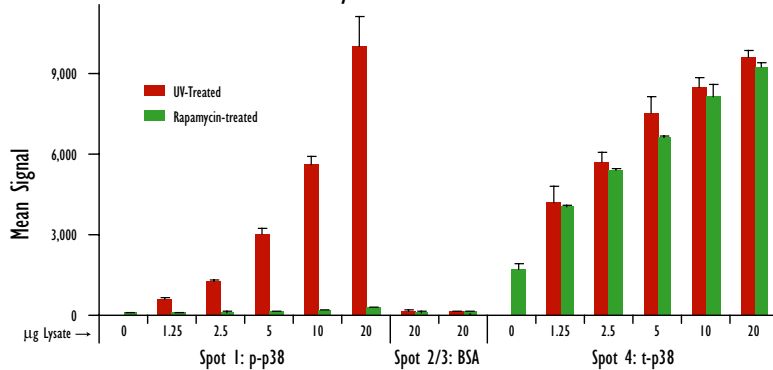
in NIH3T3 Whole Cell Lysates



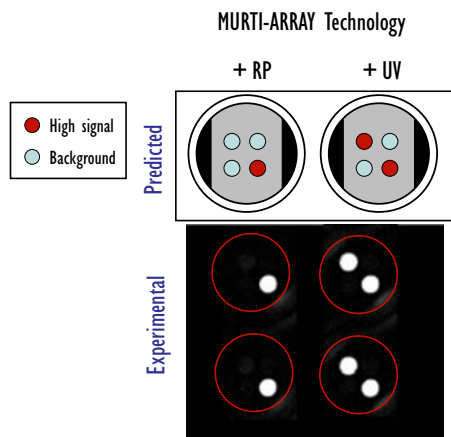
| Spot | NIH3T3 µg | UV-treated | | Untreated | | Active/ Inactive |
|--------|--------------|------------|-----|-----------|-----|---------------------|
| | | Ave. | %CV | Ave. | %CV | |
| 1:pp38 | 0.0 | | | 27 | 27 | |
| 1:pp38 | 1.25 | 847 | 2 | 34 | 13 | 25.2 |
| 1:pp38 | 2.5 | 1,706 | 5 | 47 | 35 | 36.0 |
| 1:pp38 | 5.0 | 4,071 | 2 | 66 | 10 | 62.0 |
| 1:pp38 | 10.0 | 7,821 | 3 | 81 | 10 | 97.0 |
| 1:pp38 | 20.0 | 14,750 | 5 | 99 | 6 | 148.5 |
| 2:BSA | 20.0 | 52 | 35 | 47 | 30 | |
| 3:BSA | 20.0 | 63 | 23 | 55 | 8 | |
| 4:tp38 | 0.0 | | | 1,916 | 2 | |
| 4:tp38 | 1.25 | 5,051 | 3 | 6,042 | 2 | 0.8 |
| 4:tp38 | 2.5 | 7,050 | 5 | 8,603 | 3 | 0.8 |
| 4:tp38 | 5.0 | 7,505 | 2 | 9,877 | 1 | 0.8 |
| 4:tp38 | 10.0 | 12,165 | 2 | 13,528 | 5 | 0.9 |
| 4:tp38 | 20.0 | 13,213 | 7 | 14,811 | 3 | 0.9 |



in HEK293 Whole Cell Lysates



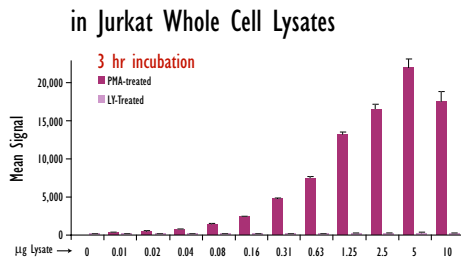
| Spot | HEK293 µg | UV-treated | | RP-treated | | Active/ Inactive |
|--------|--------------|------------|-----|------------|-----|---------------------|
| | | Ave. | %CV | Ave. | %CV | |
| 1:pp38 | 0.0 | | | 82 | 19 | |
| 1:pp38 | 1.25 | 604 | 6 | 113 | 4 | 5.3 |
| 1:pp38 | 2.5 | 1,243 | 4 | 126 | 6 | 9.9 |
| 1:pp38 | 5.0 | 3,028 | 6 | 142 | 12 | 21.4 |
| 1:pp38 | 10.0 | 5,578 | 6 | 192 | 12 | 29.1 |
| 1:pp38 | 20.0 | 10,011 | 11 | 319 | 0 | 31.4 |
| 2:BSA | 20.0 | 159 | 16 | 121 | 5 | |
| 3:BSA | 20.0 | 164 | 4 | 135 | 9 | |
| 4:tp38 | 0.0 | | | 1,721 | 12 | |
| 4:tp38 | 1.25 | 4,199 | 14 | 4,013 | 2 | 1.0 |
| 4:tp38 | 2.5 | 5,695 | 7 | 5,376 | 1 | 1.1 |
| 4:tp38 | 5.0 | 7,505 | 9 | 6,609 | 1 | 1.1 |
| 4:tp38 | 10.0 | 8,499 | 4 | 8,101 | 6 | 1.0 |
| 4:tp38 | 20.0 | 9,584 | 2 | 9,239 | 1 | 1.0 |



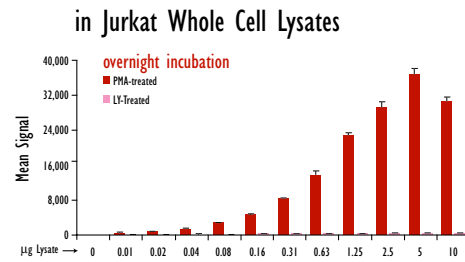
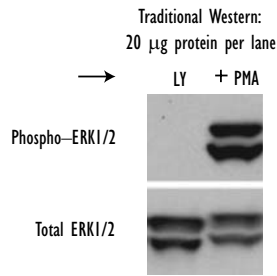
Meso Scale Discovery
A division of Meso Scale Diagnostics, LLC.

MAPK Cascade as a Signaling Convergence Point in Cancer Therapy: Simultaneous Detection of Activated p38, ERK and JNK

6 MSD Singleplex Phospho-ERK1/2 (T/Y:202/204;185/187) assay



| Jurkat µg | PMA-Treated | | LY-Treated | | S/B | Active/ Unactive |
|--------------|-------------|-----|------------|-----|-------|---------------------|
| | Ave | %CV | Ave | %CV | | |
| 0 | | | 87 | 11 | | |
| 0.01 | 280 | 3 | 89 | 12 | 3.2 | 3.1 |
| 0.02 | 449 | 5 | 81 | 6 | 5.2 | 5.6 |
| 0.039 | 716 | 5 | 68 | 11 | 8.3 | 10.6 |
| 0.078 | 1,398 | 1 | 85 | 14 | 16.1 | 16.4 |
| 0.16 | 2,404 | 1 | 92 | 7 | 27.7 | 26.2 |
| 0.31 | 4,652 | 3 | 120 | 3 | 53.1 | 38.9 |
| 0.63 | 7,376 | 4 | 141 | 1 | 85.1 | 52.2 |
| 1.25 | 13,072 | 2 | 181 | 2 | 150.8 | 72.1 |
| 2.5 | 16,483 | 4 | 219 | 1 | 190.2 | 75.4 |
| 5 | 21,853 | 5 | 257 | 7 | 252.1 | 84.9 |
| 10 | 17,330 | 8 | 244 | 4 | 200 | 71.1 |



| Jurkat µg | PMA-Treated | | LY-Treated | | S/B | Active/ Unactive |
|--------------|-------------|-----|------------|-----|-------|---------------------|
| | Ave | %CV | Ave | %CV | | |
| 0 | | | 68 | 27 | | |
| 0.01 | 449 | 4 | 74 | 5 | 6.6 | 6 |
| 0.02 | 771 | 2 | 67 | 11 | 11.3 | 11.4 |
| 0.04 | 1,334 | 3 | 82 | 17 | 19.5 | 16.3 |
| 0.08 | 2,661 | 2 | 82 | 9 | 38.9 | 32.6 |
| 0.16 | 4,537 | 5 | 100 | 5 | 66.4 | 45.4 |
| 0.31 | 8,316 | 1 | 122 | 5 | 121.7 | 68.2 |
| 0.63 | 13,688 | 6 | 177 | 6 | 200.3 | 77.5 |
| 1.25 | 22,746 | 3 | 226 | 8 | 332.9 | 100.6 |
| 2.5 | 29,147 | 4 | 314 | 3 | 426.5 | 92.9 |
| 5 | 36,659 | 3 | 349 | 6 | 536.5 | 105.1 |
| 10 | 30,455 | 3 | 347 | 1 | 445.7 | 87.7 |

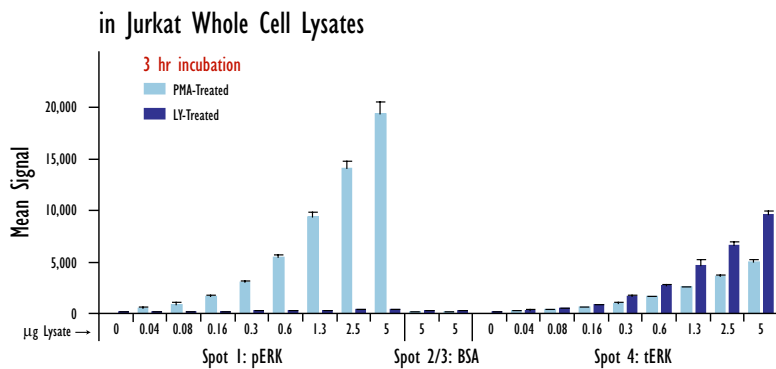
Logarithmically growing Jurkat cells were treated with LY294002 inhibitor (50 µM; 2.25 hr)(LY) or PMA (200 nM; 15 minutes)(+PMA). Whole cell lysates were added to MSD MULTI-SPOT 4 Spot 96-well plates coated with anti ERK1/2 antibody on one of the four spatially distinct electrodes per well. Phosphorylated ERK1/2 was detected with 10 nM anti phospho-ERK1/2 antibody labeled with MSD SULFO-TAG reagents.



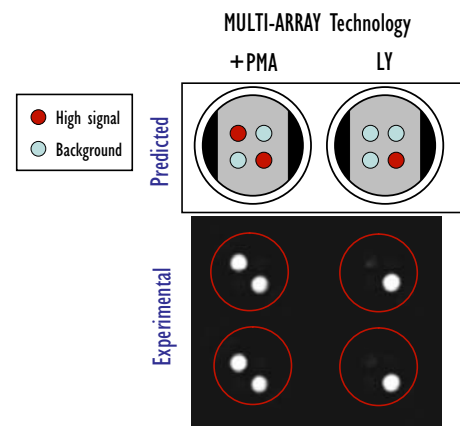
Meso Scale Discovery
A division of Meso Scale Diagnostics, LLC.

MAPK Cascade as a Signaling Convergence Point in Cancer Therapy: Simultaneous Detection of Activated p38, ERK and JNK

7 MSD Multiplexed Phospho-ERK1/2 / Total-ERK1/2 assay



| Spot | Jurkat µg | PMA-treated | | LY-Treated | | Active/Unactive |
|--------|-----------|-------------|-----|------------|-----|-----------------|
| | | Ave. | %CV | Ave. | %CV | |
| 1:pERK | 0 | | | 143 | 9 | |
| 1:pERK | 0.02 | 343 | 6 | 154 | 10 | 2.2 |
| 1:pERK | 0.04 | 511 | 4 | 134 | 5 | 3.8 |
| 1:pERK | 0.08 | 849 | 22 | 145 | 5 | 5.9 |
| 1:pERK | 0.16 | 1,660 | 3 | 148 | 5 | 11.2 |
| 1:pERK | 0.31 | 2,979 | 6 | 181 | 4 | 16.4 |
| 1:pERK | 0.63 | 5,447 | 4 | 217 | 5 | 25.1 |
| 1:pERK | 1.25 | 9,324 | 5 | 267 | 7 | 34.9 |
| 1:pERK | 2.5 | 14,091 | 5 | 317 | 1 | 44.5 |
| 1:pERK | 5.0 | 19,386 | 6 | 385 | 4 | 50.4 |
| 2:BSA | 5.0 | 154 | 7 | 215 | 5 | |
| 3:BSA | 5.0 | 161 | 4 | 237 | 3 | |
| 4:tERK | 0 | | | 151 | 9 | |
| 4:tERK | 0.02 | 197 | 6 | 205 | 5 | 1.0 |
| 4:tERK | 0.04 | 238 | 3 | 281 | 4 | 0.8 |
| 4:tERK | 0.08 | 351 | 6 | 444 | 1 | 0.8 |
| 4:tERK | 0.16 | 581 | 5 | 812 | 4 | 0.7 |
| 4:tERK | 0.31 | 973 | 2 | 1,629 | 3 | 0.6 |
| 4:tERK | 0.63 | 1,625 | 1 | 2,661 | 2 | 0.6 |
| 4:tERK | 1.25 | 2,500 | 4 | 4,590 | 13 | 0.5 |
| 4:tERK | 2.5 | 3,580 | 4 | 6,588 | 5 | 0.5 |
| 4:tERK | 5.0 | 4,999 | 3 | 9,544 | 4 | 0.5 |



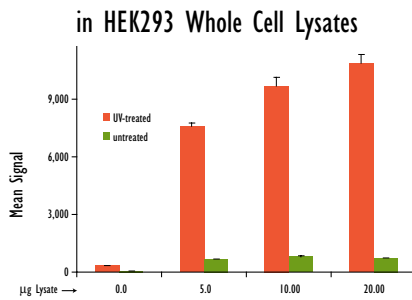
Logarithmically growing Jurkat cells were treated with LY294002 (50 nM; 2.25 hr)(-) or PMA (200 nM; 15 minutes)(+). Whole cell lysates were added to MSD MULTI-SPOT 4 Spot 96-well plates coated with anti ERK1/2 antibody on two of the four spatially distinct electrodes per well. Phosphorylated and total ERK1/2 were detected with 10 nM anti-total-ERK1/2 antibody labeled with MSD SULFO-TAG reagent.



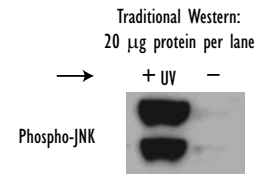
Meso Scale Discovery
A division of Meso Scale Diagnostics, LLC.

MAPK Cascade as a Signaling Convergence Point in Cancer Therapy: Simultaneous Detection of Activated p38, ERK and JNK

8 MSD Singleplex Phospho-JNK (T183/Y185) assay



| HEK293 µg | UV-Treated | | Untreated | | Active/ Unactive |
|--------------|------------|-----|-----------|-----|---------------------|
| | Ave. | %CV | Ave. | %CV | |
| 0.0 | 358 | 2 | 41 | 14 | |
| 5.0 | 7,606 | 2 | 679 | 1 | 11.2 |
| 10.0 | 9,710 | 4 | 787 | 6 | 12.3 |
| 20.0 | 10,888 | 4 | 722 | 1 | 15.1 |



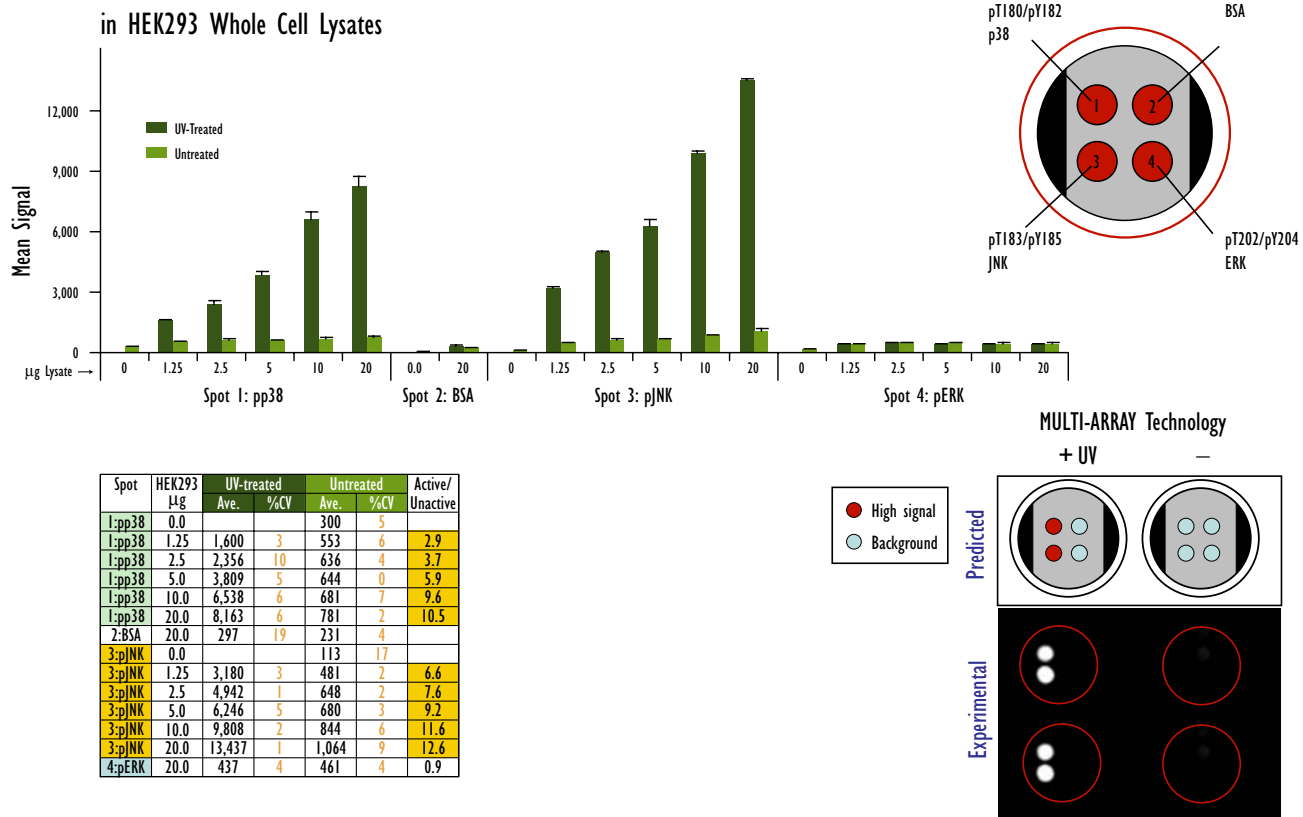
Logarithmically growing HEK293 cells were treated with +/- UV. Whole cell lysates were added to MSD MULTI-SPOT 4 Spot 96-well plates coated with capture antibody on one of the four spatially distinct electrodes per well. Phosphorylated JNK was detected with 10 nM anti-phospho-JNK antibody labeled with MSD SULFO-TAG reagents.



Meso Scale Discovery
A division of Meso Scale Diagnostics, LLC.

MAPK Cascade as a Signaling Convergence Point in Cancer Therapy: Simultaneous Detection of Activated p38, ERK and JNK

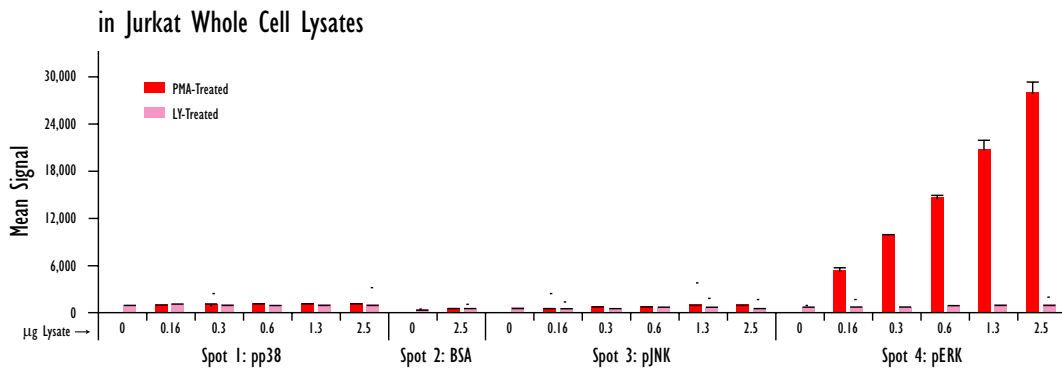
9 MSD Multiplexed MAPK Panel : p-p38 / p-ERK / p-JNK



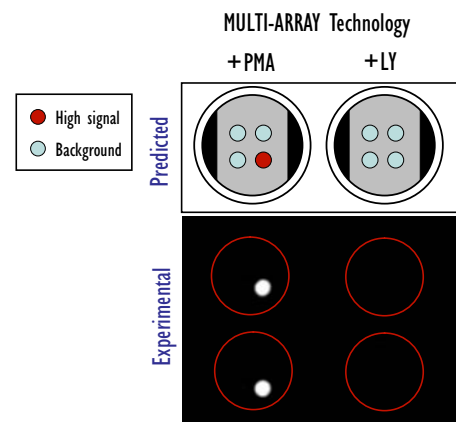
Logarithmically growing HEK293 cells were treated with +/- UV. Whole cell lysates were added to MSD MULTI-SPOT 4 Spot 96-well plates coated with capture antibody on three of the four spatially distinct electrodes per well. Phosphorylated proteins were detected with MSD SULFO-TAG detection antibodies.

MAPK Cascade as a Signaling Convergence Point in Cancer Therapy: Simultaneous Detection of Activated p38, ERK and JNK

9 MSD Multiplexed MAPK Panel : p-p38 / p-ERK / p-JNK (continued)



| Spot | Jurkat µg | PMA-treated | | LY-Treated | | Active/ Unactive |
|--------|--------------|-------------|-----|------------|-----|---------------------|
| | | Ave. | %CV | Ave. | %CV | |
| 1:pp38 | 0 | | | 808 | 9 | |
| 1:pp38 | 2.5 | 907 | 5 | 733 | 1 | 1.2 |
| 2:BSA | 2.5 | 445 | 4 | 337 | 1 | 1.3 |
| 3:pJNK | 0 | | | 354 | 11 | |
| 3:pJNK | 2.5 | 801 | 3 | 457 | 6 | 1.8 |
| 4:pERK | 0 | | | 486 | 17 | |
| 4:pERK | 0.156 | 5,220 | 8 | 526 | 12 | 9.9 |
| 4:pERK | 0.313 | 9,734 | 1 | 605 | 5 | 16.1 |
| 4:pERK | 0.625 | 14,366 | 3 | 673 | 12 | 21.3 |
| 4:pERK | 1.25 | 20,685 | 5 | 774 | 10 | 26.7 |
| 4:pERK | 2.5 | 27,770 | 5 | 766 | 15 | 36.3 |



Logarithmically growing Jurkat cells were treated with LY294002 inhibitor (50 μ M; 2.25 hr)(LY) or PMA (200 nM; 15 minutes)(+PMA). Whole cell lysates were added to MSD MULTI-SPOT 4 Spot 96-well plates coated with capture antibody on three of the four spatially distinct electrodes per well. Phosphorylated proteins were detected with MSD SULFO-TAG detection antibodies.

MAPK Cascade as a Signaling Convergence Point in Cancer Therapy: Simultaneous Detection of Activated p38, ERK and JNK

10 Conclusion

1. We present highly specific multiplexed assays for simultaneous detection of phosphorylated MAPK proteins, phospho-p38 (pT180/pY182), phospho-ERK1/2 (pT202/pY204 and pT185/pY187) and phospho-JNK (pT183/pY185).
2. These assays specifically identify phosphoproteins from human and murine cell lysates as well as recombinant purified proteins. The phospho-p38 detection assay identifies attomole levels of recombinant purified active protein with high sensitivity; signal / background ratio of about 380 (active vs inactive protein).
3. The phosphoprotein assays could be run in duplex format, thus simultaneously determining the phosphorylated and total levels of the target protein.
4. The methods we present are general, demonstrating that multiple phosphoprotein members of signaling pathways can be assayed simultaneously in a single well in short (3 hr) or long (overnight) incubation periods, using specific antibodies immobilized on MSD MULTI-SPOT plates.
5. The MULTI-ARRAY technology-based assays are useful supplements to gold-standard methods like western blots because these versatile assays are highly quantitative and easily automated, suitable for HTS, and save time and labor compared to existing techniques.

