

# Bio-Plex™ suspension array system

tech note 5394

## Bio-Plex Phosphoprotein Detection References

### Cell Culture

| Bio-Plex Assay   | Reference  |
|--|--|
| p-ERK2<br>p-IκB-α<br>p-JNK<br>p-p38 MAPK<br><br>t-ERK2<br>t-IκB-α<br>t-JNK | Bharadwaj U et al., Effects of cyclophilin A on myeloblastic cell line KG-1 derived dendritic like cells (DLC) through p38 MAP kinase activation, J Surg Res 127, 29–38 (2005)   |
| p-Akt<br>p-ATF-2<br>p-ERK1/2<br>p-GSK-3α/β<br>p-p38 MAPK                   | Comer JE et al., Direct inhibition of T-lymphocyte activation by anthrax toxins in vivo, Infect Immun 73, 8275–8281 (2005)   |
| p-ERK1/2<br>p-JNK<br>p-p38 MAPK<br><br>t-ERK2<br>t-JNK<br>t-p38 MAPK       | Li M et al., Transfection of SSTR-1 and SSTR-2 inhibits Panc-1 cell proliferation and renders Panc-1 cells responsive to somatostatin analogue, J Am Coll Surg 201, 571–578 (2005)   |
| p-ATF-2<br>p-EGFR<br>p-ERK1/2<br>p-IκB-α<br>p-JNK<br>p-p38 MAPK<br>p-Stat3 | Sakai K et al., Dimerization and the signal transduction pathway of a small in-frame deletion in the epidermal growth factor receptor, FASEB J 20, 311–313 (2006)  |
| p-ERK1/2   | Titus B et al., Endothelin axis is a target of the lung metastasis suppressor gene <i>RhoGDI2</i> , Cancer Res 65, 7320–7327 (2005)  |
| p-ERK1/2<br>p-JNK<br>p-p38 MAPK  | Wang T et al., Co-activation of ERK, NF-κB, and GADD45β in response to ionizing radiation, J Biol Chem 280, 12593–12601 (2005)   |
| p-JNK  | Wang X et al., Complete inhibition of anisomycin and UV radiation but not cytokine induced JNK and p38 activation by an aryl-substituted dihydropyrrlopyrazole quinoline and mixed lineage kinase 7 small interfering RNA, J Biol Chem 280, 19298–19305 (2005) |
| p-ERK1/2<br>p-JNK<br>p-p38 MAPK<br><br>t-ERK2<br>t-JNK                     | Yan S et al., Effects of lysophosphatidylcholine on monolayer cell permeability of human coronary artery endothelial cells, Surgery 138, 464–473 (2005)  |

## Tissue

| Bio-Plex Assay  | Reference  |
|---|--|
| p-I $\kappa$ B- $\alpha$<br>p-JNK<br>p-p38 MAPK                                   | Austin BA et al., Critical role for the oligoadenylate synthetase/RNase L pathway in response to IFN- $\beta$ during acute ocular herpes simplex virus type 1 infection, J Immunol 175, 1100–1106 (2005)                         |
| p-ATF-2<br>p-ERK1/2<br>p-I $\kappa$ B- $\alpha$<br>p-JNK<br>p-p38 MAPK<br>p-Stat3 | Turner TT et al., Testicular torsion alters the presence of specific proteins in the mouse testis as well as the phosphorylation status of specific proteins, J Androl 27, 285–293 (2006)  |
| p-ERK1/2  | Wilhelm SM et al., BAY 43-9006 exhibits broad spectrum oral antitumor activity and targets the RAF/MEK/ERK pathway and receptor tyrosine kinases involved in tumor progression and angiogenesis, Cancer Res 64, 7099–7109 (2004) |

p = Phosphoprotein assay

t = Total target assay



**Bio-Rad**  
**Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 4BIORAD **Australia** 02 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 21 3237 9400  
**Canada** 905 712 2771 **China** 86 21 6426 0808 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65  
**Germany** 089 318 84 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 455 8800 **India** 91 124 4029300/5013478 **Israel** 03 963 6050  
**Italy** 39 02 216091 **Japan** 03 5811 6270 **Korea** 82 2 3473 4460 **Mexico** 55 5200 05 20 **The Netherlands** 0318 540666 **New Zealand** 64 9415 2280  
**Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 095 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 0861 246 723  
**Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** 886 2 2578 7189/2578 7241 **United Kingdom** 020 8328 2000