

Bio-Plex Phosphoprotein Detection References

Cell Culture

Bio-Plex Assay	Reference
p-ERK2 p-IkB- α p-JNK p-p38 MAPK t-ERK2 t-IkB- α 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 p-ATF-2 p-ERK1/2 p-GSK-3 α / β 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 p-JNK p-p38 MAPK t-ERK2 t-JNK 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 p-EGFR p-ERK1/2 p-IkB- α p-JNK p-p38 MAPK 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 p-JNK p-p38 MAPK	Wang T et al., Co-activation of ERK, NF- κ B, and GADD45 β 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 p-JNK p-p38 MAPK t-ERK2 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 κ B- α p-JNK p-p38 MAPK	Austin BA et al., Critical role for the oligoadenylate synthetase/RNase L pathway in response to IFN- β during acute ocular herpes simplex virus type 1 infection, J Immunol 175, 1100–1106 (2005)
p-ATF-2 p-ERK1/2 p-I κ B- α p-JNK p-p38 MAPK 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 **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