

## Examples of known RhoA activators

| Activator*   | Treatment                                | Cell Line                                  | Response   | Type of Assay                           | Citation   |
|--|--|--|--|---|--|
| Lysophosphatidylcholine (lysophospholipid)               | 100 nM                                   | human melanocytes                          | After 1 min, 1.5 fold increase over control. Activation peaked at 5 min (5.3 fold increase) and declined to baseline by 60 min   | Rho G-LISA (Cat. # BK124)               | Scott et al., 2007. J Invest Dermatol. 127, 668  |
| Calpeptin (Cat. # CN01)                                  | 0.5 mU/mL                                | uterine myocytes                           | After 15 min treatment, 2.4 fold increase over vehicle control   | RhoA G-LISA (Cat. # BK124)              | Aguilar et al., 2011. PLoS ONE 6(6): e20903. doi:10.1371/journal.pone.0020903                  |
| Colchicine (microtubule destabilizer)                    | 10 µg/ml                                 | Swiss 3T3 cells, adherent or suspension    | Maximal activation of 2-4 fold activation after 30 min   | Rhotekin-RBD pulldown                   | Ren et al., 1999. EMBO J. 18, 578  |
| Nocodazole (microtubule destabilizer)                    | 10 µM                                    | MG63 human osteosarcoma cells & HeLa cells | Maximal activation of 2-3 fold activation after 30 min   | Actin morphology, rhotekin-RBD pulldown | Zhang et al., 1997. Mol Biol Cell. 8, 1415<br>Maddox and Burridge, 2003. J Cell Biol. 160, 255 |
| Vinblastine (microtubule destabilizer)                   | 50 µM                                    | MG63 human osteosarcoma cells              | Maximal activation of 2-4 fold activation after 30 min   | Actin morphology                        | Zhang et al., 1997. Mol Biol Cell. 8, 1415   |
| Cytochalasin D (actin filament destabilizer)             | 0.5 µg/ml                                | Swiss 3T3 cells, adherent or suspension    | Maximal activation of 1-2 fold after 60 min  | Rhotekin-RBD pulldown                   | Ren et al., 1999. EMBO J. 18, 578  |
| Sphingosine -1-phosphate (serum lipid & GPCR agonist)    | 1 µg/ml                                  | Swiss 3T3 cells, adherent or suspension    | Maximal activation of 1-2 fold after 2 min for 3T3 cells and 20 min for HUVEC cells  | Rhotekin-RBD pulldown                   | Ren et al., 1999. EMBO J. 18, 578<br>Vouret-Craviari et al., 2002. J Cell Sci. 115, 2475       |
| Serum  | 5 - 10%                                  | Swiss 3T3 cells, adherent or suspension    | Maximal activation of 2-6 fold (10%) and 1-2 fold (5%) after 1-5 min   | Rhotekin-RBD pulldown                   | Ren et al., 1999. EMBO J. 18, 578  |
| Lysophosphatidic acid (LPA) (serum lipid & GPCR agonist) | 1 µg/ml                                  | Swiss 3T3 cells, adherent or suspension    | Maximal activation of 2-6 fold after 1 min then dropping to basal after 30 min   | Rhotekin-RBD pulldown                   | Ren et al., 1999. EMBO J. 18, 578  |
| Lysophosphatidic acid (LPA) (serum lipid & GPCR agonist) | 1 µM                                     | N1E-115 neuronal cells                     | Maximal activation of 3-5 fold after 3 min   | Rho-kinase pull down assay              | Kranenburg et al., 1999. Mol Biol Cell. 10, 1851   |
| Fibronectin (extracellular matrix protein)               | Culture plate is coated with fibronectin | Swiss 3T3 cells                            | Biphasic regulation after plating cells on fibronectin coated plates. Initial period of low RhoA activity (10-20 min) followed by a 1-7 fold activation peaking at 60-90 minutes and then dropping to basal levels after 6 h | Rhotekin-RBD pulldown                   | Ren et al., 1999. EMBO J. 18, 578  |
| Bombesin   | 10 nM                                    | Swiss 3T3 cells                            | Maximal activation of 2-3 fold after 1 min which is sustained for at least 30 min  | Actin morphology                        | Ridley and Hall, 1992. Cell. 70, 389   |

## Examples of known Rac activators

| Activator*   | Treatment  | Cell Line   | Response  | Type of Assay                      | Citation  |
|--|--|---|---|------------------------------------|---|
| Epidermal Growth Factor (Cat. # CN02)  | 10 ng/ml   | Swiss 3T3 cells   | Maximal activation after 2 min  | Rac G-LISA (Cat. # BK126 or BK128) | In-house QC results                                       |
| Epidermal Growth Factor  | 50 ng/ml   | U87MG human glioblastoma cells  | Maximal activation at 5 min. With 2D cultures, 55% increase over control and with 3D cultures, 25% over control | Rac G-LISA                         | Kim et al., 2008. Mol Biol Cell. 19, 4249                 |
| MCP-1 (chemokine and ligand for GPCR CCR2)                                       | 10 ng/ml   | murine alveolar macrophages; J774 macrophages murine peritoneal macrophages | maximal activation with 4 h treatment   | Rac G-LISA; PAK-PBD pulldown       | Tanaka et al., 2010. Biochem Biophys Res Comm. 399, 677   |
| heregulin beta 1 (ligand for the ErbB3/ ErbB4 tyrosine kinase-coupled receptors) | 0-30 ng/ml; used 10 ng/ml for time course              | breast cancer cell lines MCF-7 and T-47D serum-starved for 48 h             | dose-dependent effect with activation at 2 min, maximal by 10 min (MCF) or 5 min (T-47D)                        | PAK-PBD pulldown                   | Yang et al., 2006. Molec and Cellular Biol. 26, 831       |
| Epidermal Growth Factor  | 100 ng/ml  | breast cancer cell lines MCF-7 and T-47D serum-starved for 48 h             | maximal activation by 2 min   | PAK-PBD pulldown                   | Yang et al., 2006. Molec and Cellular Biol. 26, 831       |
| hyperosmotic exposure  | 100 mM NaCl added into the isotonic medium for 10 min  | neutrophils   | 2.5 fold increase over control  | PAK-PBD pulldown                   | Lewis et al., 2002. Am J Physiol Cell. Physiol. 282, C271 |
| hyperosmotic exposure  | 200 mM sucrose added, a non-ionic osmolyte, for 10 min | neutrophils   | Est. 2.5 fold increase over control   | PAK-PBD pulldown                   | Lewis et al., 2002. Am J Physiol Cell. Physiol. 282, C271 |
| Interleukin-3  | 5 ug/ml for 5 min                                      | MC/9 cells (mouse mast cell-like cell line)                                 | 2.4 fold increase over control. maximal activation after 5 min and maintained for at least 10 min               | PAK-PBD pulldown                   | Grill and Schrader, 2002. Blood. 100, 3183                |
| Interleukin-3  | 5 ug/ml for 5 min                                      | primary bone marrow-derived mast cells (BMMCs)                              | 4.3 fold increase over control  | PAK-PBD pulldown                   | Grill and Schrader, 2002. Blood. 100, 3183                |
| colony stimulating factor-1 (CSF-1)  | 200 ng/ml, 1-10 min                                    | WEHI 274.3, myelomonocytic cell line  | maximal at 1 min (12 fold), signif drop by 5 min (8.5 fold) and at 10 min (4.5 fold) compared to control        | PAK-PBD pulldown                   | Grill and Schrader, 2002. Blood. 100, 3183                |
| Steel locus factor (SLF)   | 50 ng/ml, 1-10 min                                     | MC/9 cells  | maximal at 1 min (7 fold increase over control) and dropped by 5-10 min (4 fold increase)                       | PAK-PBD pulldown                   | Grill and Schrader, 2002. Blood. 100, 3183                |
| Steel locus factor (SLF)   | 50 ng/ml, 5 min  | primary bone marrow-derived mast cells (BMMCs)                              | 3.5 fold increase over control  | PAK-PBD pulldown                   | Grill and Schrader, 2002. Blood. 100, 3183                |
| granulocyte-macrophage colony-stimulating factor (GM-CSF)                        | 10 ug/ml for 5 min                                     | MC/9 cells (mouse mast cell-like cell line)                                 | maximal at 1 min, maintained for at least 10 min  | PAK-PBD pulldown                   | Grill and Schrader, 2002. Blood. 100, 3183                |

## Examples of known Cdc42 activators

| Activator*   | Treatment  | Cell Line   | Response                            | Type of Assay               | Citation   |
|--|--|---|-------------------------------------|-----------------------------|--|
| Epidermal Growth Factor (Cat. # CN02)  | 100 ng/ml  | Swiss 3T3 cells   | Maximal activation after 1 min      | Cdc42 G-LISA (Cat. # BK127) | In-house QC results                                      |
| heregulin beta 1 (ligand for the ErbB3/ ErbB4 tyrosine kinase-coupled receptors) | 10 ng/ml   | breast cancer cell lines MCF-7 and T-47D serum-starved for 48 h | maximal by 5 min. treated 1-60 min  | PAK-PBD pulldown            | Yang et al., 2006. Molec and Cellular Biol. 26, 831      |
| hyperosmotic exposure  | 100 mM NaCl added into the isotonic medium for 10 min  | neutrophils   | 7.2 fold increase over control      | PAK-PBD pulldown            | Lewis et al., 2002. Am J Physiol Cell Physiol. 282, C271 |
| hyperosmotic exposure  | 200 mM sucrose added, a non-ionic osmolyte, for 10 min | neutrophils   | Est. 7.2 fold increase over control | PAK-PBD pulldown            | Lewis et al., 2002. Am J Physiol Cell Physiol. 282, C271 |
| Interleukin-3  | 5 ug/ml for 5 min                                      | primary bone marrow-derived mast cells (BMMCs)                  | 2 fold increase over control        | PAK-PBD pulldown            | Grill and Schrader, 2002. Blood. 100, 3183               |