

## PAK1 PBD Agarose Beads

---

**CATALOG NUMBER:** STA-411

**STORAGE:** -20°C

**QUANTITY AND CONCENTRATION:** 800  $\mu$ L of 50% Agarose slurry, 400  $\mu$ g PAK1-PBD in 1X PBS, 50% Glycerol

**SHELF LIFE:** 1 year from receipt under proper storage conditions; avoid multiple freeze thaw cycles

### **Background**

Small GTP-binding proteins (or GTPases) are a family of proteins that serve as molecular regulators in signaling transduction pathways. Rac, a 21 kDa protein, belongs to the family of Rho GTPases regulating a variety of biological response pathways that include cell motility, cell division, gene transcription, and cell transformation. Like other small GTPases, Rac regulates molecular events by cycling between an inactive GDP-bound form and an active GTP-bound form. In its active (GTP-bound) state, Rac binds specifically to the p21-binding domain (PBD) of p21-activated protein kinase (PAK) to control downstream signaling cascades.

### **Presentation**

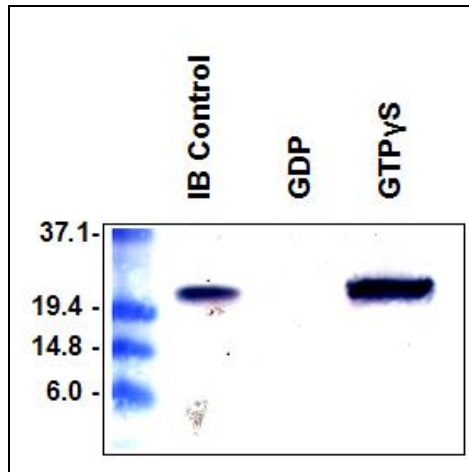
PAK PBD Agarose beads, in color, are easy to visualize, minimizing potential loss during washes and aspirations of Rac-GTP pulldown (Figure 1).



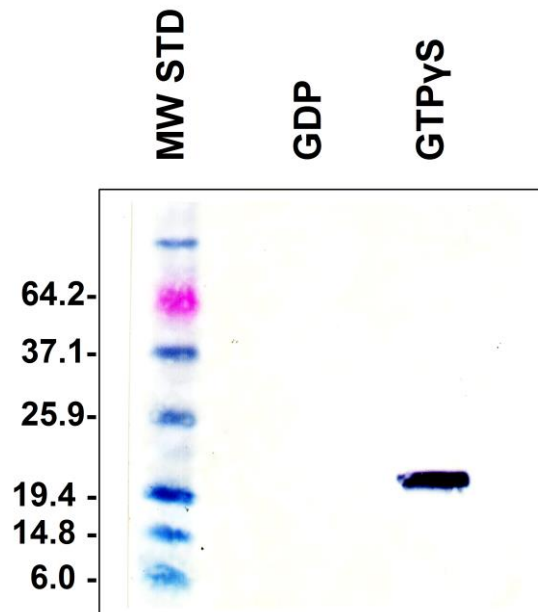
**Figure 1:** PAK-PBD Beads in Color

### Activity

Product specifically interacts and precipitates GTP-bound Rac or Cdc 42 from cell lysate (Figures 2 & 3).



**Figure 2:** Rac Activation Assay. *Lane 1*, GTPase Immunoblot Positive Control. *Lane 2*, 293 cell lysate loaded with GDP and incubated with PAK PBD Agarose beads. *Lane 3*, 293 cell lysate loaded with GTP $\gamma$ S and incubated with PAK-1 PBD Agarose beads. Samples were immunoblotted with anti-Rac antibody.



**Figure 3:** Cdc42 Activation Assay. *Lane 1*, MW Standard. *Lane 2*, 293 cell lysate loaded with GDP and incubated with PAK PBD Agarose beads. *Lane 3*, 293 cell lysate loaded with GTP $\gamma$ S and incubated with PAK-1 PBD Agarose beads. Samples were immunoblotted with anti-Cdc42 antibody.

## References

1. Raftopoulou M., and Hall A. (2004) *Dev Biol.* **265**: 23-32.
2. Bar-Sagi D., and Hall A. (2000) *Cell* **103**: 227-38.
3. Benard, V., Bohl, B. P., and Bokoch, G. M. (1999) *J. Biol. Chem.* **274**, 13198-13204.

## Recent Product Citations

1. Bernadzki, K. M. et al. (2020). Arhgef5 Binds  $\alpha$ -Dystrobrevin 1 and Regulates Neuromuscular Junction Integrity. *Front Mol Neurosci.* doi: 10.3389/fnmol.2020.00104.
2. Speranza, L. et al. (2017). Serotonin 5-HT7 receptor increases the density of dendritic spines and facilitates synaptogenesis in forebrain neurons. *J Neurochem.* doi: 10.1111/jnc.13962.
3. Bijata, M. et al. (2016). Dystroglycan controls dendritic morphogenesis of hippocampal neurons in vitro. *Front Cell Neurosci.* doi:10.3389/fncel.2015.00199.
4. Fusté, N. P. et al. (2016). Cytoplasmic cyclin D1 regulates cell invasion and metastasis through the phosphorylation of paxillin. *Nat Commun.* doi:10.1038/ncomms11581.
5. Alam, J. et al. (2014). N-acetylcysteine and the human serum components that inhibit bacterial invasion of gingival epithelial cells prevent experimental periodontitis in mice. *J Periodontal Implant Sci.* **44**:266-273.
6. Morrison, A. R. et al. (2014). Chemokine-coupled  $\beta$ 2 integrin-induced macrophage Rac2-myosin IIA interaction regulates VEGF-A mRNA stability and arteriogenesis. *J Exp Med.* **211**:1957-1968.
7. Pothula, S. et al. (2013). Regulation of Cdc42 expression and signaling is critical for promoting corneal epithelial wound healing. *Invest. Ophthalmol. Vis. Sci.* **54**: 5343-5352.
8. Cheng, J. et al.(2010).FSP-1 silencing in bone marrow cells suppresses neointima formation in vein graft. *Circ Res.* **110**:230-240.
9. Sabbatini, M. E. et al. (2010). CCK activates RhoA and Rac1 differentially through G-alpha-13 and G-alpha-q in mouse pancreatic acini. *Am. J. Physiol. Cell Physiol.* **298**:C592-C605.
10. Zhang, Q-G. et al. (2009). Estrogen attenuates ischemic oxidative damage via an estrogen receptor alpha-mediated inhibition of NADPH oxidase activation. *J. Neurosci.* **29**:13823-13836.

## Warranty

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS's sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

***This product is for RESEARCH USE ONLY; not for use in diagnostic procedures.***

## **Contact Information**

Cell Biolabs, Inc.  
7758 Arjons Drive  
San Diego, CA 92126  
Worldwide: +1 858-271-6500  
USA Toll-Free: 1-888-CBL-0505  
E-mail: [tech@cellbiolabs.com](mailto:tech@cellbiolabs.com)  
[www.cellbiolabs.com](http://www.cellbiolabs.com)

©2006-2020: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.