

RalBP1 PBD Agarose Beads

CATALOG NUMBER: STA-420

STORAGE: -20°C

QUANTITY AND CONCENTRATION: 800 μ L of 50% Agarose slurry, 400 μ g RalBP1-PBD (amino acid 397-518) 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. Ral, a 24 kDa protein of the Ras superfamily, regulates a variety of biological response pathways that include vesicle trafficking, cytoskeletal rearrangement, and migration. The Ras-like proteins RalA and RalB share 85% identity. Like other small GTPases, Ral regulates molecular events by cycling between an inactive GDP-bound form and an active GTP-bound form. In their active (GTP-bound) state, Ral A and Ral B bind specifically to the protein-binding domain (PBD) of RalBP1 to control downstream signaling cascades.

Presentation

RalBP1 PBD Agarose beads, in color, are easy to visualize, minimizing potential loss during washes and aspirations of Ral-GTP pulldown (Figure 1).

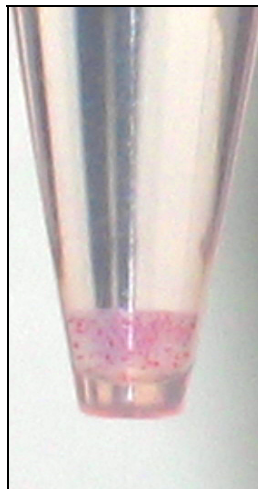


Figure 1: RalBP1-PBD Beads in Color

Activity

Product specifically interacts and precipitates GTP-bound Ral from cell lysate (Figure 2).

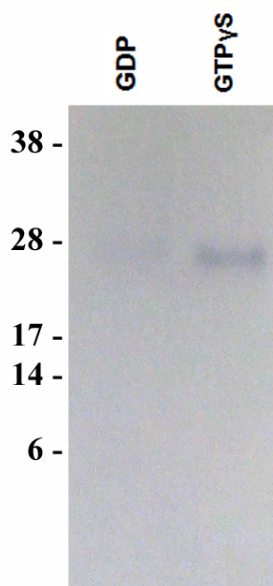


Figure 2: Ral Activation Assay. *Lane 1*, NIH 3T3 cell lysate loaded with GDP and incubated with RalBP1 PBD Agarose beads. *Lane 2*, NIH 3T3 cell lysate loaded with GTP γ S and incubated with RalBP1 PBD Agarose beads.

References

1. Wolthuis RM, Franke B, van Triest M, et al. (1998) *Mol Cell Biol* **18**:2486–2491.
2. Herrmann C, Horn G, Spaargaren M, Wittinghofer A (1996) *J Biol Chem* **271**:6794–6800.
3. Ward Y, Wang W, Woodhouse E, et al. (2001) *Mol Cell Biol* **21**:5958–5969.
4. Oxford G, Owens CR, Titus BJ, et al. (2005) *Cancer Res* **65**:7111–7120.

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
www.cellbiolabs.com

©2009: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.