

Recombinant Human Ral B

CATALOG NUMBER: STA-733 **STORAGE:** -80°C; avoid freeze/thaw cycles

QUANTITY: 10 µg/vial in Tris-HCl, pH 8.0 with 0.1 M NaCl, 1 mM DTT and 10% glycerol

SHELF LIFE: 1 year from date of receipt under proper storage conditions; for longer term storage add 0.1% BSA

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 B regulates molecular events by cycling between an inactive GDP-bound form and an active GTP-bound form. In its active (GTP-bound) state, Ral B binds specifically to the protein-binding domain (PBD) of RalBP1 to control downstream signaling cascades.

Purity and Activity

Cell Biolabs' rRal B is expressed and purified from *E. coli* as a recombinant protein with purity >90% by SDS-PAGE. It has an approximate MW of 25.8 kDa including a His-tag at the N-terminus.

Sequence

MGSSHHHHHHSSGLVPRGSHMGSHMAANKSKGQSSLALHKVIMVGGVGGKSALTQLQFMYDE
FVEDYEPTKADSYRKKVVDLGEVQIDILDTAGQEDYAAIRDNYFRSGEGFLLVFSITEHESFTAT
AEFREQLRVKAEEDKIPLLVGNKSDLEERRQVPVEEARSKAEWGVQYVETSAKTRANVDKV
FFDLMREIRTKKMSNKDKNGKSSKNKKSFKERC

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

©2012: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.