

## Recombinant Human Rap1a

**CATALOG NUMBER:** STA-735      **STORAGE:** -80°C; avoid freeze/thaw cycles

**QUANTITY:** 10 µg/vial in Tris-HCl, pH 7.5 with 0.2 M NaCl, 5 mM DTT and 40% 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. Rap, a 24 kDa protein of the Ras superfamily, regulates a variety of biological response pathways that include cell adhesion, proliferation, differentiation, and apoptosis. The Ras-like proteins Rap1 and Rap2 share 60% identity. Like other small GTPases, Rap1a regulates molecular events by cycling between an inactive GDP-bound form and an active GTP-bound form. In their active (GTP-bound) state, Rap1a binds specifically to the Rap-binding domain (RBD) of RalGDS to control downstream signaling cascades.

### **Purity and Activity**

Cell Biolabs' rRap1a is expressed and purified from *E. coli* as a recombinant protein with purity >85% by SDS-PAGE. It has an approximate MW of 22.8 kDa including a His-tag at the N-terminus.

### **Sequence**

MGSSHHHHHSSGLVPRGSHMREYKLVVLGSGGVGKSALTVQFVQGIFVEKYDPTIEDSYRKQV  
EVDCQQCMLEILDTAGTEQFTAMRDLYMKNQGQGFALVYSITAQSTFNDLQDLREQILRVKDTED  
VPMILVGNKCDLEDERVVGKEQGQNLARQWCNCAFLESSAKSKINVNEIFYDLVRQINRKTTPVE  
KKKPKKKSC

### **References**

1. Kitayama, H., Sugimoto, Y., Matsuzaki, T., Ikawa, Y., and Noda, M. (1989) *Cell* **56**, 77–84.
2. Matsubara, K., Kishida, S., Matsuura, Y., Kitayama, H., Noda, M., and Kikuchi, A. (1999) *Oncogene* **18**, 1303–1312.
3. Bos, J. L., de Rooij, J., and Reedquist, K. A. (2001) *Nature Rev.* **2**, 369–377.
4. Herrmann, C., Horn, G., Spaargaren, M., and Wittinghofer, A. (1996) *J. Biol. Chem.* **271**, 6794–6800.

### **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)

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