

RhoA Expression Vector Set

CATALOG NUMBER: STA-456

STORAGE: -80°C

QUANTITY AND CONCENTRATION: 3 vectors; each 100 µl of bacterial glycerol stock.

Components

1. pcDNA3-RhoA (Wild Type) (Part No. 345601): One tube
2. pcDNA3-RhoA T19N (Dominant Negative) (Part No. 345602): One tube
3. pcDNA3-RhoA G14V (Constitutively Active) (Part No. 345603): One tube

Background

Three members of the Rho family small GTPase, Rho, Rac, and Cdc42, have been shown to play a crucial role in regulating the organization of the actin cytoskeleton in response to extracellular stimuli. Activation of Rho, Rac, and Cdc42 in quiescent Swiss 3T3 fibroblasts induces the assembly of filamentous actin into stress fibers, lamellipodia, and filopodia, respectively. In addition to these effects on the actin cytoskeleton, it has been shown Rac and Cdc42 (and in some cells Rho) can activate JNK and p38 that leads to transcriptional activation. In fibroblast cells, Rho, Rac, and Cdc42 have each been implicated in cell cycle control.

The small GTPase gene from human is cloned in the pCMV5 expression vector (Figure 1).

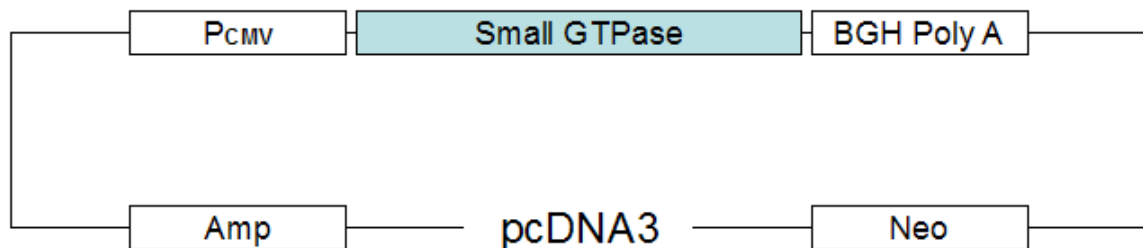


Figure 1. pcDNA3-RhoA Expression Vector Map

References

1. Morgenstern, J. P. and H Land. (1990) *Nuc. Acid Res.* 18, 3587-3596.
2. Schuck S, Manninen A, Honsho M, Fullekrug J and Simons K. (2004) *Proc Natl Acad Sci U S A.* 101, 4912-4917.
3. Machesky L. M. and Hall A. (1996) *Trends Cell Biol.* 6:304-10.

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

©2007-2012: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.