

Anti-ROCK2

Rabbit Polyclonal Antibody

Catalog # R11-63R

Lot # O2121-42

Cited Applications

WB, ELISA, IF, IHC

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the ROCK2 protein and will not cross-react with other ROCK family members

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

ROCK2 antibody was raised against a 15 amino acid synthetic peptide near the center of human ROCK2

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

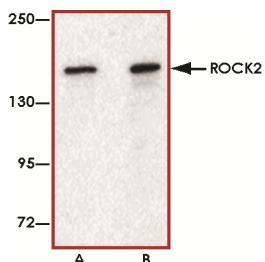
Scientific Background

ROCK2 is a ubiquitously expressed serine/threonine kinase localized in the nucleus that regulates cytokinesis, smooth muscle contraction, the formation of actin stress fibers and focal adhesions, and the activation of the c-fos serum response element (1). ROCK2 is an immediate downstream target of the small GTPase RhoA. ROCK2 may play a pivotal role in cardiovascular diseases such as vasospastic angina, ischemic stroke, and heart failure. Inhibition of ROCKs by statins or other selective inhibitors leads to the upregulation and activation of endothelial nitric oxide synthase (eNOS) and reduction of vascular inflammation and atherosclerosis (2).

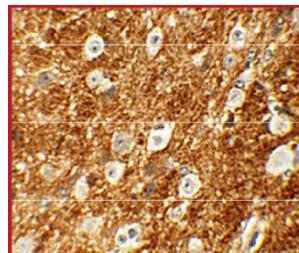
References

- Zhao, Z. et al: Rho-associated kinases play a role in endocardial cell differentiation and migration. *Dev Biol.* 2004 Nov 1;275(1):183-91.
- Noma, K. et al: Physiological role of ROCKs in the cardiovascular system. *Am J Physiol Cell Physiol.* 2006 Mar;290(3):C661-8.

Sample Data



Western blot analysis of ROCK2 in mouse brain tissue lysate with ROCK2 antibody at (A) 1 and (B) 2 ug/ml.



Immunohistochemistry of ROCK2 in mouse brain tissue with ROCK2 antibody at 5 ug/ml.

Anti-ROCK2

Rabbit Polyclonal Antibody

Catalog Number

R11-63R

Specific Lot Number

O2121-42

Purification

Affinity chromatography

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.