

## Anti-ROCK1

Rabbit Polyclonal Antibody

**Catalog # R10-63R**

Lot # O2121-41

### Cited Applications

WB, ELISA, ICC, IF

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

### Specificity

Recognizes the ROCK1 protein and will not cross-react with other ROCK family members

### Cross Reactivity

Human, Mouse and Rat

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

ROCK1 antibody was raised against a 15 amino acid synthetic peptide near the amino terminus of human ROCK1

### Formulation

PBS + 0.02% sodium azide

### Stability

1yr at -20°C from date of shipment

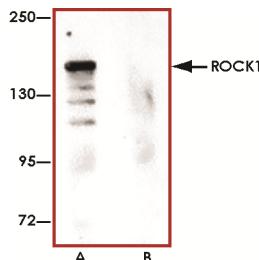
### Scientific Background

ROCK1 is a ubiquitously expressed serine/threonine kinase that is a downstream target of the small GTPase RhoA. ROCK1 is involved in diverse cellular functions, including smooth muscle contraction, actin cytoskeleton organization, cell adhesion and motility, and gene expression (1). ROCK1 contributes to the development of cardiac fibrosis and induction of fibrogenic cytokines in cardiomyocytes in response to pathological stimuli. ROCK1 knockout mice exhibit reduced perivascular and interstitial fibrosis which is associated with reduced expression of a variety of extracellular matrix (ECM) proteins and fibrogenic cytokines (2).

### References

- Zhao, Y M. et al: Rho-associated kinases play a role in endocardial cell differentiation and migration. *Dev Biol.* 2004 Nov 1;275(1):183-91.
- Zhang, C. et al: Targeted deletion of ROCK1 protects the heart against pressure overload by inhibiting reactive fibrosis. *FASEB J.* 2006 May; 20(7):916-25.

### Sample Data



Western blot analysis of ROCK1 in 293 cell lysate with ROCK1 antibody at 1 ug/ml in (A) the absence and (B) the presence of blocking peptide.



Immunocytochemistry of ROCK1 in 293 cells with ROCK1 antibody at 10 ug/ml.

## Anti-ROCK1

Rabbit Polyclonal Antibody

Catalog Number      R10-63R

Specific Lot Number      O2121-41

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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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