

## Anti-PAK1

Rabbit Polyclonal Antibody

**Catalog # P02-63R**

Lot # H283-5

### Cited Applications

- Western blot (1:1000)

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

### Specificity

Recognizes the PAK1 protein

### Cross Reactivity

- Western blot PAK1 from human and mouse cells

PAK1 from other species may also be detectable

### Host

Rabbit

### Immunogen

Synthetic peptide corresponding to amino acids 210-214 of human PAK1

### Formulation

PBS pH 7.4, 0.02% sodium azide, 50% glycerol

### Stability

Store at 4°C (add 0.1% NaN<sub>3</sub>) for several months, and at -20°C for longer periods. 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.

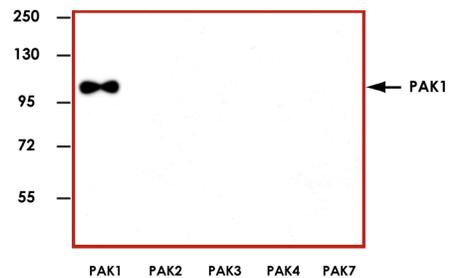
### Scientific Background

PAK1 is a member of the p21-activated kinases (PAKs) which have been implicated in the regulation of cell morphology, motility and transformation. These serine/threonine kinases are activated by and are effectors of small GTPases, CDC42 and RAC. PAK1 belongs to the Group I PAKs which also includes PAK2 and PAK3 (1). PAK1 is a key regulator of the actin cytoskeleton, adhesion and cell motility. Inactive dimeric PAK1 is mainly cytosolic and interaction with the activators Cdc42-GTP and Rac1-GTP stimulates the kinase at the sites of cellular protrusions forming adhesions to the extracellular matrix (2).

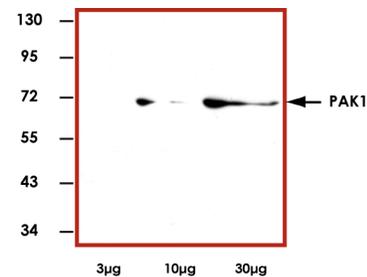
### References

1. Jaffer, Z M. et al: p21-activated kinases: three more join the Pak. Int J Biochem Cell Biol. 2002 Jul;34(7):713-7.
2. Parrini, M C. et al: Spatiotemporal regulation of the Pak1 kinase. Biochem Soc Trans. 2005 Aug;33(Pt 4):646-8.

### Sample Data



Representative western blot with Anti-PAK1 (1:1000) using 20 ng of human recombinant PAK1, PAK2, PAK3, PAK4, and PAK7.



Representative western blot with Anti-PAK1 (1:1000) using 3 µg, 10 µg, and 30 µg of Jurkat cell lysate.

## Anti-PAK1

Rabbit Polyclonal Antibody

Catalog Number

P02-63R

Specific Lot Number

H283-5

Purification

Affinity chromatography

Concentration

1.0 µg/µL

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.**