

## Anti-PAK4

Rabbit Polyclonal Antibody

**Catalog # P05-163R**

Lot # O2109-22

### Cited Applications

WB, ELISA, ICC, IF

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

### Specificity

Recognizes the PAK4 protein

### Cross Reactivity

Human

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

PAK4 antibody was raised against a 13 amino acid synthetic peptide from near the center of human PAK4

### Formulation

PBS + 0.02% sodium azide

### Stability

1yr at -20°C from date of shipment

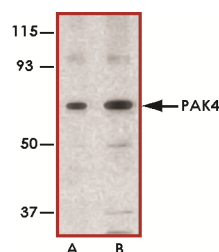
### Scientific Background

PAK4 is a recently identified 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, cdc 42 and Rac. PAK4 belongs to the Group II PAKs which also includes PAK5 and PAK6. PAK4 has been shown to regulate cell morphology and cytoskeletal organization in mammalian cells.

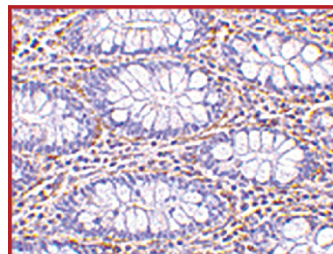
### References

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. Qu, J. et al: Activated PAK4 regulates cell adhesion and anchorage-independent growth. *Mol Cell Biol.* 2001 May;21(10):3523-33.

### Sample Data



Western blot analysis of PAK4 in SW480 lysate with PAK4 antibody at (A) 1 and (B) 2 ug/ml.



Immunohistochemistry of PAK4 in human colon tissue with PAK4 antibody at 10 ug/ml.

## Anti-PAK4

Rabbit Polyclonal Antibody

Catalog Number

P05-163R

Specific Lot Number

O2109-22

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.**