

## Anti-PAK6

Rabbit Polyclonal Antibody

Catalog # P06-63R

Lot # O2109-20

### Cited Applications

WB, ELISA, IF, IHC

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

### Specificity

Recognizes the PAK6 protein

### Cross Reactivity

Human

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

PAK6 antibody was raised against a 13 amino acid synthetic peptide from near the center of human PAK6

### Formulation

PBS + 0.02% sodium azide

### Stability

1yr at -20°C from date of shipment

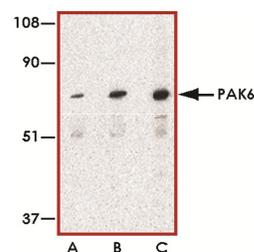
### Scientific Background

PAK6 is a Ser/Thr protein kinase that is a member of the PAK (p21-activated kinase) family that are implicated in the regulation of a number of cellular processes, including cytoskeleton rearrangement, apoptosis and the MAP kinase signaling pathway. PAK6 is highly expressed in the testis and prostate tissues and interacts with the androgen receptor (AR). In response to androgen, PAK6 cotranslocate into the nucleus with AR (1). PAK6 is weakly expressed in normal prostate epithelium but the expression is increased in primary and metastatic prostate cancer cells and is further increased in tumors that relapse after androgen deprivation therapy (2).

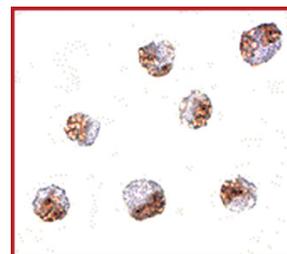
### References

1. Yang F, et al: Androgen receptor specifically interacts with a novel p21-activated kinase, PAK6. J. Biol. Chem. 276: 15345-15353, 2001.
2. Kaur R, et al: Increased PAK6 expression in prostate cancer and identification of PAK6 associated proteins. Prostate. 2008 Oct 1;68(14):1510-6.

### Sample Data



Western blot analysis of PAK6 in Raji lysate with PAK6 antibody at (A) 1, (B) 2, and (C) 4 ug/ml.



Immunocytochemistry of PAK6 in Raji cells with PAK6 antibody at 10 ug/ml.

## Anti-PAK6

Rabbit Polyclonal Antibody

Catalog Number

P06-63R

Specific Lot Number

O2109-20

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.**