



Anti-PAK6

(p21-activated kinase 6)

CATALOG NO.: 54512

BACKGROUND:

The p21-activated kinases (PAKs) are serine-threonine kinases that bind to the active forms of Cdc42 and Rac (1). They are divided into two groups, the first of which include PAK1, 2 and 3, and can be activated by Cdc42/Rac binding. Group 1 PAKs contain an autoinhibitory domain whose activity is regulated by Cdc42/Rac binding. The group 1 PAKs are known to be involved in cellular processes such as gene transcription, apoptosis, and cell morphology and motility. Much less is known about the second group, which includes PAK4, 5 and 6. These proteins are not activated by Cdc42/Rac binding. PAK6 was initially identified as an androgen receptor in a yeast two hybrid screen and was found to be highly expressed in testis and prostate tissues (2). Later experiments have shown it to be activated by MAP kinase kinase 6 and p38 MAP kinase (3), suggesting that PAK6 may play a role in the cellular response to stress-related signals.

SOURCE & REACTIVITY:

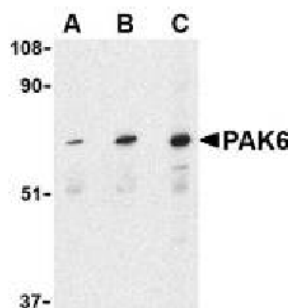
Rabbit polyclonal anti-PAK6 was raised against a 13 amino acid peptide from near the center of human PAK6 (Genbank accession No. NP_064553). Anti-PAK6 is human specific.

APPLICATION:

The following concentration ranges are recommended starting points for this product.

WB: 1.0-2.0 µg/ml

Positive Control: Raji cell lysate



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

Immunocytochemistry of PAK6 in Raji cells with anti-PAK6 at 10 µg/ml.



This product is for in vitro research purposes only.

RELATED PRODUCTS:

Anti-PAK4, Catalog No. **54514**
Anti-PAK5, Catalog No. **54513**

STORAGE:

The antibody is supplied as purified IgG, 50 µg in 250 µl of 1X PBS containing 0.02% sodium azide. Store at 4 °C for up to one year. Avoid repeated freezing and thawing.

REFERENCES:

1. Jaffer ZM, et al (2002) *Int. J. Biochem. Cell Biol.* 34:713-7.
2. Yang F, et al (2001) *J. Biol. Chem.* 276:15345-53.
3. Kaur R, et al (2005) *J. Biol. Chem.* 280:3323-30.