

Anti-phospho-PAK1/2/3 (Thr423)

Rabbit Polyclonal Antibody

Catalog # P02-653R

Lot # B3216-70

Cited Applications

ELISA, WB

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

Specificity

Recognizes the PAK (P21-GTPase Activated Protein Kinase)

Cross Reactivity

Human

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

The antibody was produced against human PAK1/2/3 phospho peptide corresponding to a region in the protein conjugated to KLH.

Formulation

0.02 M Potassium Phosphate, 0.15 M Sodium Chloride, pH 7.2 + 0.01% (w/v) Sodium Azide.

Stability

1yr at -20°C from date of shipment

Scientific Background

In mammals, there are several identified isoforms of p21-activated protein kinases or PAKs. PAKs 1, 2 and 3 are members of the serine/threonine p21-activating kinases that serve as targets for the small GTP binding proteins Cdc42 and RAC1. They have been implicated in wide range of biological activities, which include cytoskeletal reorganization and nuclear signaling. CDC42 and RAC1 induce autophosphorylation of PAK2, which stimulates sustained phosphorylation of other substrates [2]. Autophosphorylation of Thr402 is essential for activation of PAK (Jakobi et al., 2000).

References:

1. Martin, G. A. et al: A novel serine kinase activated by rac1/CDC42Hs-dependent autophosphorylation is related to PAK65 and STE20. EMBO J. 14: 1970-1978, 1995.
2. Jakobi R. et al: Substrates enhance autophosphorylation and activation of p21-activated protein kinase gamma-PAK in the absence of activation loop phosphorylation. Eur J Biochem 267:4414-4421, 2000.

Anti-phospho-PAK1/2/3 (Thr423)

Rabbit Polyclonal Antibody

Catalog #	P02-653R
Lot #	B3216-70
Purification	Immunoaffinity 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
www.signalchem.com

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