

Recombinant PAK2 protein

Catalog No: 31188

Expressed In: Baculovirus

Quantity: 10 µg

Source: Human

Buffer Contents: 10 µg of Recombinant PAK2 protein in 50 mM Tris-HCl, pH 8.0; 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20% glycerol. Protein is supplied at 0.403 µg/µl.

Background: The p21 activated kinases (PAK) are critical effectors that link Rho GTPases to cytoskeleton reorganization and nuclear signaling. The PAK proteins are a family of serine/threonine kinases that serve as targets for the small GTP binding proteins, CDC42 and RAC1, and have been implicated in a wide range of biological activities. The protein encoded by this gene is activated by proteolytic cleavage during caspase-mediated apoptosis, and may play a role in regulating the apoptotic events in the dying cell.

Protein Details: Recombinant human PAK2 protein was produced using baculovirus infected Sf9 cells. The protein was made against amino acids D3-R534, accession number NM_002577 and N-terminally fused to GST-HIS₆-Thrombin cleavage site. Purified by GSH-agarose affinity purification.

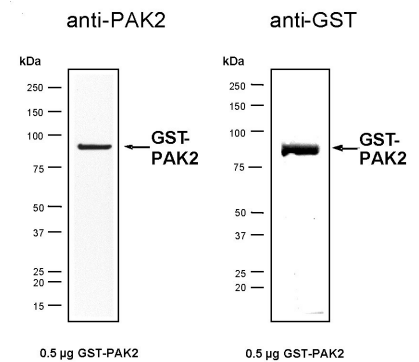
Application Notes: Recombinant PAK2 is suitable for kinase assays and Western blot. The molecular weight of the protein is ~87.954 kDa. The activity of the protein is ~ 228 pmol/µg min.

Recommended kinase reaction conditions: 60 mM HEPES-NaOH, pH 7.5, 3 mM MgCl₂, 3 mM MnCl₂, 3 µM Na-orthovanadate, 1.2 mM DTT, ATP (variable), 2.5 µg/50 µl PEG20.000, Substrate: tetra (LRRWSLG), 5 µg / 50 µl, Recombinant PAK2 : 50 ng/50 µl.

Kinase activity may vary depending on the substrate and reaction conditions used.

Storage and Guarantee: This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



Recombinant PAK2 protein tested by SDS-PAGE gel.

0.5 µg of PAK2 protein was analyzed using SDS/PAGE followed by Western blot with either an anti-PAK2 or anti-GST antibody.