

PAK-GST Protein
(human p21 activated kinase PBD)
Cat. # PAK01

Upon arrival store at 4°C (desiccated)
See datasheet for storage after reconstitution

Material

The Rac/Cdc42 binding domain of the human p21 activated kinase 1 protein (PAK) has been overexpressed as a GST-tagged recombinant protein in a bacterial expression system. The recombinant protein (amino acids 67-150) includes the highly conserved PBD region (also referred to as the CRIB region) and sequences required for the high affinity interaction with GTP-Rac and GTP-Cdc42 proteins. The recombinant protein is tagged with GST (28 kDa) at its amino terminus and has an approximate molecular weight of 34 kDa. PAK-GST protein (250 µg protein per tube) is supplied as a white lyophilized powder.

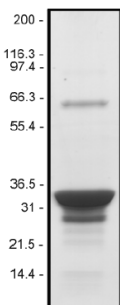
Storage and Reconstitution

Briefly centrifuge to collect the product at the bottom of the tube. The protein should be reconstituted to 3 mg/ml by the addition of 83 µl of distilled water. When reconstituted, the protein will be in the following buffer: 50 mM Tris pH 7.5, 50 mM NaCl, 1.0% (w/v) dextran and 5% (w/v) sucrose. For storage, the reconstituted protein should be aliquoted into experiment sized amounts, snap frozen in liquid nitrogen and stored at -70°C. Under these conditions the protein is stable for 6 months. **To maintain high biological activity, the protein should not be exposed to repeated freeze thaw cycles.** The lyophilized protein is stable for 6 months if stored desiccated to <10% humidity at 4°C.

Purity

Protein purity is determined by scanning densitometry of Coomassie Blue stained protein on a 12% SDS polyacrylamide gel. PAK-GST protein was determined to be 88% pure (see Figure 1).

Figure 1. PAK-GST Protein Purity Determination. A 20 µg sample of PAK-GST protein (molecular weight approx. 34 kDa) was separated by electrophoresis in a 12% SDS-PAGE system and stained with Coomassie Blue. Protein quantitation was performed using the Precision Red Protein Assay Reagent (Cat. # ADV02). Mark12™ molecular weight markers are from Invitrogen.



Biological Activity Assay

PAK-GST protein specifically recognizes and binds the active "GTP-bound" forms of the Rac and Cdc42 proteins (1). It has a much lower affinity for the inactive "GDP-bound" forms of Rac and Cdc42. When coupled to a colored glutathione sepharose matrix, the PAK-GST protein beads become a convenient tool for assaying the activity of the Rac and Cdc42 proteins. A standard biological assay for PAK-GST protein consists of a Rac protein pulldown from human platelet extracts loaded with either GTPγS or GDP.

Reagents

1. PAK-GST protein (Cat. # PAK01)
2. Loading buffer (150 mM EDTA)
3. Stop buffer (600 mM MgCl₂)
4. Wash buffer (25 mM Tris pH 7.5, 30 mM MgCl₂, 40 mM NaCl)
5. Cell lysis buffer (50 mM Tris pH 7.5, 10 mM MgCl₂, 0.3 M NaCl, 2% IGEPAL)
6. GTPγS (20 mM solution)
7. GDP (100 mM solution)
8. Human platelet extract prepared in 50 mM PIPES pH 7.0, 130 mM NaCl, 1 mM PMSF, 1 mM DTT, 5 µg/ml leupeptin, 0.5% Triton X-100
9. Protease inhibitor cocktail (Cat. # PIC02)
10. Anti-Rac polyclonal antibody (Cat. # ARC03)
11. Glutathione beads

Equipment

1. Microfuge at 4°C
2. SDS-PAGE and Western blot apparatus

Method

1. Dilute the platelet extract to 1 mg/ml with cell lysis buffer containing protease inhibitors.
2. Centrifuge the extract at 14k rpm at 4°C for 5 min to pellet cell membranes and insoluble material.
3. Aliquot 500 µg (500 µl) of the extract into two experimental tubes.
4. Add 1/10th the volume of loading buffer to each tube (final conc. 15 mM EDTA).
5. Add 1/100th the volume of GDP to one tube (final conc. 1.0 mM GDP).
6. Add 1/100th the volume of GTPγS to the other tube (final conc. 0.2 mM GTPγS).
7. Incubate both tubes at room temperature for 15 min.
8. Stop the reaction by adding 1/10th the volume of stop buffer to each tube (final conc. 60 mM MgCl₂).
9. Resuspend PAK-GST protein to 3 mg/ml and add 20 µg (6.7 µl) protein to each reaction tube.

10. Gently rotate the tubes at 4°C for 30 minutes.
11. Add 10 μ l of glutathione beads (equilibrated in cell lysis buffer) to each reaction and incubate on the rotator for a further 30 minutes at 4°C.
12. Centrifuge the beads at 8k rpm at 4°C for 1 min.
13. Remove the supernatant and wash the beads in 500 μ l of wash buffer.
14. Pellet the beads and resuspend in 20 μ l of SDS sample buffer.
15. The protein samples can now be analyzed by Western blot procedure using a Rac1 monoclonal antibody (Cat. # ARC03). NOTE: The PAK01 can also be used in a Cdc42 activation assay using anti-Cdc42 monoclonal antibody (Cat. # ACD03).
16. Typical assay results are shown in Figure 2.

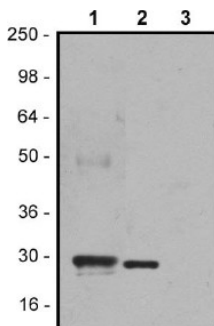


Figure 2. Selective Binding of PAK-GST Protein to the GTP-bound Form of Rac1 *In Vitro*. Human platelet extracts (500 μ g) were loaded with either GTPyS (lane 2), or GDP (lane 3) as described in the method. The extracts were then incubated with 20 μ g of PAK-GST protein, subsequently the PAK-GST was captured by addition of glutathione beads. The protein-bead complexes were recovered by centrifugation and subjected to Western blot analysis using a Rac1 specific polyclonal antibody. Lane 1 shows 50 ng of recombinant Rac-His control protein (Note: His-Rac1 runs slightly higher than endogenous Rac1 due to the presence of the 6x-His tag). SeeBlue™ molecular weight markers are from Invitrogen.

Product Uses

- Measurement of the GTP/GDP ratio of Rac or Cdc42 *in vitro*.
- Quantitation of GTP-Rac/Cdc42 from tissue and tissue culture cell lysates.

References

1. Manser, E. et al. 1994. *Nature*. **367**: 40-46.

Product Citations/Related Products

For the latest citations and related products please visit www.cytoskeleton.com.