

His-Rac1 L61 Mutant Protein Constitutively Active Cat. # R6101

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

Material

The constitutively active form of the human Rac1 protein has been produced in a bacterial expression system. The protein has a glutamine to leucine substitution at amino acid 61, creating a constitutively active mutant protein that will not hydrolyze GTP. The recombinant protein contains six histidine residues (His-tag) at its amino terminus. The approximate molecular weight of His-L61 Rac1 protein is 28 kDa. His-Rac1 L61 protein is supplied as a white lyophilized powder.

Storage and Reconstitution

Before reconstitution, briefly centrifuge to collect the product to the bottom of the tube. The protein should be reconstituted to 1 mg/ml by the addition of 10 μ l of distilled water. When reconstituted, the protein will be in the following buffer: 10 mM Tris pH 7.5, 10 mM NaCl, 0.1 mM MgCl₂, 0.5% (w/v) sucrose, and 0.1% (w/v) dextran. In order to maintain high biological activity of the protein, it is strongly recommended that the protein solution be supplemented with DTT to 1 mM final concentration, aliquoted into "experiment sized" amounts, snap frozen in liquid nitrogen and stored at -70°C. The protein is stable for six months under these conditions. The protein must not be exposed to repeated freeze-thaw cycles. The lyophilized protein is stable at 4°C desiccated (<10% humidity) for six months.

Purity

Protein purity is determined by scanning densitometry of Coomassie Blue stained protein on a 12% polyacrylamide gel. His-Rac1 L61 protein was determined to be 90% pure (see Figure 1). The higher molecular weight band in Fig. 1 is a doublet of R6101.

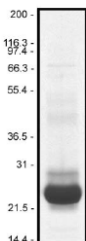


Figure 1. His-Rac1 L61 Protein Purity Determination. A 10 μ g sample of recombinant His-Rac1 L61 protein (molecular weight approx. 28 kDa) was separated by electrophoresis in a 12% SDS-PAGE system, and stained with Coomassie Blue. Protein quantitation was performed using the Precision Red Protein Assay Reagent (Cat. # ADV02). Mark12 molecular weight markers are from Invitrogen.

Biological Activity Assay

His-Rac1 L61 mutant protein can bind GTP but its intrinsic GTPase activity has been eliminated, resulting in a constitutively active protein. A standard biological assay for His-Rac1 L61 activity consists of a pull-down assay using PAK-PBD beads (Cat. # PAK02). The PAK (p21 Activated Kinase CRIB domain) protein is an effector of Rac1, and will specifically bind to active GTP bound Rac1. Stringent quality control ensures that >80% of His-Rac1 L61 protein can be pulled down using this method.

Reagents

1. Recombinant His-Rac1 L61 constitutively active protein (Cat. # R6101)
2. Recombinant His-Rac1 wild-type protein (Cat. # RC01)
3. PAK-PBD beads (Cat. # PAK02)
4. Cell lysis buffer (50 mM Tris pH 7.5, 10 mM MgCl₂, 0.3 M NaCl, 2% IGEPAAL)
5. Loading buffer (150 mM EDTA)
6. Stop buffer (600 mM MgCl₂)
7. Wash buffer (25 mM Tris pH 7.5, 30 mM MgCl₂, 40 mM NaCl)
8. GTPyS (20 mM solution) (Cat. # BS01)
9. GDP (100 mM solution)
10. BSA (10 mg/ml)
11. Anti-Rac1 monoclonal antibody (Cat. # ARC03)

Equipment

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

Method

1. Dilute His-Rac1 L61 constitutively active protein to 0.1 μ g/ μ l with cold cell lysis buffer.
2. Dilute His-Rac1 wild-type protein to 0.1 μ g/ μ l with cold cell lysis buffer.
3. Resuspend PAK-PBD beads to 1 μ g/ μ l by the addition of 500 μ l cold distilled water.
4. Add 23 μ l of cell lysis buffer and 2 μ l of loading buffer into two microfuge tubes on ice.
5. Add 2 μ l (200 ng) of His-Rac1 wild-type protein into both tubes.
6. Add 3 μ l of GTPyS to one tube and 3 μ l of GDP to the other tube. Incubate the loading reactions at room temperature for 15 min.
7. Repeat the nucleotide loading steps 4 through 6 with His-Rac1 L61 constitutively active protein.
8. Stop all reactions with the addition of 4 μ l of stop buffer and place on ice.
9. Add 215 μ l of cold cell lysis buffer and 20 μ l of BSA to each

10. tube on ice.
11. Add 10 μ l (10 μ g) of PAK-PBD beads to each tube and rotate for 30 min at 4°C
12. Pellet the beads at 6800 x g in a microfuge 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 as before and resuspend in 20 μ l of SDS sample buffer.
15. The bead and supernatant samples can now be analyzed by Western blot using a Rac1 specific monoclonal antibody (Cat. # ARC03).
16. Typical assay results are shown in Figure 2.

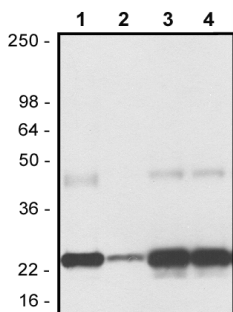


Figure 2. Binding of Wild-type and Constitutively active His-Rac1 to PAK-PBD beads *in vitro*. 200 ng of wild-type and constitutively active His-Rac1 protein were loaded with either GTP γ S (lanes 1 and 3) or GDP (lanes 2 and 4), subjected to a pull-down assay with 10 μ g of PAK-PBD beads and analyzed by Western blot using a Rac1 specific monoclonal antibody as described in the method. Lanes 1 and 2, wild-type His-Rac1. Lanes 3 and 4, constitutively active His-Rac1. Note: GTP γ S and GDP samples look identical in the constitutively active His-Rac1 pull-down due to extremely poor nucleotide exchange. See Blue molecular weight markers are from Invitrogen.

Product Uses

- Identification of Rac1 binding proteins
- Study of Rac1 function *in vivo* by the introduction of constitutively active His-Rac1 into live cells

Product Citations/Related Products

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