

**His-Cdc42 L61 Mutant Protein
Constitutively Active
Cat. # C6101**

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

Material

The constitutively active form of the human Cdc42 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-Cdc42 L61 protein is 25 kDa. His-Cdc42 L61 protein is supplied as a lyophilized white powder.

Storage and Reconstitution

The protein should be reconstituted to 1 mg/ml by the addition of 10 µl of distilled water. The protein will be in the following buffer; 10 mM Tris pH 7.5, 10 mM NaCl, 0.1 mM MgCl₂, 0.5% sucrose and 0.1% 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 6 months under these conditions. The protein must not be exposed to repeated freeze-thaw cycles. The lyophilized protein is stable at 4°C for 1 year.

Purity

Protein purity is determined by scanning densitometry of Coomassie Blue stained protein on a 12% polyacrylamide gel. His-Cdc42 L61 protein was determined to be 70% pure (see Figure 1).

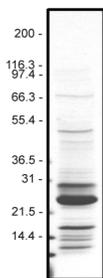


Figure 1. His-Cdc42 L61 Protein Purity Determination. A 10 µg sample of recombinant His-Cdc42 L61 protein (molecular weight approx. 25 kDa) was separated by electrophoresis in a 12% SDS-PAGE system, and stained with Coomassie Blue. Protein quantitation was performed using the Precision Red Protein Assay Reagent (Cat. # ADV02). Mark12 molecular weight markers are from Invitrogen.

Biological Activity Assay

His-Cdc42 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-Cdc42 L61 activity consists of a pulldown assay using PAK-PBD beads (Cat. # PAK02). The PAK (p21 Activated Kinase CRIB domain) protein is an effector of Cdc42 and will specifically bind to active GTP-Cdc42. Stringent quality control ensures that > 80% of His-Cdc42 L61 protein can be pulled down using this method.

Reagents

1. Recombinant His-Cdc42 L61 constitutively active protein (Cat. # C6101)
2. Recombinant His-Cdc42 wild-type protein (Cat. # CD01)
3. PAK-PBD beads (Cat. # PAK02)
4. Loading buffer (150 mM EDTA)
5. Stop buffer (600 mM MgCl₂)
6. Wash buffer (25 mM Tris pH 7.5, 30 mM MgCl₂, 40 mM NaCl)
7. Cell lysis buffer (50 mM Tris pH 7.5, 10 mM MgCl₂, 0.3M NaCl, 2% IGEPAL)
8. GTPγS (20 mM solution) (Cat. # BS01)
9. GDP (100 mM solution)
10. BSA (10 mg/ml)
11. Anti-Cdc42 polyclonal antibody (Cat. # ACD02)

Equipment

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

Method

1. Dilute His-Cdc42 L61 constitutively active protein to 0.1 µg/µl with cold Cell lysis buffer.
2. Dilute His-Cdc42 wild-type protein to 0.1 µg/µl with cold Cell lysis buffer.
3. Resuspend PAK-PBD beads to 1 µg/µl by the addition of 500 µl distilled water.
4. Add 23 µl of Cell lysis buffer and 2 µl of Loading buffer into two microfuge tubes on ice.
5. Add 2 µl (200 ng) of His-Cdc42 wild-type protein into both tubes.
6. Add 3 µl of GTPγS to one tube and 3 µl 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-Cdc42 L61 constitutively active protein.
8. Stop all reactions with the addition of 4 µl of Stop buffer and place on ice.
9. Add 215 µl of cold Cell lysis buffer and 20 µl of BSA to each tube on ice.

10. Add 10 μ l (10 μ g) of PAK-PBD beads to each tube and rotate for 30 min at 4°C
11. Pellet the beads at 8k rpm in a microfuge at 4°C for 1 min.
12. Remove the supernatant and wash the beads in 500 μ l of Wash buffer.
13. Pellet the beads as before and resuspend in 20 μ l of SDS sample buffer.
14. The bead and supernatant samples can now be analyzed by Western blot using a Cdc42 specific polyclonal antibody (Cat. # ACD02).
15. Typical assay results are shown in Figure 2.

Product Uses

- Identification of Cdc42 binding proteins
- Study of Cdc42 function *in vivo* by the introduction of constitutively active His-Cdc42 into live cells

Product Citations/Related Products

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

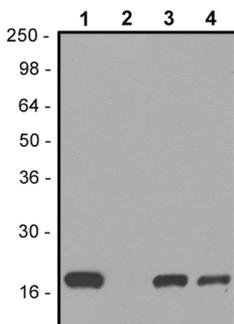


Figure 2. Binding of Wild-type and Constitutively active His-Cdc42 to PAK-PBD beads *in vitro*. 200 ng of wild-type and constitutively active His-Cdc42 protein were loaded with either GTP γ S (lanes 1 and 3) or GDP (lanes 2 and 4), subjected to a pull-down assay with 10 μ g of PAK-PBD beads and analyzed by Western blot using a Cdc42 specific polyclonal antibody as described in the method. Lanes 1 and 2, wild-type His-Cdc42. Lanes 3 and 4, constitutively active His-Cdc42. Note: GTP γ S and GDP samples look identical in the constitutively active His-Cdc42 pull down due to extremely poor nucleotide exchange. SeeBlue molecular weight markers are from Invitrogen.