

**Cdc42 N17-GST Protein**  
**Dominant Negative**  
**Cat. # C17G01**

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

#### Material

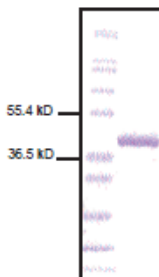
The dominant negative form of the human Cdc42 protein has been produced in a bacterial expression system. This protein has a threonine to asparagine substitution at amino acid 17. The recombinant protein is tagged with GST (28 kDa) at its amino terminus. The approximate molecular weight of the Cdc42 N17-GST protein is 45 kDa. Cdc42 N17-GST 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 1 mg/ml by the addition of 25  $\mu$ l of distilled water. The protein will then be in the following buffer: 10 mM Tris pH 7.5, 10 mM NaCl, 0.3 mM  $MgCl_2$ , 1.0% sucrose, 0.2% 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  $-70^\circ C$ . The protein is stable for 6 months if stored at  $-70^\circ C$ . The protein should not be exposed to repeated freeze-thaw cycles. The lyophilized protein is stable for 1 year if stored desiccated to <10% humidity at  $4^\circ C$ .

#### Purity

Protein purity is determined by scanning densitometry of Coomassie Blue stained protein on a 12% polyacrylamide gel. Cdc42 N17-GST protein was determined to be >90% pure (see Figure 1).



**Figure 1. Cdc42 N17-GST Protein Purity Determination.** A 10  $\mu$ g sample of recombinant Cdc42 N17-GST protein (molecular weight approx. 45 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.

#### Product Citations/Related Products

For the latest citations and related products please visit [www.cytoskeleton.com](http://www.cytoskeleton.com).