



The Protein  
Experts

Cytoskeleton, Inc

# Datasheet

v. 1.0

### **p50Rho GAP protein: Catalytic domain (human recombinant)**

Cat. # GAS01

**Upon arrival store at 4°C (desiccated)**

**See datasheet for storage after reconstitution**

## Material

The catalytic domain of human p50RhoGAP protein has been produced in a bacterial expression system. The protein consists of the RhoGAP domain of p50 RhoGAP (amino acids 198-439, approx. 29 kDa) and an amino terminal GST protein tag (approx. 28 kDa). The protein is supplied as a lyophilized powder. Each tube contains 50 µg of protein.

#### **Storage and Reconstitution**

The protein should be reconstituted to 1 mg/ml by the addition of 50  $\mu$ l of distilled water. The protein will be in the following buffer; 2 mM Tris pH 7.5, 0.5 mM MgCl<sub>2</sub>, 0.5% sucrose, 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" aliquots and snap frozen in liquid nitrogen. The protein can be stored a -70°C for 6 months. The protein must not be exposed to repeated freeze-thaw cycles. The lyophilized protein is stable at 4°C or -70°C for 1 year when stored dessicated (<10% humidity).

## Purity

Protein purity is determined by scanning densitometry of Coomassie blue stained protein on a 4-20% polyacrylamide gradient gel. The protein was determined to be >85% pure. (see Figure 1).

**Figure 1.** p50RhoGAP catalytic domain Protein Purity Determination. A 10 ug sample of recombinant GAS01 protein (molecular weight approx. 57 kD) was separated by electrophoresis in a 4-20% SDS-PAGE system. The protein was stained with coomassie blue. Protein quantitation was determined using the Precision Red Protein Assay Reagent (Cat.# ADV02).



## Biological Activity Assay

Biological activity of p50 RhoGAP catalytic domain protein is determined by the ability of the protein to catalyse GTP hydrolysis by RhoA and Rac1 proteins.

## Reagents

1. Recombinant p50 RhoGAP (Cat.# GAS01)
  2. Recombinant RhoA-His protein (Cat.# RH01)
  3. Recombinant Rac1-His protein (Cat.# RC01)
  4. GTP stock (100 mM) (Cat.# BST06-001)
  5. 2X reaction buffer (80 mM PIPES pH 7.5, 20 mM EDTA, 200 mM NaCl)
  6. CytoPhos reagent (Cat.# BK054)

All reagents are available in biochemical kit Cat # BK105

## Equipment

1. Spectrophotometer set to 650 nm
  2. Corning 96-well half area plates (Cat # 3686) or other plates with low protein binding surface.

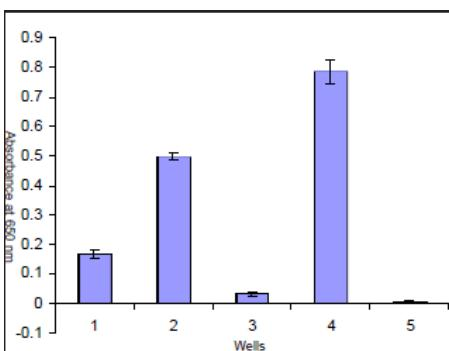
## Method

1. Resuspend all protein reagents in distilled water to give 1 mg/ml final concentration.
  2. Add 15  $\mu$ l 2X reaction buffer to each of 5 wells in a half area microtiter plate.
  3. Add the following to each well; well # 1 = 5  $\mu$ l Rac1 protein, well # 2 = 5  $\mu$ l Rac1 protein plus 5  $\mu$ l of GAS01, well # 3 = 5  $\mu$ l RhoA protein, well # 4 = 5  $\mu$ l RhoA protein plus 5  $\mu$ l of GAS01, well # 5 = 5  $\mu$ l GAS01.
  4. Make up the volume of each reaction to 30  $\mu$ l using distilled water.
  5. Initiate all reactions by adding GTP to 200  $\mu$ M final concentration and incubate at 37°C for 20 minutes.
  6. Stop the reaction by adding 120  $\mu$ l of CytoPhos reagent into each well.
  7. Allow the green color to develop for 10 minutes.
  8. Read reactions at 650 nm. An example of standard results are given in Figure 2, increased OD650 readings indicate enhanced GTP hydrolysis in those samples.

NOTE: The assay as described here is a qualitative analysis of GAP activity. Specific GAP activity in nmoles/min/mg of small G-protein can be determined by taking timepoints from a scaled up GAP reaction. Readings can be quantitated against a standard phosphate curve. The standard curve can be performed as follows:

- a. Prepare a 0.1 mM KH<sub>2</sub>PO<sub>4</sub> solution in distilled water and add 1, 2, 5 and 20 µl to individual wells of a half area plate.

- b. Bring the volumes to 30  $\mu$ l with distilled water and add 120  $\mu$ l of CytoPhos reagent into each well.
- c. Proceed with color development as described above.

**Figure 2.**
**Rac1 & RhoA GAP Assay**


Rac1 (wells 1 & 2) and RhoA (wells 3 & 4) proteins were incubated at 37°C in the absence (wells 1 & 3) or presence (wells 2 & 4) of p50RhoGAP catalytic domain protein. Each 30  $\mu$ l reaction contains 5  $\mu$ g of either Rac1 or RhoA protein and +/- 5  $\mu$ g of p50RhoGAP catalytic domain protein. Well 5 contains p50RhoGAP catalytic domain protein only. All reactions were incubated at 37°C for 20 minutes followed by the addition of 120  $\mu$ l of CytoPhos reagent. After a 10 minute incubation the reactions were read at 650nm. Increased OD650 indicates enhanced GTP hydrolysis.

**Product Uses**

- Study of GAP activity with different GTPases.
- Identification of GAP binding proteins
- Positive control for other RhoGAP studies.

**References**

1. J. Biol. Chem. (1994) 269: 1137-1142.

**Product Citations/Related Products**

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