



## His-Rac2 Protein

### Wild-type

(Human recombinant)

Cat. # RC02

Lot # 001 Amount 1 x 100 µg

Store at 4°C (desiccated) or at -70°C when reconstituted

### Material

Wild-type human Rac2 protein has been produced in a bacterial expression system. The recombinant protein contains six histidine residues (His-tag) at its amino terminus. The molecular weight of the His-Rac2 is approximately 22 kDa. His-Rac2 protein is supplied as a white lyophilized powder.

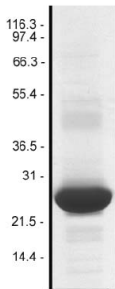
### Storage and Reconstitution

Before reconstitution, briefly centrifuge to collect the product at the bottom of the tube. The protein should be reconstituted to 5 mg/ml with the addition of Milli-Q water. When reconstituted, the protein will be in the following buffer: 50 mM Tris pH 7.5, 0.5 mM MgCl<sub>2</sub>, 50 mM NaCl, 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 if stored at -70°C. **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 gradient gel. His-Rac2 protein was determined to be >90% pure. (see Figure 1).

**Figure 1. His-Rac2 Protein Purity Determination.** A 20 µg sample of recombinant His-Rac2 protein (molecular weight approx. 22 kDa) was separated by electrophoresis in a 4-20% SDS-PAGE system and stained with Coomassie Blue. Protein quantitation was determined using the Precision Red Protein Assay Reagent (Cat. # ADV02). Mark12 molecular weight markers are from Invitrogen.



## Biological Activity Assay

The biological activity of His-Rac2 can be determined from its ability to catalyze the hydrolysis of GTP (GTPase activity). The RhoGAP Assay Biochem Kit (Cat. # BK105) is used to monitor GTP hydrolysis by His-Rac2 in the presence of p50 RhoGAP. Stringent quality control ensures that the hydrolysis of GTP by His-Rac2 is enhanced three fold in the presence of the p50 RhoGAP.

### Reagents

- 1) Recombinant His-Rac2 protein (Cat. # RC02)
- 2) RhoGAP Assay Biochem Kit (Cat. # BK105)

### Equipment

- 1) Microplate spectrometer capable of reading at 650 nm. Cytoskeleton Inc. recommends the SpectroMax250 from Molecular Devices Inc.
- 2) Corning 96-well half area plate (Cat. # 3696) or other plate with low protein binding surface.

### Method

- 1) Resuspend the His-Rac2 protein as described in the reconstitution section for a 250 µM solution. Dilute one aliquot to 50 µM with cold Milli-Q water. Keep on ice.
- 2) Prepare two reaction mixes according to the RhoGAP Assay Biochem Kit. Reaction mix #1 contains p50 RhoGAP (32 µg/100 µl) and will be used to monitor p50 RhoGAP activated GTPase activity. Reaction mix #2 contain no GAP protein and will be used to monitor intrinsic GTPase activity.
- 3) Prepare and dilute a GTP stock solution to 800 µM in cold Milli-Q water. Keep on ice.
- 4) Add the following proteins into duplicate wells of a 96 well plate on ice:

Wells A1 and A2: 5 µl of 50 µM His-Rac2

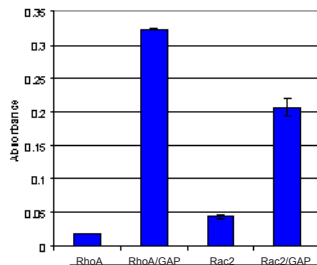
Wells B1 and B2: 5 µl of 50 µM His RhoA

Wells C1 and C2: 5 µl of 50 µM His-Rac2

Wells D1 and D2: 5 µl of 50 µM His RhoA

- 5) Pipette 25 µl of reaction mix #1 into wells A1, A1, B1 and B2.
- 6) Pipette 25 µl of reaction mix #2 into wells C1, C2, D1 and D2.
- 7) Using a multichannel pipette, add 10 µl of 800 µM GTP to each well and incubate at 37°C for 20 min. Shake the plate for 5 s to ensure complete solution mixing.
- 8) After 20 min, remove the plate and add 120 µl of Cytophos reagent (included in BK105) to each well and incubate at room temperature for 10 min.
- 9) Read the absorbance at 650 nm. A typical GAP assay result is shown in Figure 2.

**Figure 2. GTPase Activity of Recombinant Rac2 and RhoA.** Recombinant His-Rac2 and His-RhoA were assayed for GTPase activity using the RhoGAP Assay Biochem Kit (Cat. # BK105) as described. Each reaction contains +/- 5 µg His-Rac2, +/- 5 µg His-RhoA, +/- 8 µg p50 RhoGAP and 200 µM GTP. Reactions were incubated at 37°C for 20 min followed by the addition of Cytophos reagent for 10 min to determine the phosphate generated by the hydrolysis of GTP. His-Rac2 shows a three fold increase in GTP hydrolysis in the presence of p50 RhoGAP. His-RhoA activity is shown for comparison.



### Product Uses

- \* Control for the measurement of the GTP/GDP ratio of Rac2 *in vitro*
- \* Identification of Rac2 binding proteins.
- \* Study of Rac2 function *in vivo* by the introduction of His-Rac2 into live cells
- \* Quantitation standard for activated Rac2 in tissue culture lysates

### Related Products

Cytoskeleton Inc. is the leading supplier of purified small G-proteins, visit our web site or call for information on the small G-proteins currently available. These include the small G-protein Activation Assay Kits, and a variety of affinity reagents for small G-protein activation assays:

- |   |               |
|---|---------------|
| * RhoA Activation Assay, ELISA format, Luminescence | Cat. # BK121  |
| * RhoA Activation Assay, ELISA format, Absorbance   | Cat. # BK124  |
| * Rac Activation Assay Kit                          | Cat. # BK035  |
| * Cdc42 Activation Assay Kit                        | Cat. # BK034  |
| * RhoA Activation Assay Kit                         | Cat. # BK036  |
|   |               |
| * EasyRad Phosphate Assay Kit                       | Cat. # BK055  |
| * GEF Assay Kit                                     | Cat. # BK100  |
| * GAP Assay Kit                                     | Cat. # BK105  |
| * His-Rac1 protein: constitutively active           | Cat. # R6101  |
|   |               |
| * GST-Rac1 protein: wild-type                       | Cat. # RCG01  |
|   |               |
| * GST-Rac1 protein: constitutively active           | Cat. # R61G01 |
|   |               |
| * GST-Rac1 protein: dominant negative               | Cat. # R17G01 |
| * PAK-1 PBD beads                                   | Cat. # PAK02  |