

## R-Ras Protein: wild-type

(Human recombinant)

Cat. # CS-RS05

Lot # 012 1 x 300 µg

Upon arrival store at 4°C (desiccated)

See datasheet for storage after reconstitution

### Material

The wild-type human Ras-related (R-Ras) protein (Spaargaren et al. 1994) has been produced in a bacterial expression system. The recombinant protein contains six histidine residues at its amino terminus (His-tag). The molecular weight of 6xHis tagged R-Ras is approximately 25 kDa and it is supplied as a white lyophilized powder.

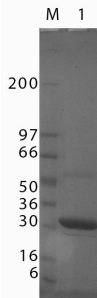
Spaargaren M, Martin GA, McCormick F, Fernandez-Sarabia MJ, Bischoff JR. The Ras-related protein R-ras interacts directly with Raf-1 in a GTP-dependent manner. *Biochem J.* 1994;300 ( Pt 2) (Pt 2):303-307. doi:10.1042/bj3000303

### 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 60 µl of Milli-Q water (300 µg size). When reconstituted, the protein will be in the following buffer: 50 mM Tris pH 7.5, 50 mM NaCl, 0.5 mM MgCl<sub>2</sub>, 5% (w/v) sucrose, and 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 should not be exposed to repeated freeze-thaw cycles.** The lyophilized protein is stable at 4°C desiccated (<10% humidity) for one year.

### Purity

Protein purity is determined by scanning densitometry of Coomassie Blue-stained protein on a 4-20% polyacrylamide gradient gel. His tagged R-Ras protein was determined to be >90% pure. (see Figure 1).



**Figure 1. R-Ras Protein Purity Determination.** A 10 µg sample of recombinant R-Ras protein (molecular weight approx. 25 kDa) was separated by electrophoresis in a 4-20% SDS-PAGE system and stained with Coomassie Blue. Protein quantitation was deter-

mined using the Precision Red Protein Assay Reagent (Cat. # ADV02). Mark12 molecular weight markers are from Life Technologies Inc.

### Biological Activity Assay

The biological activity of R-Ras can be determined from the ability of the RasGRF1 exchange domain (Ras-GRF1-ExD) to catalyze the exchange of GDP for GTP on R-Ras. A standard biological assay for monitoring the biological activity of R-Ras is an exchange assay utilizing the 2X Exchange Buffer from the RhoGEF exchange assay kit (Cat.# BK100) and the human RasGRF GEF domain (Cat.# CS-GE03).

### Reagents

1. Recombinant R-Ras protein (Cat.# CS-RS05)
2. Recombinant RAS-GRF GEF protein (Cat.# CS-GE03)
3. 2X Exchange Buffer (40 mM Tris pH 7.5, 100 mM NaCl, 20 mM MgCl<sub>2</sub>, 0.1 mg/ml BSA, 1.5 µM N-MAR-GTP)
4. Dilution Buffer (20 mM Tris pH 7.5, 50 mM NaCl, 10 mM MgCl<sub>2</sub>, 0.1 mg/ml BSA)

### Equipment

1. Fluorescence spectrophotometer (λ<sub>exc</sub>=485 nm, λ<sub>em</sub>=515 nm)
2. Corning 96-well half area plates (Cat. # 3686) or other plate with low protein binding surface.

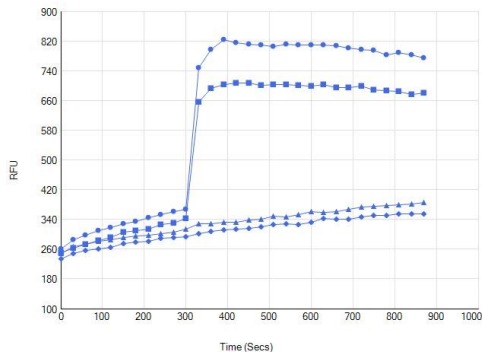
### Method

1. Dilute Ras-GRF GEF protein (Cat.# CS-GE03) to 1 µM (0.06 mg/ml) with Dilution Buffer.
2. Dilute R-Ras to 50 µM (1.1 mg/ml) with Dilution Buffer.
3. Dissolve lyophilized 2X Exchange Buffer in 5 ml Milli-Q water and keep at room temperature.
5. Set up the plate reader for kinetic fluorescence measurements (Excitation wavelength at 485 nm and emission wavelength at 515nm) with readings every 30 seconds for 30 minutes.
6. Add the following components together and mix well by gentle pipetting:

Exchange reaction mix	96 well black plate
2X Exchange Buffer	50 µl
dH <sub>2</sub> O	26 µl
50 µM R-Ras	4 µl

5. Pipette 20 µl of 1 µM Ras-GRF GEF protein or Dilution Buffer into their respective wells and immediately pipette up and down twice and begin reading the fluorescence.
6. Once the readings are complete and the plate reader file has been saved, the exchange rate can be calculated by reducing the data to Vmax with the software that accompanies the plate reader.

**Figure 2. EDTA mediated N-MAR-GTP exchange in to R-Ras.**



Legend: R-Ras protein was prepared in Exchange buffer to 100 µg/ml and 100 µl was pipetted in to duplicate wells of a 96-well half-area plate. To initiate the N-MAR-GTP exchange reaction, 10 mM EDTA was added to two wells at T = 300 s. Data was obtained using a Molecular Device Spectrofluor M2. See increase in fluorescence at T = 300s which relates to exchange of GTP on R-Ras for N-MAR-GTP reporter.

#### Product Uses

- Study of R-Ras exchange activity with different GEFs
- Identification of R-Ras exchange factors (GEFs)
- Positive control for GEF studies
- Biochemical characterization of R-Ras protein interactions
- Western blot standard

#### Product Citations/Related Products

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