

## RHEB Protein

Full-length recombinant protein expressed in Sf9 cells

**Catalog # R42-30H**

Lot # 1098-3

### Product Description

Recombinant full-length human RHEB was expressed by baculovirus in Sf9 cells using an N-terminal His tag. The gene accession number is [NM\\_005614](#).

### Gene Aliases

RHEB2

### Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 2mM DTT, 25% glycerol.

### Storage and Stability

Store product at  $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

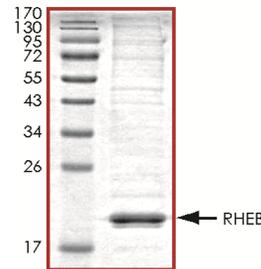
### Scientific Background

RHEB is a member of the small GTPase superfamily and encodes a lipid-anchored, cell membrane protein with five repeats of the RAS-related GTP-binding region. RHEB is vital in the regulation of growth and cell cycle progression. RHEB has GTPase activity and shuttles between a GDP-bound form and a GTP-bound form. Drosophila RHEB is a direct target of Tsc2 GTPase-activating protein (GAP) activity both in vivo and in vitro. Point mutations in the Drosophila RHEB gene inhibit cell growth while overexpression promotes cell growth. RHEB functions downstream of Tsc1 and Tsc2 in the mTOR signaling pathway.

### References

1. Stocker, H. et al: Rheb is an essential regulator of S6K in controlling cell growth in Drosophila. *Nature Cell Biol.* 5: 559-565, 2003.
2. Zhang, Y. et al: Rheb is a direct target of the tuberous sclerosis tumour suppressor proteins. *Nature Cell Biol.* 5: 578-581, 2003.

### Purity



The purity was determined to be **>90%** by densitometry. Approx. MW **21 kDa**.

## RHEB Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog Number	R42-30H
Specific Lot Number	1098-3
Purity	>90%
Concentration	0.2 µg/µl
Stability	1yr At $-70^{\circ}\text{C}$ from date of shipment
Storage & Shipping	Store product at $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**