

## Rab2<sup>His</sup>-GST

human, recombinant, *E. coli*

Cat. No.	Amount
PR-112	50 µg

For *in vitro* use only  
Quality guaranteed for 12 months  
Store at -80°C

### Avoid freeze / thaw cycles

### Form

Liquid. Supplied as a 1 mg/ml solution in 25 mM HEPES, pH 7.2, 40 mM NaCl, 5 mM DTT, 2 mM MgCl<sub>2</sub> and 10 mM GDP.

### Activity

GTPγS binding: >750 mmol/mol.

### Molecular weight

53,5 kDa.

### Purity

≥90% by SDS-PAGE.

### Description

Rab proteins are small-molecular-weight guanosine triphosphatases (GTPases) that control vesicular traffic in eukaryotic cells. The small GTPase Rab2 is a resident of pre-Golgi intermediates and is required for protein transport from the endoplasmic reticulum to the Golgi complex.

The GST-Tag facilitates the protein's application in typical GST pull-down assays.

### Selected References:

Tisdale E.J. and Balch W.E. (1996) Rab2 is essential for the maturation of pre-Golgi intermediates. *J. Biol. Chem.* **271**:29372.

### Associated products available from Jena Bioscience

**Non-hydrolyzable GTP-analogs** (such as GTPγS, GppCp, GppNHp, NPE-caged-GTP...)

**Fluorescent GTPs** (such as MANT, ANT, TNP, or GTP labeled with a dye of your choice)

**Recombinant G-protein signaling reagents** (including GAPs, GEFs, heterotrimeric G-proteins, and many more...)

For further information please visit [www.jenabioscience.com](http://www.jenabioscience.com).