

## Rab1B<sup>GST-His</sup>

rat, recombinant, *E. coli*

Cat. No.	Amount
PR-114	50 µg

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

### Avoid freeze / thaw cycles

### Form

Liquid. Supplied in 50 mM Tris, pH 8.0, 200 mM NaCl, 2 mM MgCl<sub>2</sub>, 10 µM GDP and 2 mM β-mercaptoethanol.

### Activity

100 pmol of protein can bind > 80 pmol of GDP.

### Purity

≥90% by SDS-PAGE.

### Description

Rab1B is a small GTPase that belongs to the Ras superfamily. Rab proteins play an important role in various aspects of membrane traffic, including cargo selection, vesicle budding, vesicle motility, tethering, docking, and fusion. For example, Rab1A (cat.# PR-113), Rab1B, and Rab2 (cat.# PR-111, PR-112) are located in the ER to Golgi intermediate compartment and cis Golgi cisternae and are required for ER to Golgi and intra Golgi transport. The GST-Tag facilitates the protein's application in typical GST pull-down assays.

### Selected References:

- Stenmark H. and Olkkonen V.M. (2001) The Rab GTPase family. *Genome Biol.* **2**:30071.  
Plutner *et al.* (1991) Rab1b regulates vesicular transport between the endoplasmic reticulum and successive Golgi compartments. *J. Cell Biol.* **115**:31.  
Jin *et al.* (1996) Rab1a and multiple other Rab proteins are associated with the transcytotic pathway in rat liver. *J. Biol. Chem.* **271**:30105.

### 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).