

Rab4B^{GST-His}

human, recombinant, *E. coli*

Cat. No.	Amount
PR-183	

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

Avoid freeze / thaw cycles

Activity

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

Form

Liquid. Supplied in 50 mM Tris, pH 8.0, 100 mM NaCl, 10 mM MgCl₂, 10 μM GDP and 2 mM mercaptoethanol.

Purity

≥ 90% by SDS-PAGE.

Description

Rab4B 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.

Rab4 is associated with early endosomes and regulates recycling vesicle formation. Together with Rab5 it acts to control influx and efflux out of early endosomes.

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.
- Mohrmann et al. (2002) Rab4 function in membrane recycling from early endosomes depends on a membrane to cytoplasm cycle. *J. Biol. Chem.* **277**:32029.
- Somsel Rodman J. and Wandinger-Ness A. (2000) Rab GTPases coordinate endocytosis. *J. Cell Sci.* **113**:183.
- Chamberlain et al. (2004) The p85 alpha Subunit of Phosphatidylinositol 3'-Kinase Binds to and Stimulates the GTPase Activity of Rab Proteins. *J. Biol. Chem.* **279**:48607.