

Rab4A^{GST-His}

human, recombinant, *E. coli*

Cat. No.	Amount
PR-116	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, 10 mM MgCl₂ and 2 mM β-mercaptoethanol.

Activity

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

Purity

>90% by SDS-PAGE.

Description

Rab4A 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. The monomeric GTPase Rab4 is associated with early endosomes and regulates recycling vesicle formation. Together with Rab5 it act 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:

- Chamberlain et al. (2008) Disrupted RabGAP Function of the p85 Subunit of Phosphatidylinositol 3-Kinase Results in Cell Transformation. *J. Biol. Chem.* **283**:15861-15868
- 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 p85α Subunit of Phosphatidylinositol 3'-Kinase Binds to and Stimulates the GTPase Activity of Rab Proteins. *J. Biol. Chem.* **279**:48607.

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.