

**Rab4B**

Ras-associated, small GTP-binding protein  
human, recombinant, *E. coli*

Cat. No.	Amount
PR-183	50 µg

**For general laboratory use.**

**Shipping:** shipped on dry ice

**Storage Conditions:** store at -80 °C

**Additional Storage Conditions:** avoid freeze/thaw cycles

**Shelf Life:** 12 months

**Accession number:** NM\_004578

**Purity:** > 90 % (SDS-PAGE)

**Form:** liquid (Supplied in 50 mM Tris-HCl pH 8.0, 100 mM NaCl, 10 mM MgCl<sub>2</sub>, 10 µM GDP and 2 mM mercaptoethanol)

**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.

**Activity:**

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

**Selected References:**

Stenmark *et al.* (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 *et al.* (2000) Rab GTPases coordinate endocytosis. *J. Cell Sci.* **113**:183.

Chamberlain *et al.* (2004) The p85alpha Subunit of Phosphatidylinositol 3-kinase Binds to and Stimulates the GTPase Activity of Rab Proteins. *J. Biol. Chem.* **279** (47):48607.