

**Rab19**

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

Cat. No.	Amount
PR-189	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: NP_035356

Purity: > 90 % (SDS-PAGE)

Form: liquid (Supplied in 50 mM Tris-HCl pH 8.0, 100 mM NaCl, 10 mM MgCl₂, 10 µM GDP and 2 mM mercaptoethanol)

Description:

Rab19 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. Rab 19 is expressed at high levels in intestine, lung and spleen, and at a lower level in kidney.

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.