

**Rab24^{GST-His}**

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

Cat. No.	Amount
PR-120	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

Molecular Weight: 52 kDa

Accession number: NM_009000

Purity: > 90 % (SDS-PAGE)

Form: liquid (Supplied in 25 mM HEPES pH 7.2, 40 mM NaCl, 2 mM MgCl₂, 10 µM GDP and 3 mM DTE)

Description:

N-terminal tagged Rab24 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. Rab24 possesses several unusual characteristics that distinguish it from other Rab proteins. Rab24 exists predominantly in the GTP state when expressed in cultured cells. The low GTPase activity is related to the presence of serine instead of glutamine at the position cognate to Ras Gln-61. Rab24 was found in the endoplasmic reticulum/cis-Golgi region and on late endosomal structures. The localization of Rab24 may indicate its involvement in autophagy-related processes. The GST-Tag facilitates the protein's application in typical GST pull-down assays.

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.

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

Erdman *et al.* (2000) Rab24 Is an Atypical Member of the Rab GTPase Family. Deficient GTPase activity, GDP dissociation inhibitor interaction, and prenylation of Rab24 expressed in cultured cells. *J. Biol. Chem.* **275**:3848.

Olkkonen *et al.* (1993) Molecular cloning and subcellular localization of three GTP-binding proteins of the rab subfamily. *J. Cell Sci.* **106**:1249.