

**Rab23**

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

Cat. No.	Amount
PR-191	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_033025

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:

Rab23 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 23 is expressed at high levels in brain and localized to plasma membrane and endosomes. It regulates ciliary and sonic hedgehog signalling pathways and is involved in embryogenesis.

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.

Yamaguchi *et al.* (2002) A GDP/GTP exchange protein for the Rab3 small G protein family up-regulates a postdocking step of synaptic exocytosis in central synapses. *Proc. Natl. Acad. Sci. USA.* 99:14536.