

**Rab22A**

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

Cat. No.	Amount
PR-190	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\_065724

**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:**

Rab22A 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. Rab22A is found in early endosomes and trans Golgi. It interacts with EEA 1 and controls endosomal membrane trafficking.

**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.