

Rab3C^{GST-His}

bovine, recombinant, *E. coli*

Cat. No.	Amount
PR-182	50 µg

For *in vitro* use only
Quality guaranteed for 12 months
Store at -80°C

Avoid freeze / thaw cycles

Activity

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

Form

Liquid. Supplied in 50 mM Tris, pH 8.0, 100 mM NaCl, 10 mM MgCl₂, 10 µM GDP and 2 mM mercaptoethanol.

Purity

≥ 90% by SDS-PAGE.

Description

Rab3C 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.

The Rab3C recycling is coupled with synaptic and secretory vesicle trafficking. It facilitates translocation and docking of the synaptic vesicles to the active zone.

The GST-Tag facilitates the protein's application in typical GST pull-down assays.

Selected References:

- Stenmark H. and Olkkonen V.M. (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.