

**Rab27A^{GST-His}**

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

Cat. No.	Amount
PR-122	50 µg

For in vitro use only!

Shipping: shipped on dry ice

Storage Conditions: store at -80 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Molecular Weight: 54 kDa

Accession number: P23640

Purity: > 90 % (SDS-PAGE)

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

Description:

N-terminal tagged Rab27 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 monomeric GTPase Rab27 subfamily regulates the exocytosis of these cell-specific store organelles. Rab27 acts through organelle-specific effector proteins, such as granuphilin in pancreatic beta cells and melanophilin in melanocytes. The Rab27 and effector complex then interacts with proteins that are essential for membrane transport and fusion. 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.

Izumi *et al.* (2003) The roles of Rab27 and its effectors in the regulated secretory pathways. *Cell Struct. Funct.* **28**:465.