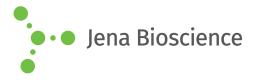
# **DATA SHEET**





## Rab31

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

Cat. No.	Amount
PR-193	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: 26 kDa

Accession number: NP\_006859

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:**

Rab31 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 small Rab31 GTPase is associated with early endosomes and trans Golgi. It regulates biosynthetic pathways from endosomes to Golgi. It plays an important role in phagosome maturation.

### 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.

