

**Rab2<sup>GST-His</sup>**

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

Cat. No.	Amount
PR-112	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:** 53.5 kDa

**Accession number:** NM\_002865

**Purity:** > 90 % (SDS-PAGE)

**Form:** liquid (Supplied in 25 mM HEPES pH 7.2, 40 mM NaCl, 5 mM DTT, 2 mM MgCl<sub>2</sub> and 10 mM GDP)

**Description:**

Rab proteins are small-molecular-weight guanosine triphosphatases (GTPases) that control vesicular traffic in eukaryotic cells. The small GTPase Rab2 is a resident of pre-Golgi intermediates and is required for protein transport from the endoplasmic reticulum to the Golgi complex. The GST-Tag facilitates the protein's application in typical GST pull-down assays.

**Activity:**

GTPyS binding: >750 mmol/mol.

**Selected References:**

Tisdale *et al.* (1996) Rab2 is essential for the maturation of pre-Golgi intermediates. *J. Biol. Chem.* **271**:29372.