

Rab2

human, recombinant, *E. coli*

Cat. No.	Amount
PR-111	50 μ g

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

Avoid freeze / thaw cycles

Form

Liquid. Supplied as a 1 mg/ml solution in 25 mM HEPES, pH 7.2, 40 mM NaCl, 5 mM DTT, 2 mM MgCl₂ and 10 mM GDP.

Activity

GTP γ S binding: >750 mmol/mol.

Molecular weight

23.5 kDa.

Purity

\geq 90% by SDS-PAGE.

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.

Selected References:

Tisdale E.J. and Balch W.E. (1996) Rab2 is essential for the maturation of pre-Golgi intermediates. *J. Biol. Chem.* **271**:29372.

Associated products available from Jena Bioscience

Non-hydrolyzable GTP-analogs (such as GTP γ S, GppCp, GppNHp, NPE-caged-GTP...)

Fluorescent GTPs (such as MANT, ANT, TNP, or GTP labeled with a dye of your choice)

Recombinant G-protein signaling reagents (including GAPs, GEFs, heterotrimeric G-proteins, and many more...)

For further information please visit www.jenabioscience.com.