

REP-1^{His}

Rab escort Protein

rat, recombinant, *Spodoptera frugiperda*

Cat. No.	Amount
PR-105	50 µg

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

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 25 mM HEPES pH 7.2, 40 mM NaCl and 5 mM DTT.

Activity

1 pmol of REP-1 will mediate transfer of 2 pmol of geranylgeraniol by RabGGTase onto 1 pmol of Rab7 in 3 minutes.

Purity

> 90% by SDS-PAGE

Description

C-terminal His-tagged REP-1 (Rab escort protein-1) binds to a wide range of RabGGTase with nanomolar affinities and supports the prenylation of RabGTPase mediated by RabGGTase.

Selected References:

- Seabra *et al.* (1995) Deficient geranylgeranylation of Ram/Rab27 in choroideremia. *J. Biol. Chem.* **270**:24420.
Alexandrov *et al.* (1994) Rab escort protein-1 is a multifunctional protein that accompanies newly prenylated rab proteins to their target membranes. *EMBO J.* **13**:5262.
Watanabe *et al.* (2008) Inhibitors of Protein Geranylgeranyltransferase I and Rab Geranylgeranyltransferase Identified from a Library of Allenoate-derived Compounds. *J. Biol. Chem.* **283**(15):9571-9579.

Associated products available from Jena Bioscience

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

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

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

For further information please visit www.jenabioscience.com.