

REP-1^{His}

(Rab escort protein)

Rat, Recombinant, *Spodoptera frugiperda*

Cat. No.	Amount
PR-105	50 µg

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

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.

AVOID FREEZE/THAW CYCLES.

For in vitro use only!

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

Store: -80 °C

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.

Associated products available from Jena Bioscience

Non-hydrolyzable GTP-analogs (such as GTPγS, Gppppp, GppNHp, NPE-caged-GTP...)
Activity:

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

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

For detailed information please view the sections on

<http://www.jenabioscience.com>