

REP-2^{His}

Rab escort Protein

human, recombinant, *Saccharomyces cerevisiae*

Cat. No.	Amount
PR-104	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-2 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-2 (Rab escort protein-2) binds to a wide range of RabGGTase with nanomolar affinities and supports the prenylation of RabGTPase mediated by RabGGTase, with an exception of Rab27.

Selected References:

Cremers *et al.* (1994) REP-2, a Rab escort protein encoded by the choroideremia-like gene. *J. Biol. Chem.* **269**:2111.

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.