

REP-2^{His}

(Rab Escort Protein)

human, Recombinant, *Saccharomyces cerevisiae*

Cat. No.	Amount
PR-104	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-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.

AVOID FREEZE/THAW CYCLES.

For in vitro use only!

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

Store: -80 °C

Selected References:

Creemers *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

Non-hydrolyzable GTP-analogs (such as GTPγS, 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, GTPases, and many more...

For detailed information please view the sections on

Activity:

<http://www.jenabioscience.com>