

**REP-1^{His}**

Rab Escort Protein
fish, recombinant, *E. coli*

Cat. No.	Amount
PR-105	50 µg

For general laboratory use.

Shipping: shipped on dry ice

Storage Conditions: store at -80 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Purity: > 90 % (SDS-PAGE)

Form: liquid (Supplied in 25 mM HEPES pH 7.2, 40 mM NaCl and 5 mM DTT)

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.

Activity:

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

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 Alkenoate-derived Compounds. *J. Biol. Chem.* **283** (15):9571-9579.