

GGTase-I^{GST}

Protein geranylgeranyltransferase type I, α - and β -subunit

rat, recombinant, *E. coli*

Cat. No.	Amount
PR-360	50 μ g

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, 25 mM NaCl and 5 mM DTT.

Activity

1 pmol of GGTase-I will transfer 5 pmol of Farnesyl to RhoA in 15 min at 37°C.

Molecular Weight

α 44 kDa and β 38 kDa

Purity

> 90% by SDS-PAGE

Description

GGTase-I (Geranylgeranyltransferase-I) catalyzes the transfer of the farnesyl and geranylgeranyl groups from farnesyl and geranylgeranyl-diphosphate to proteins containing a C-terminal CaaX motif, where 'C' is a conserved cysteine that is the site of farnesyl modification, 'a' is usually an aliphatic amino acid, and 'X' is leucine or phenylalanine. Farnesyltransferase (FT) and GGTase-I are closely related, sharing a common α subunit and 30% identity in their β subunits.

The GST-Tag facilitates the protein's application in typical GST pull-down assays.

Selected References:

Lackner *et al.* (2005) Chemical genetics identifies Rab geranylgeranyl transferase as an apoptotic target of farnesyl transferase inhibitors. *Cancer Cell*. **7**:325.

Yokoyama *et al.* (1993) Purification of a mammalian protein geranylgeranyltransferase. Formation and catalytic properties of an enzyme-geranylgeranyl pyrophosphate complex. *J. Biol. Chem.* **268**:4055.

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.