

GGTase-I^{GST} (Geranyl-Geranyl-Transferase) Rat, Recombinant, *E. coli*

Cat. No.	Amount
PR-360	50 µg

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

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.

AVOID FREEZE/THAW CYCLES!

For in vitro use only!

Purity: > 90% by SDS-PAGE

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

Store: -80 °C

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

Non-hydrolyzable GTP-analogs (such as GTP γ S, 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

<http://www.jenabioscience.com>