

## FTase<sup>GST</sup>

### Protein farnesyltransferase, $\alpha$ - and $\beta$ -subunit rat, recombinant, *E. coli*

Cat. No.	Amount
PR-361	50 $\mu$ g

For *in vitro* use only  
Quality guaranteed for 12 months  
Store at -80°C

#### Avoid freeze / thaw cycles

#### Form

Liquid. Supplied in 100 mM HEPES pH 7.2, 100 mM NaCl and 5 mM DTT.

#### Activity

1 pmol of FTase will transfer 1 pmol of farnesyl to H-Ras in 15 min at 37°C.

#### Molecular Weight

$\alpha$ -subunit: 44 kDa and  $\beta$ -subunit: 35 kDa

#### Purity

> 90% by SDS-PAGE

#### Description

FTase catalyzes the transfer of the farnesyl group from farnesyl 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 methionine, serine, glutamine, or alanine. 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.  
Zimmerman *et al.* (1998) High-level expression of rat farnesyl:proteintransferase in *Escherichia coli* as a translationally coupled heterodimer. *Protein Express. Purif.* **14**:395.

#### 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 $\gamma$ 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](http://www.jenabioscience.com).