

FTase

(Protein Farnesyltransferase, α - and β -subunit)

Rat, Recombinant, *E. coli*

Cat. No.	Amount
PR-102	50 μ g

Liquid. Supplied as a 1 mg/ml solution in 100 mM HEPES, pH 7.2, 100 mM NaCl, and 5 mM DTT.

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.

AVOID FREEZE/THAW CYCLES.

For in vitro use only!

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

Purity: \geq 90% by SDS-PAGE.

Molecular weight:

α -subunit: 44 kDa and β -subunit: 35 kDa

Store: -80 °

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

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>