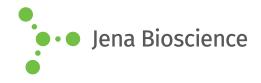
# **DATA SHEET**





## **■ FTase**<sub>W102T\_Y154T</sub>GST

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

Cat. No.	Amount
PR-952	20 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

Molecular Weight: GST-α: 73 kDa, β: 48 kDa

Purity: > 90 % (SDS-PAGE)

Form: liquid (Supplied in 25 mM HEPES pH 7.2, 40 mM NaCl and 2 mM

DTT)

#### **Description:**

FTase is one of the three mammalian protein prenyltransferases. The double mutant of FTase was engineered to efficiently transfer biotin-geranylpyrophosphate (BGPP, Cat.# LI-015) onto its native protein substrates. This approach allows the detection of femtomolar amounts of prenylatable proteins in eukaryotic cells and tissues as well as effective inhibitor screening.

#### Activity:

1 pmol of FTase double mutant will transfer 0.4 pmol of BGPP to H-Ras in 15 min at 37  $^{\circ}\text{C}.$ 

#### Selected References:

Hougland et al. (2009) Getting a handle on protein prenylation. Nat. Chem. Biol. 5 (4):197.

Nguyen et al. (2009) Analysis of the eukaryotic prenylome by isoprenoid affinity tagging. Nat. Chem. Biol. 5 (4):227.