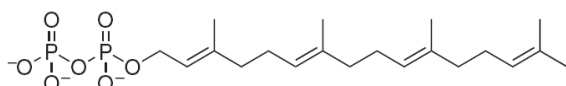


## GGPP

### Geranylgeranyl pyrophosphate, ammonium salt

### Natural lipid donor for protein geranylgeranyltransferases (RabGGTase, GGTase-I)

Cat. No.	Amount
LI-017	200 µg



For *in vitro* use only  
Quality guaranteed for 12 months  
Store at -20°C

#### Avoid freeze-thaw cycles

#### Form

Liquid, 1 mg/ml supplied as a solution in methanol:  
NH<sub>4</sub>OH (7:3)

#### Molecular weight

450.44 g/mol (free acid)

#### Molecular Formula

C<sub>20</sub>H<sub>36</sub>O<sub>7</sub>P<sub>2</sub> (free acid)

#### Purity

>95% by TLC

#### CAS Number

6699-20-3 (free acid)

#### Description

GGPP is the natural prenyl donor for protein geranylgeranyltransferases (RabGGTase, GGTase-I), therefore it's suitable as control in prenylation assays.

#### Selected References:

Lane *et al.* (2006) Structural biology of protein farnesyltransferase and geranylgeranyltransferase type I. *Journal of Lipid Research* **47**:681