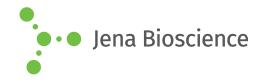
DATA SHEET





B-GPP

Biotin-labeled geranyl pyrophosphate, triethylammonium salt

| Cat. No. | Amount |
|----------|--------|
| LI-015 | 20 μl |

Structural formula of B-GPP

For general laboratory use.

Shipping: shipped on gel packs **Storage Conditions:** store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Molecular Formula: C₂₀H₃₅N₃O₉P₂S **Molecular Weight:** 555.52 g/mol

Purity: mass identification conforms (ESI-MS)

Form: liquid (Supplied as 1 mM solution in 25 mM (NH₄)₂CO₃)

Concentration: 1 mM

Description:

B-GPP is a biotin-labeled analog of geranyl pyrophosphate (GPP). It serves as lipid donor for the native RabGGTase and mutants of GGTase-I and FTase. B-GPP allows for straightforward mammalian prenylome labeling and thus, not only identifying protein prenyl-transferase substrates but efficiently screening for new inhibitors.

Selected References:

Kovtun *et al.* (2010) Towards the construction of expressed proteomes using a Leishmania tarentolae based cell-free expression system. *PLoS One.* **5** (12):e14388.

Nguyen et al. (2009) Analysis of the eukaryotic prenylome by isoprenoid affinity tagging. Nature Chemical Biology. **5 (4)**:227.