

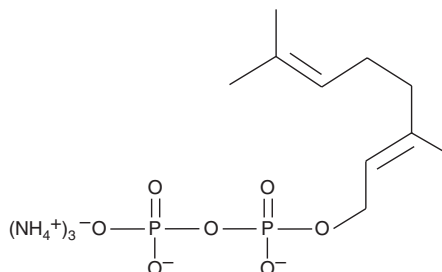
PRODUCT INFORMATION



Geranyl Pyrophosphate (triammonium salt)

Item No. 63320

CAS Registry No.: 116057-55-7
Formal Name: 3E,7-dimethyl-2,6-octadienyl-diphosphoric acid, triammonium salt
Synonyms: GDP, Geranyl Diphosphate, GPP
MF: C₁₀H₂₀O₇P₂ · 3NH₃
FW: 365.3
Purity: ≥90%
Supplied as: A solution in methanol
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Geranyl pyrophosphate (triammonium salt) is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol under a gentle stream of nitrogen and immediately add the solvent of choice. A stock solution may be made by dissolving the geranyl pyrophosphate (triammonium salt) in the solvent of choice. Geranyl pyrophosphate (triammonium salt) is slightly soluble in water.

Description

Geranyl pyrophosphate is an intermediate in the mevalonate pathway. It is formed from dimethylallyl pyrophosphate (DMAPP; Item No. 63180) and isopentenyl pyrophosphate by geranyl pyrophosphate synthase.¹ Geranyl pyrophosphate is used in the biosynthesis of farnesyl pyrophosphate (Item No. 63250), geranylgeranyl pyrophosphate (Item No. 63330), cholesterol, terpenes, and terpenoids.

Reference

1. Dorsey, J.K., Dorsey, J.A. and Porter, J.W. The purification and properties of pig liver geranyl pyrophosphate synthetase. *J. Biol. Chem.* **241(22)**, 5353-5360 (1966).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/08/2020

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM