

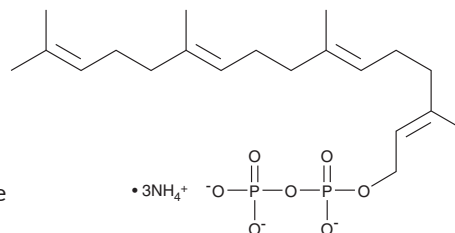
# PRODUCT INFORMATION



## Geranylgeranyl Pyrophosphate (ammonium salt)

Item No. 63330

**CAS Registry No.:** 313263-08-0  
**Formal Name:** P-[(2E,6E,10E)-3,7,11,15-tetramethyl-2,6,10,14-hexadecatetraen-1-yl] ester-diphosphoric acid, triammonium salt  
**Synonym:** GGPP  
**MF:** C<sub>20</sub>H<sub>33</sub>O<sub>7</sub>P<sub>2</sub> • 3NH<sub>4</sub>  
**FW:** 501.5  
**Purity:** ≥95%  
**Supplied as:** A 1 mg/ml solution in 10 mM ammonium hydroxide  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Description

GGPP is an intermediate in the HMG-CoA reductase pathway derived directly from farnesyl pyrophosphate and used in the biosynthesis of terpenes and terpenoids.<sup>1</sup> It also serves as a substrate in the prenylation of a variety of critical intracellular proteins including small GTPases.<sup>2</sup> This post-translational modification is necessary for correct localization of proteins to intracellular membranes for proper functionality and has become a focus of anticancer drug discovery.<sup>3,4</sup>

### References

1. Vance, D.E. Cholesterol and related derivatives. *Biochemistry*. Zubay, G., editor, 2<sup>nd</sup> edition., Macmillan Publishing Company (1988).
2. McTaggart, S.J. Isoprenylated proteins. *Cell Mol. Life Sci.* **63(3)**, 255-267 (2006).
3. Baron, R.A. and Seabra, M.C. Rab geranylgeranylation occurs preferentially via the pre-formed REP-RGGT complex and is regulated by geranylgeranyl pyrophosphate. *Biochem. J.* **415(1)**, 67-75 (2008).
4. Maynor, M., Scott, S.A., Rickert, E.L., et al. Synthesis and evaluation of 3- and 7-substituted geranylgeranyl pyrophosphate analogs. *Bioorg. Med. Chem. Lett.* **18(6)**, 1889-1892 (2008).

#### 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, 07/29/2025

#### 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