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:	$C_{20}H_{33}O_7P_2 \bullet 3NH_4$	
FW:	501.5	0 0
Purity:	≥95%	• 3NH4+ -O_P_O_P_O
Supplied as:	A solution in 10 mM ammonium hydroxide	
Storage:	-20°C	0 0
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.¹ It is also serves as a substrate in the prenylation of a variety of critical intracellular proteins including small GTPases.² 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.^{3,4}

References

- 1. Vance, D.E. Cholesterol and related derivatives. Biochemistry. Zubay, G., editor, 2nd 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.

SAFFTY 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.

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, 04/14/2023

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