

PRODUCT INFORMATION

Isopentenyl Pyrophosphate (ammonium salt)

Item No. 63480

Formal Name: 3-methylbut-3-en-1-yl
diphosphate, triammonium salt

Synonym: IPP

MF: $C_5H_9O_7P_2 \cdot 3NH_4$

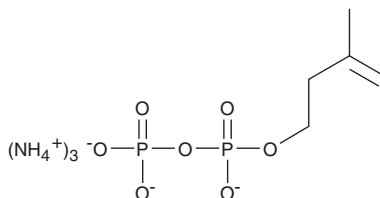
FW: 297.2

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^\circ C$

Stability: ≥ 4 years



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

Laboratory Procedures

Isopentenyl pyrophosphate (IPP) (ammonium salt) is supplied as a solid. A stock solution may be made by dissolving the IPP (ammonium salt) in the solvent of choice, which should be purged with an inert gas. IPP (ammonium salt) is soluble in methanol.

Description

IPP is a precursor in the biosynthesis of isoprenoids.¹ It is formed from acetyl-CoA (Item No. 16160) via the mevalonate pathway, and it can be isomerized to dimethylallyl diphosphate (DMAPP; Item No. 63180) by IPP isomerase. IPP can also be formed from pyruvate and glyceraldehyde-3-phosphate via the nonmevalonate pathway.

Reference

1. Chatzivasileiou, A.O., Ward, V., Edgar, S.M., et al. Two-step pathway for isoprenoid synthesis. *Proc. Natl. Acad. Sci. USA* **116**(2), 506-511 (2019).

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, 11/23/2022

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