

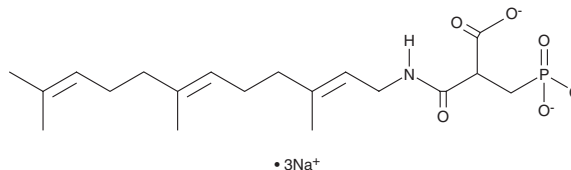
# PRODUCT INFORMATION



## FPT Inhibitor I

Item No. 45256

**CAS Registry No.:** 162934-54-5  
**Formal Name:** (E,E)-3-oxo-2-(phosphonomethyl)-3-[(3,7,11-trimethyl-2,6,10-dodecatrienyl)amino]-propanoic acid, trisodium salt  
**MF:** C<sub>19</sub>H<sub>29</sub>NO<sub>6</sub>P • 3Na  
**FW:** 467.4  
**Purity:** ≥85%  
**Supplied as:** A solid  
**Storage:** -80°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

FPT inhibitor I is supplied as a solid. A stock solution may be made by dissolving the FPT inhibitor I in water. FPT inhibitor I is slightly soluble (0.1-1 mg/ml) in water. We do not recommend storing the aqueous solution for more than one day.

### Description

FPT inhibitor I is a farnesyltransferase (FTase) inhibitor ( $IC_{50} = 0.083 \mu M$ ).<sup>1</sup> It is selective for FTase over geranylgeranyltransferase I (GGTase I), geranylgeranyltransferase II (GGTase II), squalene synthetase (SQS), and mevalonate kinase (MK;  $IC_{50}$ s = 26, 47, 516, and >200  $\mu M$ , respectively). FPT inhibitor I decreases FTase-mediated farnesylation of H-Ras in a cell-free assay.

### Reference

1. Manne, V., Ricca, C.S., Brown, J.G., *et al.* Ras farnesylation as a target for novel antitumor agents: Potent and selective farnesyl diphosphate analog inhibitors of farnesyltransferase. *Drug Develop. Res.* **34(2)**, 121-137 (1995).

#### 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, 02/26/2026

#### 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