

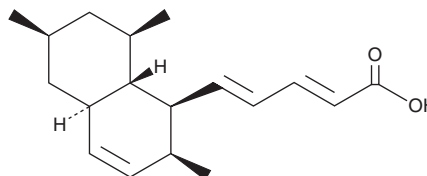
PRODUCT INFORMATION



Tanzawaic Acid B

Item No. 40909

CAS Registry No.: 169181-33-3
Formal Name: (2E,4E)-5-[(1S,2S,4aR,6S,8R,8aS)-1,2,4a,5,6,7,8,8a-octahydro-2,6,8-trimethyl-1-naphthalenyl]-2,4-pentadienoic acid
Synonym: (+)-Tanzawaic Acid B
MF: C₁₈H₂₆O₂
FW: 274.4
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4years
Item Origin: Bacterium/*Penicillium* sp.



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

Laboratory Procedures

Tanzawaic acid B is supplied as a solid. A stock solution may be made by dissolving the tanzawaic acid B in the solvent of choice, which should be purged with an inert gas. Tanzawaic acid B is soluble in ethanol, methanol, dichloromethane, and DMSO.

Description

Tanzawaic acid B is a fungal metabolite that has been found in *Penicillium* and has diverse biological activities.^{1,2} It inhibits protein tyrosine phosphatase 1B (PTP1B) in a cell-free assay (IC₅₀ = 8.2 μM) and LPS-induced production of nitric oxide (NO) in BV-2 microglia (IC₅₀ = 42.5 μM).¹ Tanzawaic acid B is active against chloroquine-resistant and chloroquine-sensitive *P. falciparum* strains (IC₅₀s = 78.5 and 79.2 μM, respectively).²

References

1. Quang, T.H., N.T.T., N., Ko, W., et al. Tanzawaic acid derivatives from a marine isolate of *Penicillium* sp. (SF-6013) with anti-inflammatory and PTP1B inhibitory activities. *Bioorg. Med. Chem. Lett.* **24**(24), 5787-5791 (2014).
2. Matsuo, H., Hokari, R., Ishiyama, A., et al. Hatsusamides A and B: Two new metabolites produced by the deep-sea-derived fungal strain *Penicillium steckii* FKJ-0213. *Mar Drugs* **18**(10), 513 (2020).

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, 05/16/2024

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