

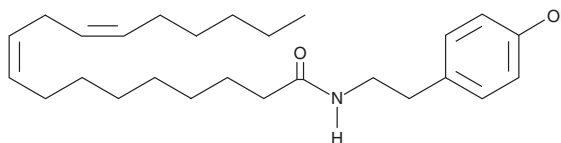
PRODUCT INFORMATION



Termitomycamide E

Item No. 11191

CAS Registry No.: 1177258-62-6
Formal Name: N-[2-(4-hydroxyphenyl)ethyl]-
9Z,12Z-octadecadienamide
Synonym: 2'-Deoxotermitomycamide A
MF: C₂₆H₄₁NO₂
FW: 399.6
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 225, 279, 349 nm



Laboratory Procedures

For long term storage, we suggest that termitomycamide E be stored as supplied at -20°C. It should be stable for at least two years.

Termitomycamide E is supplied as a crystalline solid. A stock solution may be made by dissolving the termitomycamide E in the solvent of choice. Termitomycamide E is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of termitomycamide E in these solvents is approximately 30 mg/ml.

Termitomycamide E is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, termitomycamide E should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Termitomycamide E has a solubility of approximately 0.05 mg/ml in a 1:20 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Endoplasmic reticulum (ER) stress, caused by accumulation of misfolded proteins and a disruption of calcium homeostasis, has been linked to several neuronal diseases including, Parkinson's, Alzheimer's, and prion diseases. A screen for protective activity against ER stress-dependent cell death identified termitomycamides, extracts from the fruiting bodies of *T. titanicus*, a notably large edible mushroom that is cultivated symbiotically in the nests of termites. Termitomycamide E is a fatty acid isolated from *T. titanicus* that at a dose of 0.1 µg/ml protects against ER stress-dependent cell death in Neuro2a cells induced by tunicamycin.¹

References

1. Choi, J.-H., Maeda, K., Nagai, K., *et al.* Termitomycamides A to E, fatty acid amides isolated from the mushroom *Termitomyces titanicus*, suppress endoplasmic reticulum stress. *Org. Lett.* **12(21)**, 5012-5015 (2010).

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, 08/19/2015

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