

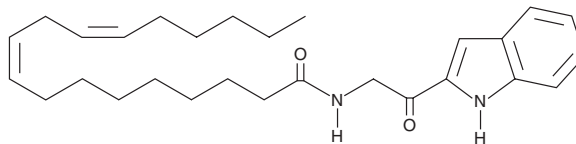
PRODUCT INFORMATION



Termitomycamide B

Item No. 11190

CAS Registry No.: 1254277-89-8
Formal Name: N-[2-(1H-indol-2-yl)-2-oxoethyl]-
9Z,12Z-octadecadienamide
MF: C₂₈H₄₀N₂O₂
FW: 436.6
Purity: ≥97%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 206, 240, 297 nm



Laboratory Procedures

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

Termitomycamide B is supplied as a crystalline solid. A stock solution may be made by dissolving the termitomycamide B in the solvent of choice. Termitomycamide B is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of termitomycamide B in these solvents is approximately 20 mg/ml.

Termitomycamide B is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. 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 B 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

Reference

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