

PRODUCT INFORMATION



Nanangenine G

Item No. 33522

Formal Name: 4,5-dihydroxy-3,4-bis(hydroxymethyl)-4a,8,8-trimethyl-1R,4R,4aR,5R,6,7,8,8aS-octahydronaphthalen-1-yl hexanoate

MF: C₂₁H₃₆O₆

FW: 384.5

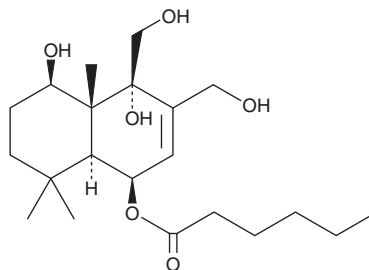
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Aspergillus nanangensis*



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

Laboratory Procedures

Nanangenine G is supplied as a solid. A stock solution may be made by dissolving the nanangenine G in the solvent of choice, which should be purged with an inert gas. Nanangenine G is soluble in methanol and DMSO.

Description

Nanangenine G is a drimane sesquiterpene that has been found in *Aspergillus*.¹ It is cytotoxic to NS-1 mouse myeloma cells (IC₅₀ = 47 µg/ml).

Reference

1. Lacey, H.J., Gilchrist, C.L.M., Crombie, A., et al. Nanangenines: Drimane sesquiterpenoids as the dominant metabolite cohort of a novel Australian fungus, *Aspergillus nanangensis*. *Beilstein J. Org. Chem.* **15**, 2631-2643 (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, 08/25/2023

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