

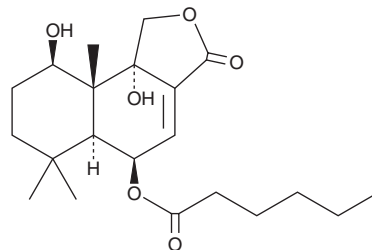
PRODUCT INFORMATION



Isonanangenine B

Item No. 33518

CAS Registry No.: 1623153-14-9
Formal Name: (5R,5aS,9R,9aR,9bS)-1,3,5,5a,6,7,8,9,9a,9b-decahydro-9,9b-dihydroxy-6,6,9a-trimethyl-3-oxonaphtho[1,2-c]furan-5-yl ester, hexanoic acid
Synonym: SF002-96-1
MF: C₂₁H₃₂O₆
FW: 380.5
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

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

Description

Isonanangenine B is a drimane sesquiterpene lactone that has been found in *Aspergillus*.^{1,2} It decreases survivin mRNA expression in a reporter assay when used at concentrations of 13.6, 18.42, and 26.31 μM. Isonanangenine B (18.42 μM) inhibits STAT3 and NF-κB binding to the survivin promoter in chromatin immunoprecipitation (ChIP) assays. It also induces apoptosis in COLO 320DM cells.

References

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).
2. Felix, S., Sandjo, L.P., Opatz, T., *et al.* SF002-96-1, a new drimane sesquiterpene lactone from an *Aspergillus* species, inhibits survivin expression. *Beilstein J. Org. Chem.* **9**, 2866-2876 (2013).

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