

PRODUCT INFORMATION



20-O-Acetylingenol-3-angelate

Item No. 39904

CAS Registry No.: 82425-35-2
Formal Name: (2Z)-2-methyl-2-butenic acid, (1aR,2S,5R,5aS,6S,8aS,9R,10aR)-4-[(acetyloxy)methyl]-1a,2,5,5a,6,9,10,10a-octahydro-5,5a-dihydroxy-1,1,7,9-tetramethyl-11-oxo-1H-2,8a-methanocyclopenta[a]cyclopropa[e]cyclodecen-6-yl ester

Synonyms: AAI, 3-O-Angeloyl-20-O-acetyl Ingenol, PEP008

MF: C₂₇H₃₆O₇

FW: 472.6

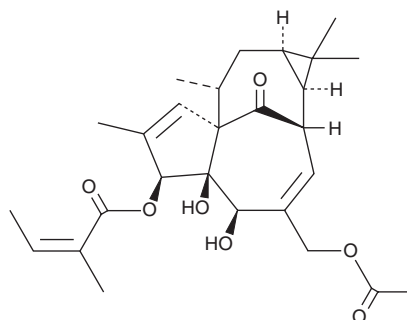
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/*Euphorbia pekinensis* Rupr.



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

Laboratory Procedures

20-O-Acetylingenol-3-angelate is supplied as a solid. A stock solution may be made by dissolving the 20-O-acetylingenol-3-angelate in the solvent of choice, which should be purged with an inert gas. 20-O-Acetylingenol-3-angelate is sparingly soluble (1-10 mg/ml) in chloroform and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

20-O-Acetylingenol-3-angelate is a diterpene ester that has been found in *E. peplus* and a derivative of ingenol-3-angelate (Item No. 16207) that has anticancer activity.^{1,2} It inhibits the growth of K562, HL-60, KT-1, MCF-7/adr, A549, and NIH3T3 cells when used at a concentration of 12.5 μM. 20-O-Acetylingenol-3-angelate (31.25-500 nM) induces cell cycle arrest at the G₂/M phase, apoptosis, and necrosis in K562 cells.

References

1. Mason, S.A., Cozzi, S.-J., Pierce, C.J., *et al.* The induction of senescence-like growth arrest by protein kinase C-activating diterpene esters in solid tumor cells. *Invest. New Drugs* **28**(5), 575-586 (2010).
2. Liu, M., Chen, F., Yu, R., *et al.* Synthesis and cytotoxicity against K562 cells of 3-O-angeloyl-20-O-acetyl ingenol, a derivative of ingenol mebutate. *Int. J. Mol. Sci.* **17**(8), 1348 (2016).

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