

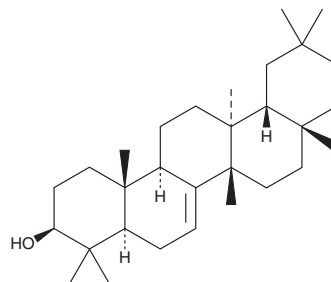
PRODUCT INFORMATION



Multiflorenol

Item No. 29709

CAS Registry No.: 2270-62-4
Formal Name: (14β)-13α-methyl-26-norolean-7-en-3β-ol
MF: C₃₀H₅₀O
FW: 426.7
Purity: ≥85%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Melicope octandra*



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

Laboratory Procedures

Multiflorenol is supplied as a solid. A stock solution may be made by dissolving the multiflorenol in the solvent of choice, which should be purged with an inert gas. Multiflorenol is soluble in organic solvents such as methanol.

Description

Multiflorenol is a triterpene that has been found in *T. kirilowii* seeds.¹ It inhibits *in vitro* activation of Epstein-Barr virus early antigen (EBV-EA) induced by the tumor promoter phorbol 12-myristate 13-acetate (TPA; Item No. 10008014) in a concentration-dependent manner.

Reference

1. Akihisa, T., Tokuda, H., Ichiishi, E., *et al.* Anti-tumor promoting effects of multiflorane-type triterpenoids and cytotoxic activity of karounidiol against human cancer cell lines. *Cancer Lett.* **173**(1), 9-14 (2001).

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, 01/26/2024

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