

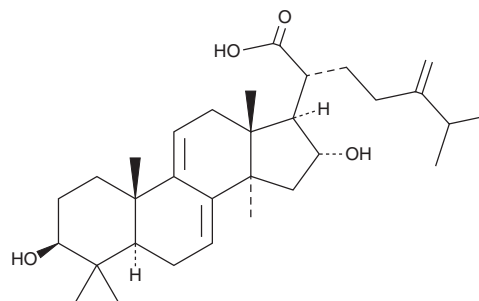
PRODUCT INFORMATION



Dehydrotumulosic Acid

Item No. 34938

CAS Registry No.: 6754-16-1
Formal Name: 3 β ,16 α -dihydroxy-24-methylene-lanosta-7,9(11)-dien-21-oic acid
MF: C₃₁H₄₈O₄
FW: 484.7
Purity: \geq 98%
UV/Vis.: λ_{max} : 244 nm
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Fungus/*Poria cocos*



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

Laboratory Procedures

Dehydrotumulosic acid is supplied as a solid. A stock solution may be made by dissolving the dehydrotumulosic acid in the solvent of choice, which should be purged with an inert gas. Dehydrotumulosic acid is soluble in ethanol.

Description

Dehydrotumulosic acid is a triterpene that has been found in *P. cocos* and has anticancer and anti-inflammatory activities.^{1,2} It enhances the cytotoxicity of vincristine (Item No. 11764) against multidrug-resistant KBV200 cells when used at concentrations of 12.5 or 25 $\mu\text{g}/\text{ml}$.¹ Dehydrotumulosic acid (50 mg/kg) reduces serotonin-induced hind paw edema, as well as ear edema induced by phorbol 12-myristate 13-acetate (TPA; Item No. 10008014), in mice.²

References

1. Shan, H., Qinglin, Z., Fengjun, X., *et al.* Reversal of multidrug resistance of KBV200 cells by triterpenoids isolated from *Poria cocos*. *Planta Med.* **78(5)**, 428-433 (2012).
2. Giner, E.M., Máñez, S., Recio, M.C., *et al.* *In vivo* studies on the anti-inflammatory activity of pachymic and dehydrotumulosic acids. *Planta Med.* **66(3)**, 221-227 (2000).

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