

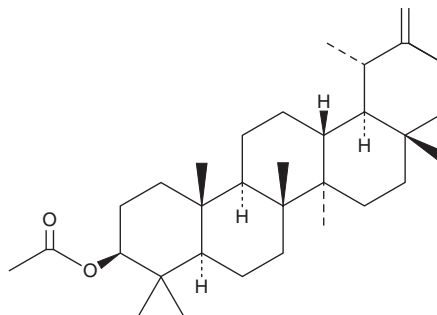
PRODUCT INFORMATION



Taraxasterol Acetate

Item No. 36131

CAS Registry No.: 6426-43-3
Formal Name: (3 β ,18 α ,19 α)-urs-20(30)-en-3-ol, 3-acetate
Synonyms: NSC 401400, Taraxasteryl Acetate
MF: C₃₂H₅₂O₂
FW: 468.8
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Inula britannica*



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

Laboratory Procedures

Taraxasterol acetate is supplied as a solid. A stock solution may be made by dissolving the taraxasterol acetate in the solvent of choice, which should be purged with an inert gas. Taraxasterol acetate is soluble in the organic solvent DMSO (warmed/sonicated) at a concentration of approximately 0.1 mg/ml.

Description

Taraxasterol acetate is a pentacyclic triterpene that has been found in *P. sagittalis* and has diverse biological activities.^{1,2} It reduces the production of reactive oxygen species (ROS) and Hsp72 induced by phorbol 12-myristate 13-acetate (PMA; Item No. 10008014) in isolated human neutrophils when used at concentrations of 10 and 100 μ M and increases Hsp72 production in unstimulated isolated human neutrophils at 100 μ M.¹ Taraxasterol acetate (50 μ g/ml) reduces protein levels of the E3 ubiquitin ligase RNF31 and increases levels of the autophagy marker LC3-I/II in HCT116 and SW480 cells.² Topical administration of taraxasterol acetate (46 μ g/ear) reduces ear edema induced by arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607), croton oil, or PMA in mice.¹

References

1. Pérez-García, F., Marín, E., Parella, T., *et al.* Activity of taraxasteryl acetate on inflammation and heat shock protein synthesis. *Phytomedicine* **12**(4), 278-284 (2005).
2. Tang, C.-T., Yang, J., Liu, Z.-D., *et al.* Taraxasterol acetate targets RNF31 to inhibit RNF31/p53 axis-driven cell proliferation in colorectal cancer. *Cell Death Discov.* **7**(1), 66 (2021).

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, 12/19/2022

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