

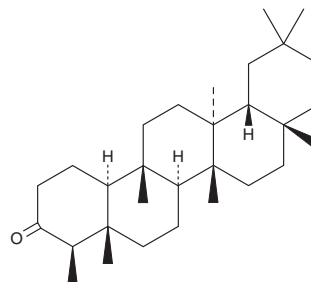
PRODUCT INFORMATION



Friedelin

Item No. 43506

CAS Registry No.: 559-74-0
Formal Name: (4 β ,5 β ,8 α ,9 β ,10 α ,13 α ,14 β)-5,9,13-trimethyl-24,25,26-trinoroleanan-3-one
Synonyms: Friedelanone, (-)-Friedelin, NSC 55141
MF: C₃₀H₅₀O
FW: 426.7
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/(*Imperata cylindrica* (L.) P. Beauv.)



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

Laboratory Procedures

Friedelin is supplied as a solid. A stock solution may be made by dissolving the friedelin in the solvent of choice, which should be purged with an inert gas. Friedelin is slightly soluble (0.1-1 mg/ml) in chloroform.

Description

Friedelin is a phytosterol that has been found in *L. eriocalyx* and has diverse biological activities.^{1,2} It scavenges DPPH (Item No. 14805) radicals in a cell-free assay (IC₅₀ = 21.1 mM).¹ Friedelin is toxic to mosquito larvae and adults (LC₅₀s = 456.99 and 891.23 μ g/ml, respectively).² It decreases the proliferation of *P. falciparum* (IC₅₀ = 250.7 ng/ml). *In vivo*, friedelin (40 mg/kg per day) reduces serum levels of glutamate dehydrogenase (GDH), alanine transaminase (ALT), and lactate dehydrogenase (Ldh) and increases hepatic levels of superoxide dismutase (SOD), catalase (Cat), glutathione (GSH), and glutathione peroxidase (Gpx) in a rat model of liver damage induced by carbon tetrachloride (CCl₄).

References

1. Manguro, L.O.A., Owuor, P.O., and Ochung, A.A. Isolation, characterization and biological activities of phytoconstituents from *Lonchocarpus eriocalyx* harms leaves. *Trends Phytochem. Res.* **2(3)**, 135-146 (2018).
2. Sunil, C., Duraipandiyar, V., Ignacimuthu, S., et al. Antioxidant, free radical scavenging and liver protective effects of friedelin isolated from *Azima tetraacantha* Lam. leaves. *Food Chem.* **139(1-4)**, 860-865 (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, 06/27/2025

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