PRODUCT INFORMATION



Epifriedelanol

Item No. 36148

CAS Registry No.: Formal Name:	16844-71-6 (3β,4β,5β,8α,9β,10α,13α,14β)-5,9,13- trimethyl-4,25,26-trinoroleanan-3-ol	
Synonyms:	Friedelan- 3β -ol, β -Friedelinol, <i>epi</i> -Friedelanol, Longan Triterpane A	H
MF:	C ₃₀ H ₅₂ O	
FW:	428.7	
Purity:	≥85%	
Supplied as:	A solid	но
Storage:	-20°C	
Stability:	≥4 years	•

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

Laboratory Procedures

Epifriedelanol is supplied as a solid. A stock solution may be made by dissolving the epifriedelanol in the solvent of choice, which should be purged with an inert gas. Epifriedelanol is slightly soluble in DMSO, dimethyl formamide, and acetonitrile.

Description

Epifriedelanol is a triterpene that has been found in U. davidiana root bark and has anticancer and senescence modifying activities.¹ It increases levels of P-glycoprotein (P-gp) and multidrug resistance protein 2 (MDR2) and enhances doxorubicin-induced cytotoxicity in doxorubicin-resistant K562/ADM cells when used at concentrations ranging from 5 to 20 μ M. It also inhibits doxorubicin-induced senescence in isolated human dermal fibroblasts (HDFs), as well as replicative senescence in HDFs and human umbilical vein endothelial cells (HUVECs).²

References

- 1. Li, Y., Liao, Z., Wei, X., et al. Epifriedelanol enhances adriamycin-induced cytotoxicity towards K562/ADM cells by down regulating of P-gp and MRP2. Xenobiotica 1-28 (2022).
- 2. Yang, H.H., Son, J.-K., Jung, B., et al. Epifriedelanol from the root bark of Ulmus davidiana inhibits cellular senescence in human primary cells. Planta Med. 77(5), 441-449 (2011).

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.

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