

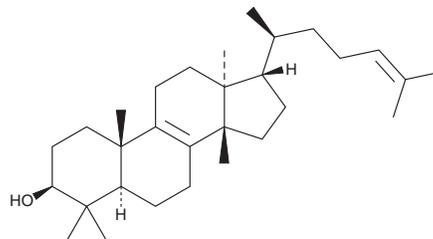
PRODUCT INFORMATION



Tirucallol

Item No. 35623

CAS Registry No.: 514-46-5
Formal Name: (3 β ,13 α ,14 β ,17 α ,20S)-lanosta-8,24-dien-3-ol
Synonyms: NSC 403164, (+)-Tirucallol
MF: C₃₀H₅₀O
FW: 426.7
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/*Euphorbia ebracteolata* Hayata



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

Laboratory Procedures

Tirucallol is supplied as a solid. A stock solution may be made by dissolving the tirucallol in the solvent of choice, which should be purged with an inert gas. Tirucallol is soluble in organic solvents such as DMSO. The solubility of tirucallol in these solvents is approximately 10 mM. It is also soluble in acetone.

Description

Tirucallol is a triterpenoid that has been found in *Camellia* and has reverse transcriptase inhibitory and anti-inflammatory activities.^{1,2} It inhibits HIV-1 reverse transcriptase (IC₅₀ = 3.8 μ M).¹ Tirucallol (25, 50, and 100 μ M) reduces nitric oxide (NO) and prostaglandin E₂ (PGE₂; Item No. 14010) production in LPS-stimulated isolated mouse peritoneal macrophages.² Topical application of tirucallol (0.25, 0.5, and 1 mg/ear) reduces ear edema induced by phorbol 12-myristate 13-acetate (TPA; Item No. 10008014) in mice.

References

1. Akihisa, T., Ogihara, J., Kato, J., *et al.* Inhibitory effects of triterpenoids and sterols on human immunodeficiency virus-1 reverse transcriptase. *Lipids* **36**(5), 507-512 (2001).
2. Fernandez-Arche, A., Saenz, M.T., Arroyo, M., *et al.* Topical anti-inflammatory effect of tirucallol, a triterpene isolated from *Euphorbia lactea* latex. *Phytomedicine* **17**(2), 146-148 (2010).

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, 12/12/2022

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