

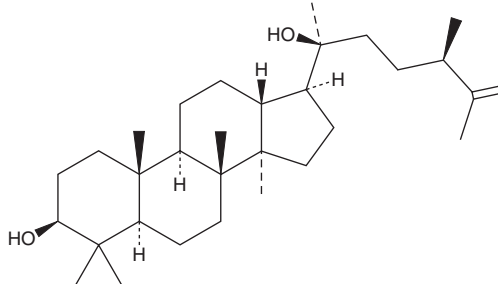
PRODUCT INFORMATION



Carnaubadiol

Item No. 29728

CAS Registry No.: 4547-30-2
Formal Name: 24R-methyl-dammar-25-ene-3 β ,20-diol
MF: C₃₁H₅₄O₂
FW: 458.8
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/*Copernicia prunifera*



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

Laboratory Procedures

Carnaubadiol is supplied as a solid. A stock solution may be made by dissolving the carnaubadiol in the solvent of choice, which should be purged with an inert gas. Carnaubadiol is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

Carnaubadiol is a triterpene that has been found in *E. corollata* and has antiprotozoal activities.^{1,2} It is active against *L. infantum* promastigotes and amastigotes, as well as *T. cruzi* trypomastigotes (IC₅₀s = 46.2, 7.8, and 14.9 μ M, respectively).²

References

1. Piatak, D.M. and Reimann, K.A. Plant investigations IV. Corolladiol A new triterpene from *euphorbia corollata*. *Tetrahedron Lett.* **13(44)**, 4525-4528 (1972).
2. de Almeida, B.C., Araújo, B.Q., Carvalho, A.A., et al. Antiprotozoal activity of extracts and isolated triterpenoids of 'carnauba' (*Copernicia prunifera*) wax from Brazil. *Pharm. Biol.* **54(12)**, 3280-3284 (2016).

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, 10/21/2022

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