

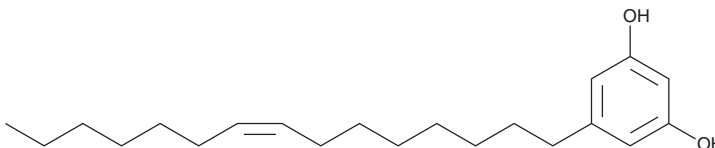
PRODUCT INFORMATION



Cardol monoene

Item No. 30033

CAS Registry No.: 22910-86-7
Formal Name: 5-(8Z)-8-pentadecenyl-1,3-benzenediol
Synonym: Bilobol
MF: $C_{21}H_{34}O_2$
FW: 318.5
Purity: $\geq 95\%$
Supplied as: A 5 mg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥ 2 years
Item Origin: Synthetic



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

Description

Cardol monoene is a phenol that has been found in cashew nut shell liquid and has diverse biological activities.¹⁻⁴ It inhibits α -glucosidase in a cell-free assay ($IC_{50} = 14.32 \text{ nM}$).² Cardol ($15 \text{ }\mu\text{g/ml}$) inhibits LPS-induced IL-6 and IL-8 production in HepG2 cells.⁴ It inhibits *T. brucei* and *P. falciparum* replication in cell-free assays ($EC_{50}\text{s} = 0.7$ and $12.4 \text{ }\mu\text{g/ml}$, respectively).³

References

1. Tyman, J.H. and Kiong, L.S. Long chain phenols: Part XI. Composition of natural cashew nutshell liquid (*Anacardium occidentale*) from various sources. *Lipids* **13**(8), 525-532 (1978).
2. He, Q., Ma, P., and Torshizi, R. The impact of some natural phenolic compounds on α -glucosidase and sorbitol dehydrogenase enzymes, and anti-leukemia cancer potential, spin density distributions, and in silico studies. *J. Oleo Sci.* **71**(6), 863-873 (2022).
3. Siheri, W., Ebiloma, G.U., Igoli, J.O., et al. Isolation of a novel flavanonol and alkylresorcinol with highly potent anti-trypanosomal activity from Libyan propolis. *Molecules* **24**(6), (2019).
4. Xu, J., Li, Y., Yang, X., et al. Bilobol inhibits the lipopolysaccharide-induced expression and distribution of RhoA in HepG2 human hepatocellular carcinoma cells. *Oncol. Lett.* **10**(2), 962-966 (2015).

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, 08/08/2025

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