

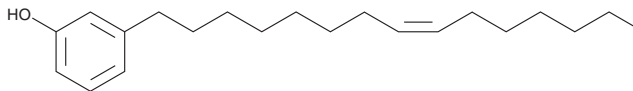
PRODUCT INFORMATION



Cardanol monoene

Item No. 23154

CAS Registry No.: 501-26-8
Formal Name: 3-8Z-pentadecen-1-yl-phenol
Synonym: Cardanol C15:1
MF: C₂₁H₃₄O
FW: 302.5
Purity: ≥98%
UV/Vis.: λ_{max}: 276 nm
Supplied as: A neat oil
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

Cardanol monoene is supplied as a neat oil. A stock solution may be made by dissolving the oil in the solvent of choice. Cardanol monoene is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of cardanol monoene in these solvents is approximately 22, 15, and 20 mg/ml, respectively.

Cardanol monoene is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, cardanol monoene should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Cardanol monoene has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Cardanol monoene is a phenol found in cashew nut shell liquid that reversibly inhibits tyrosinase with an IC₅₀ value of 56 μM *in vitro*.¹ It halts the cell cycle at the S phase and inhibits proliferation dose- and time-dependently in M14 melanoma cells (IC₅₀s = 23.15 and 12.30 μM after 24 and 48 hours of treatment, respectively).² Cardanol monoene also induces apoptosis *via* the intrinsic pathway following the accumulation of reactive oxygen species (ROS). A mixture of cardanol mono-, di-, and triene is used to synthesize cardanol-metal complexes that inhibit uropathogenic *E. coli* biofilm formation.³

References

1. Yu, X.-P., Su, W.-C., Wang, Q., *et al.* Inhibitory mechanism of cardanols on tyrosinase. *Proc. Biochem.* **51(12)**, 2330-2237 (2016).
2. Su, W.-C., Lin, Y.-F., Wang, Y.-X., *et al.* Mitochondria-associated apoptosis in human melanoma cells induced by cardanol monoene from cashew nut shell liquid. *J. Agric. Food Chem.* **65(28)**, 5620-5631 (2017).
3. Lalitha, K., Prasad, Y.S., Sridharan, V., *et al.* Intrinsic hydrophobic antibacterial thin film from renewable resources: Application in the development of anti-biofilm urinary catheters. *ACS Sus. Chem. Eng. (Univ. of Berlin)* **5(1)**, 436 (2017).

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, 09/20/2017

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