

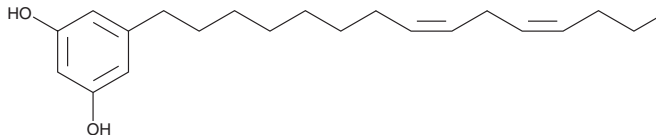
PRODUCT INFORMATION



Cardol diene

Item No. 23373

CAS Registry No.: 79473-25-9
Formal Name: 5-(8Z,11Z)-8,11-pentadecadien-1-yl-1,3-benzenediol
MF: C₂₁H₃₂O₂
FW: 316.5
Purity: ≥98%
UV/Vis.: λ_{max}: 279 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

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

Cardol diene is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, cardol diene should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Cardol diene 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

Cardol diene is a phenol found in cashew nut shell liquid.¹ It is schistosomicidal, killing 50, 100, 100, and 100% of *S. mansoni* worms after 24 hours when used at concentrations of 25, 50, 100, or 200 μM, respectively.² Cardol diene has an LC₅₀ value of 32.2 μM after 24 and 48 hours against *S. mansoni* worms. It has been used as a starting material for the synthesis of *bis*-benzoxazines.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. Alvarenga, T.A., de Oliveira, P.F., de Souza, J.M., et al. Schistosomicidal activity of alkyl-phenols from the cashew *Anacardium occidentale* against *Schistosoma mansoni* adult worms. *J. Agric. Food Chem.* **64(46)**, 8821-8827 (2016).
3. Attanasi, O.A., Behalo, M.S., Favi, G., et al. Solvent free synthesis of novel mono- and *bis*-benzoxazines from cashew nut shell liquid components. *Curr. Org. Chem.* **16(21)**, 2613-2621 (2012).

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, 10/18/2017

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