

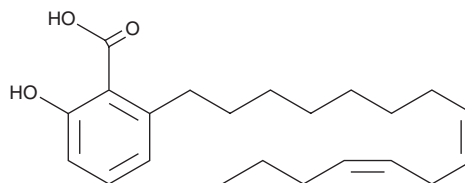
PRODUCT INFORMATION



Anacardic Acid diene

Item No. 22663

CAS Registry No.: 103904-74-1
Formal Name: 2-hydroxy-6-8Z,11Z-pentadecadien-1-yl-benzoic acid
Synonym: Anacardic Acid 15:2
MF: C₂₂H₃₂O₃
FW: 344.5
Purity: ≥95%
UV/Vis.: λ_{max}: 244, 313 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

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

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

Anacardic acid diene is a polyunsaturated form of anacardic acid (Item No. 13144) that has been found in cashew nut shell liquid.¹ It has antibacterial activity against methicillin resistant *S. aureus* (MRSA) and *S. mutans* (MICs = 12.5 and 6.25 µg/ml, respectively).² Anacardic acid diene has schistosomicidal activity against adult *S. mansoni* worms when used at a concentration of 100 µM.¹ It also inhibits soybean lipoxygenase-1 in a time-dependent manner.³

References

1. 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).
2. Kubo, I., Nihei, K., and Tsujimoto, K. Antibacterial action of anacardic acids against methicillin resistant *Staphylococcus aureus* (MRSA). *J. Agric. Food Chem.* **51(26)**, 7624-7628 (2003).
3. Ha, T.J. and Kubo, I. Lipoxygenase inhibitory activity of anacardic acids. *J. Agric. Food Chem.* **53(11)**, 4350-4354 (2005).

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, 09/12/2019

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