

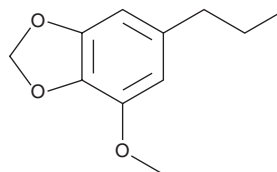
PRODUCT INFORMATION



Dihydromyristicin

Item No. 39272

CAS Registry No.: 52811-28-6
Formal Name: 4-methoxy-6-propyl-1,3-benzodioxole
MF: C₁₁H₁₄O₃
FW: 194.2
Purity: ≥95%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Dihydromyristicin is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO purged with an inert gas can be used. Dihydromyristicin is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

Description

Dihydromyristicin is a derivative of the alkenylbenzene myristicin.¹ It induces glutathione S-transferase (GST) activity in the liver and small intestinal mucosa, but not the forestomach, lungs, large intestinal mucosa, or bladder, in mice when administered at a dose of 20 mg/animal.

Reference

1. Zheng, G.-Q., Kenney, P.M., and Lam, L.K.T. Myristicin: A potential cancer chemopreventive agent from parsley leaf oil. *J. Agric. Food Chem.* **40(1)**, 107-110 (1992).

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, 01/11/2024

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