

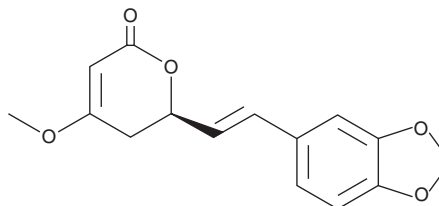
PRODUCT INFORMATION



Methysticin

Item No. 20258

CAS Registry No.: 495-85-2
Formal Name: (6R)-6-[(1E)-2-(1,3-benzodioxol-5-yl)ethenyl]-5,6-dihydro-4-methoxy-2H-pyran-2-one
Synonyms: Kavatin, (+)-Methysticin, NSC 112158
MF: C₁₅H₁₄O₅
FW: 274.3
Purity: ≥98%
UV/Vis.: λ_{max}: 208, 225, 265, 306 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Methysticin is supplied as a crystalline solid. A stock solution may be made by dissolving the methysticin in the solvent of choice, which should be purged with an inert gas. Methysticin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of methysticin in these solvents is approximately 1, 5, and 2.5 mg/ml, respectively.

Methysticin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, methysticin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Methysticin has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Methysticin is a natural kavalactone from the kava plant, *P. methysticum*. It enhances the binding of bicuculline methochloride at the γ -aminobutyric acid (GABA) receptor GABA_A at 0.1 μ M.¹ Methysticin inhibits the cytochrome P450 (CYP) isoforms CYP2C9 and CYP2D6.²

References

1. Boonen, G. and Häberlein, H. Influence of genuine kavapyrone enantiomers on the GABA_A binding site. *Planta Med.* **64(6)**, 504-506 (1998).
2. Murray, M. Toxicological actions of plant-derived and anthropogenic methylenedioxyphenyl-substituted chemicals in mammals and insects. *J. Toxicol. Environ. Health B. Crit. Rev.* **15(6)**, 365-395 (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, 09/29/2022

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