

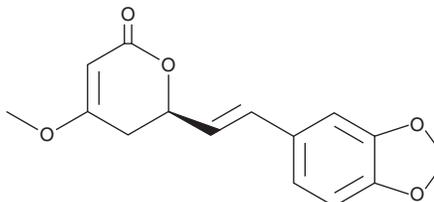
PRODUCT INFORMATION



Methysticin

Item No. 33882

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}: 227, 268 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

Methysticin (Item No. 33882) is an analytical reference standard categorized as a kavalactone.¹ This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 20258).

Reference

1. Volgin, A., Yang, L., Amstislavskaya, T., *et al.* DARK classics in chemical neuroscience: Kava. *ACS Chem. Neurosci.* **11**(23), 3893-3904 (2020).

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, 07/06/2023

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