

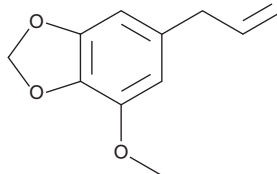
PRODUCT INFORMATION



Myristicin

Item No. 15201

CAS Registry No.: 607-91-0
Formal Name: 4-methoxy-6-(2-propen-1-yl)-1,3-benzodioxole
MF: $C_{11}H_{12}O_3$
FW: 192.2
Purity: $\geq 98\%$
UV/Vis.: $\lambda_{\max}$: 212 nm
Supplied as: A 10 mg/ml 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.

Description

Myristicin (Item No. 15201) is an alkenylbenzene present in small amounts in the essential oil of nutmeg that is reported to act as a serotonin receptor antagonist, a weak monamine oxidase inhibitor, and to produce hallucinogenic effects.¹ Abuse of myristicin has led to fatal poisoning, which has prompted improved methods for HPLC determination of myristicin in human plasma.² This product is intended for forensic and research purposes.

References

1. Kalbhen, D.A. Nutmeg as a narcotic. A contribution to the chemistry and pharmacology of nutmeg. *Angew. Chem. Int. Ed. Engl.* **10(6)**, 370-374 (1971).
2. Dawidowicz, A.L. and Dybowski, M.P. Simple and rapid determination of myristicin in human serum. *Forensic Toxicol.* **31(1)**, 119-123 (2013).

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, 01/14/2026

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