

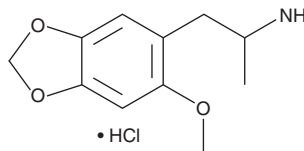
PRODUCT INFORMATION



MMDA-2 (hydrochloride)

Item No. 17659

CAS Registry No.: 64778-82-1
Formal Name: 6-methoxy- α -methyl-1,3-benzodioxole-5-ethanamine, monohydrochloride
MF: C₁₁H₁₅NO₃ • HCl
FW: 245.7
Purity: ≥95%
UV/Vis.: λ_{max} : 238, 300 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

MMDA-2 (hydrochloride) (Item No. 17659) is an analytical reference standard that is classified as an amphetamine. It only weakly induces the release of serotonin or dopamine from rat brain synaptosomes and does not produce amphetamine-like responses in drug discrimination studies in rats.^{1,2} Instead, MMDA-2 is thought to act as a 5-HT_{2A} agonist, which may account for hallucinogenic effects.³ This product is intended for forensic and research applications.

References

1. McKenna, D.J., Guan, X.-M., and Shulgin, A.T. 3,4-Methylenedioxyamphetamine (MDA) analogues exhibit differential effects on synaptosomal release of ³H-dopamine and ³H-5-hydroxytryptamine. *Pharmacol. Biochem. Behav.* **38(3)**, 505-512 (1991).
2. Glennon, R.A., Yousif, M., Naiman, N., *et al.* Methcathinone: A new and potent amphetamine-like agent. *Pharmacol. Biochem. Behav.* **26(3)**, 547-551 (1987).
3. Clare, B.W. QSAR of benzene derivatives: Comparison of classical descriptors, quantum theoretic parameters and flip regression, exemplified by phenylalkylamine hallucinogens. *J. Comput. Aided Mol. Des.* **16(8-9)**, 611-633 (2002).

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