

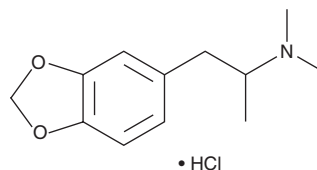
PRODUCT INFORMATION



MDDMA (hydrochloride)

Item No. 14217

CAS Registry No.: 74341-79-0
Formal Name: N,N,α-trimethyl-1,3-benzodioxole-5-ethanamine, monohydrochloride
Synonym: 3,4-Methylenedioxydimethylamphetamine
MF: C₁₂H₁₇NO₂ • HCl
FW: 243.7
Purity: ≥98%
UV/Vis.: λ_{max}: 236, 287 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

MDDMA is a centrally active substituted methylenedioxyphenethylamine that has appeared in illicit drug samples.^{1,2} It is structurally similar to 3,4-MDMA (Item No. 13971) with an additional methyl group that is predicted to result in relatively reduced central activity compared to 3,4-MDMA.³ This product is intended for research and forensic applications.

References

1. Braun, U., Shulgin, A.T., and Braun, G. Centrally active N-substituted analogs of 3,4-methylenedioxyphenylisopropylamine (3,4-methylenedioxyamphetamine). *J. Pharm. Sci.* **69(2)**, 192-195 (1980).
2. Thigpen, A., DeRuiter, J., and Clark, C.R. GC-MS studies on the regioisomeric 2,3- and 3,4-methylenedioxyphenethylamines related to MDEA, MDMMA, and MBDB. *J. Chromatogr. Sci.* **45(5)**, 229-235 (2007).
3. Dal Cason, T.A. An evaluation of the potential for clandestine manufacture of 3,4-methylenedioxyamphetamine (MDA) analogs and homologs. *J. Forensic Sci.* **35(3)**, 675-697 (1990).

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, 06/15/2023

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