

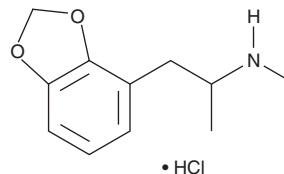
PRODUCT INFORMATION



2,3-MDMA (hydrochloride)

Item No. 13970

CAS Registry No.: 168968-01-2
Formal Name: N, α -dimethyl-1,3-benzodioxole-4-ethanamine, monohydrochloride
Synonym: 2,3-Methylenedioxyamphetamine
MF: C₁₁H₁₅NO₂ • HCl
FW: 229.7
Purity: \geq 98%
UV/Vis.: λ_{\max} : 234, 283 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 5 years



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

Description

2,3-MDMA is an aromatic positional isomer of 3,4-MDMA (ecstasy, Item No. 13971), a euphoric entactogen that is illegal in most countries and is regulated (Schedule I) in the United States.^{1,2} While 2,3-MDMA exhibits a similar potency to 3,4-MDMA for inhibition of the noradrenalin transporter (IC₅₀s = 6.2 and 6.6 μ M, respectively), it is less potent than 3,4-MDMA for inhibition of the serotonin transporter (IC₅₀s = 82 μ M and 34.8 μ M, respectively).³ This compound is intended for forensic and research purposes only.

References

1. Casale, J.F., Hays, P.A., and Klein, R.F.X. Synthesis and characterization of the 2,3-methylenedioxyamphetamines. *J. Forensic Sci.* **40(3)**, 391-400 (1995).
2. Chakraborty, K., Neogi, R., and Basu, D. Club drugs: Review of the 'rave' with a note of concern for the Indian scenario. *Indian J. Med. Res.* **133(6)**, 594-604 (2011).
3. Montgomery, T., Buon, C., Eibauer, S., et al. Comparative potencies of 3,4-methylenedioxyamphetamine (MDMA) analogues as inhibitors of [³H]noradrenaline and [³H]5-HT transport in mammalian cell lines. *Br. J. Pharmacol.* **152(7)**, 1121-1130 (2007).

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, 07/25/2023

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