

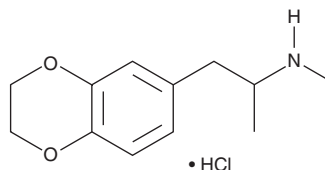
PRODUCT INFORMATION



3,4-EDMA (hydrochloride)

Item No. 15164

CAS Registry No.: 2748304-41-6
Formal Name: 2,3-dihydro-N, α -dimethyl-1,4-benzodioxin-6-ethanamine, monohydrochloride
Synonyms: 3,4-EDO-N-methylamphetamine, 3,4-Ethylenedioxy-N-methylamphetamine
MF: C₁₂H₁₇NO₂ • HCl
FW: 243.7
Purity: \geq 98%
UV/Vis.: λ_{max} : 284 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

3,4-MDMA (Item No. 13971) is a euphoric entactogen that is regulated as a Schedule I compound in the United States. 3,4-EDMA is an analog of 3,4-MDMA that stimulates the release of serotonin (Item No. 14332) and dopamine from rat brain synaptosomes at 1 μ M.¹ This product is intended for forensic and research applications.

Reference

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).

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, 06/13/2023

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