

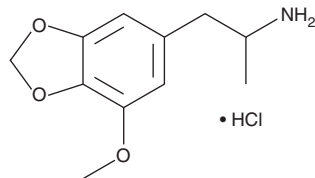
# PRODUCT INFORMATION



## MMDA (hydrochloride)

Item No. 17658

**Formal Name:** 7-methoxy- $\alpha$ -methyl-1,3-benzodioxole-5-ethanamine, monohydrochloride  
**Synonyms:** 3-MeO MDA, 3-methoxy MDA  
**MF:** C<sub>11</sub>H<sub>15</sub>NO<sub>3</sub> • HCl  
**FW:** 245.7  
**Purity:** ≥98%  
**Supplied as:** A neat solid  
**Storage:** -20°C  
**Stability:** ≥5 years



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

### Description

3,4-MDA (Item No. ISO60189 | 11554) is a psychotropic amphetamine derivative that acts as a serotonergic 5-HT<sub>2A</sub> receptor agonist and releases monoamines by interacting with monoamine transporters.<sup>1-3</sup> MMDA is a substituted form of 3,4-MDA bearing a methoxy moiety at the 3 position of the phenyl group. It is at least as potent as 3,4-MDA at promoting monoamine release from synaptosomes.<sup>4</sup> This product is intended for forensic and research applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

### References

1. Lyon, R.A., Glennon, R.A., and Titeler, M. 3,4-Methylenedioxyamphetamine (MDMA): Stereoselective interactions at brain 5-HT<sub>1</sub> and 5-HT<sub>2</sub> receptors. *Psychopharmacology* **88(4)**, 525-526 (1986).
2. Baggott, M.J., Siegrist, J.D., Galloway, G.P., *et al.* Investigating the mechanisms of hallucinogen-induced visions using 3,4-methylenedioxyamphetamine (MDA): A randomized controlled trial in humans. *PLoS One* **5(12)**, 1-13 (2010).
3. Kaizaki, A., Tanaka, S., Tsujikawa, K., *et al.* Recreational drugs, 3,4-methylenedioxyamphetamine (MDMA), 3,4-methylenedioxyamphetamine (MDA) and diphenylprolinol, inhibit neurite outgrowth in PC12 cells. *J. Toxicol. Sci.* **35(3)**, 375-381 (2010).
4. 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, 07/01/2019

#### 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