

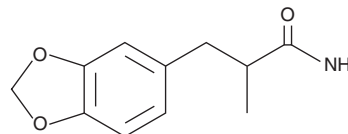
# PRODUCT INFORMATION



## MDA 2-amido analog

Item No. 9002291

**CAS Registry No.:** 858215-05-1  
**Formal Name:**  $\alpha$ -methyl-1,3-benzodioxole-5-propanamide  
**Synonyms:** 3-(1,3-Benzodioxol-5-yl)-2-methylpropanamide,  $\alpha$ -methyl-3,4-Methylenedioxyphenylpropanamide, MMDPPA  
**MF:**  $C_{11}H_{13}NO_3$   
**FW:** 207.2  
**Purity:**  $\geq 98\%$   
**Supplied as:** A neat solid  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 1$  year



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

### Description

3,4-Methylenedioxyamphetamine (MDA; Item Nos. ISO60189 | 11554) is a psychedelic drug that is regulated in the United States. MDA 2-amido analog is an intermediate produced in the synthesis of MDA from helional (also known as ocean propanal, aquanal, or tropional).<sup>1</sup> Also known as MMDPPA, MDA 2-amido analog can also be used as a direct precursor in the synthesis of MDA.<sup>2</sup> The physiological and toxicological properties of this compound are not known. 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. Kovacs, E.J.I. and Kirby, D.A. Manufacture of 3,4-methylenedioxyamphetamine from helional using Beckmann and Hofmann rearrangements. *J. Clan. Lab. Inv. Chem.* **23(1)**, 5-14 (2013).
2. Dal Cason, T.A., Corbett, C.A., Poole, P.K., et al. An unusual clandestine laboratory synthesis of 3,4-methylenedioxyamphetamine (MDA). *Forensic Sci. Int.* **223**, 279-291 (2012).

#### 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, 07/11/2019

#### 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