

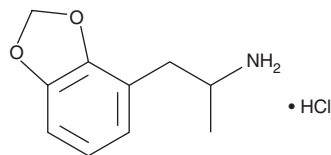
# PRODUCT INFORMATION



## 2,3-MDA (hydrochloride)

Item No. 13969

**CAS Registry No.:** 86029-48-3  
**Formal Name:**  $\alpha$ -methyl-1,3-Benzodioxole-4-ethanamine, monohydrochloride  
**Synonym:** 2,3-Methylenedioxyamphetamine  
**MF:**  $C_{10}H_{13}NO_2 \cdot HCl$   
**FW:** 215.7  
**Purity:**  $\geq 98\%$   
**UV/Vis.:**  $\lambda_{max}$ : 234, 283 nm  
**Supplied as:** A crystalline solid  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 5$  years



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

### Description

3,4-Methylenedioxyamphetamine (MDA) is psychedelic and entactogenic drug that is scheduled as a controlled substance in the United States.<sup>1</sup> 2,3-MDA is a structural isomer of MDA which differs only in the placement of the methylenedioxy group. The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

### Reference

1. 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).

#### 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