

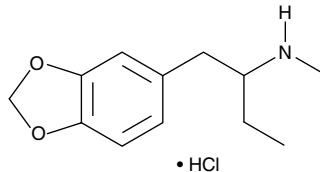
# Product Information



## MBDB (hydrochloride)

Item No. 14218

**CAS Registry No.:** 128767-12-4  
**Formal Name:**  $\alpha$ -ethyl-N-methyl-1,3-benzodioxole-5-ethanamine, monohydrochloride  
**Synonym:** N-Methyl-1,3-benzodioxolylbutanamine  
**MF:**  $C_{12}H_{17}NO_2 \cdot HCl$   
**FW:** 243.7  
**Purity:**  $\geq 95\%$   
**Stability:**  $\geq 2$  years at  $-20^\circ C$   
**Supplied as:** A crystalline solid  
**UV/Vis.:**  $\lambda_{max}$ : 236, 288 nm



### Laboratory Procedures

For long term storage, we suggest that MBDB (hydrochloride) be stored as supplied at  $-20^\circ C$ . It should be stable for at least two years.

MBDB (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the MBDB (hydrochloride) in the solvent of choice. MBDB (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of MBDB (hydrochloride) in these solvents is approximately 16, 10, and 10 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of MBDB (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of MBDB (hydrochloride) in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

MBDB is an analog of methylenedioxymethamphetamine (MDMA, Item No. 13971), differing only by the substitution of an ethyl for methyl group at the  $\alpha$ -carbon. Like MDMA, MBDB is an entactogen that potently inhibits monoamine uptake.<sup>1,2</sup> This product is intended for forensic and research purposes.

### References

- Nichols, D.E., Brewster, W.K., Johnson, M.P., *et al.* Nonneurotoxic tetralin and indan analogues of 3,4-(methylenedioxy)amphetamine (MDA). *J. Med. Chem.* **33**(2), 703-710 (1990).
- Nagai, F., Nonaka, R., and Satoh Hisashi Kamimura, K. The effects of non-medically used psychoactive drugs on monoamine neurotransmission in rat brain. *Eur. J. Pharmacol.* **559**(2-3), 132-137 (2007).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/14218](http://www.caymanchem.com/catalog/14218)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 11/20/2014

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)