

PRODUCT INFORMATION



MBDB (hydrochloride)

Item No. 14218

CAS Registry No.: 128767-12-4

Formal Name: a-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

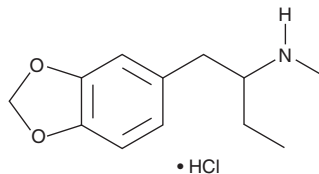
Purity: $\geq 95\%$

UV/Vis.: λ_{max} : 236, 288 nm

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 5 years



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

Description

MBDB is an analog of methylenedioxymethamphetamine (MDMA, Item No. 13971), differing only by the substitution of an ethyl for methyl group at the α -carbon. Like MDMA, MBDB is an entactogen that potently inhibits monoamine uptake.^{1,2} This product is intended for forensic and research purposes.

References

1. 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).
2. 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).

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, 06/16/2023

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