

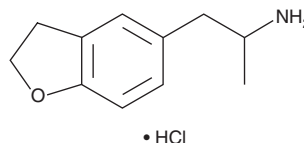
PRODUCT INFORMATION



5-APDB (hydrochloride)

Item No. 11456

CAS Registry No.: 152623-94-4
Formal Name: 2,3-dihydro- α -methyl-5-benzofuranethanamine, monohydrochloride
Synonyms: 5-(2-Aminopropyl)-2,3-dihydrobenzofuran, EMA-4, 3-desoxy-MDA
MF: C₁₁H₁₅NO • HCl
FW: 213.7
Purity: \geq 98%
Stability: \geq 2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max} : 231, 287 nm



Laboratory Procedures

For long term storage, we suggest that 5-APDB (hydrochloride) be stored as supplied at -20°C. It should be stable for at least two years.

5-APDB (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the 5-APDB (hydrochloride) in the solvent of choice. 5-APDB (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of 5-APDB (hydrochloride) in these solvents is approximately 20 mg/ml.

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 5-APDB (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 5-APDB (hydrochloride) in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

5-APDB is an entactogen belonging to the amphetamine and the phenethylamine classes. It is an analog of MDA where the heterocyclic 3-position oxygen from the 3,4-methylenedioxy ring has been replaced by a methylene group. 5-APDB can inhibit serotonin, dopamine and norepinephrine reuptake with IC₅₀ values of 0.13, 7.1, and 3.2 μ M, respectively, and generalizes closely to the non-stimulant MDMA analogs, MBDB and MMAI, in drug discrimination studies.¹ This product is intended for forensic purposes.

Reference

1. Monte, A.P., Marona-Lewicka, D., Cozzi, N.V., *et al.* Synthesis and pharmacological examination of benzofuran, indan, and tetralin analogues of 3,4-(methylenedioxy)amphetamine. *J. Med. Chem.* **36**, 3700-3706 (1993).

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, 10/08/2015

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