

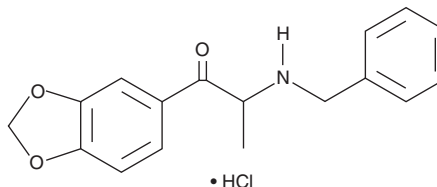
PRODUCT INFORMATION



3,4-Methylenedioxy-N-benzylcathinone (hydrochloride)

Item No. 9001330

CAS Registry No.: 1823274-68-5
Formal Name: 1-(1,3-benzodioxol-5-yl)-2-[(phenylmethyl)amino]-1-propanone, monohydrochloride
Synonym: BMDP
MF: C₁₇H₁₇NO₃ • HCl
FW: 319.8
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 204, 234, 281, 319 nm



Laboratory Procedures

For long term storage, we suggest that 3,4-methylenedioxy-N-benzylcathinone (BMDP) (hydrochloride) be stored as supplied at -20°C. It should be stable for at least two years.

BMDP (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the BMDP (hydrochloride) in the solvent of choice. BMDP (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of BMDP (hydrochloride) in ethanol is approximately 3 mg/ml and approximately 5 mg/ml in DMSO and DMF.

BMDP (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BMDP (hydrochloride) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. BMDP (hydrochloride) has a solubility of approximately 0.13 mg/ml in a 1:7 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Substituted cathinones are psychoactive compounds sold in packets labeled as bath salts or plant food.^{1,2} BMDP is a substituted cathinone that is structurally analogous to methylone, a narcotic that is regulated in the United States. This compound has a phenyl group added to the amino methyl group, substantially decreasing its solubility in aqueous solutions. The physiological and toxicological properties of this compound are not known. This product is intended for research and forensic applications.

References

1. Coppola, M. and Mondola, R. 3,4-Methylenedioxypyrovalerone (MDPV): Chemistry, pharmacology and toxicology of a new designer drug of abuse marketed online. *Toxicol. Lett.* **208(1)**, 12-15 (2012).
2. Prosser, J.M. and Nelson, L.S. The toxicology of bath salts: A review of synthetic cathinones. *J. Med. Toxicol.* **8(1)**, 33-42 (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, 01/04/2016

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