

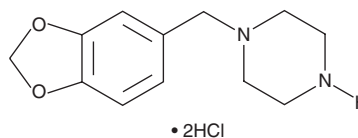
PRODUCT INFORMATION



MDBP (hydrochloride)

Item No. 19215

CAS Registry No.: 38063-96-6
Formal Name: 1-(1,3-benzodioxol-5-ylmethyl)-piperazine, dihydrochloride
Synonyms: 1-(3,4-Methylenedioxybenzyl)piperazine, 1-Piperonylpiperazine
MF: C₁₂H₁₆N₂O₂ • 2HCl
FW: 293.2
Purity: ≥95%
UV/Vis.: λ_{max}: 238, 285 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

MDBP (hydrochloride) (Item No. 19215) is an analytical reference standard that is structurally classified as a piperazine. Its metabolism and toxicological analysis in rat urine using GC-MS has been reported.¹ This product is intended for research and forensic purposes.

Reference

1. Stasck, R.F., Maurer, H.H. New designer drug 1-(3,4-methylenedioxybenzyl)piperazine (MDBP): Studies on its metabolism and toxicological detection in rat urine using gas chromatography/mass spectrometry. *J. Mass. Spectrom.* **30**, 255-261 (2004).

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, 06/19/2023

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