

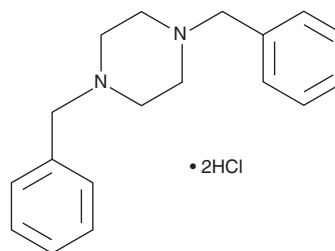
PRODUCT INFORMATION



1,4-Dibenzylpiperazine (hydrochloride)

Item No. 11206

CAS Registry No.: 2298-55-7
Formal Name: 1,4-bis(phenylmethyl)-piperazine, dihydrochloride
Synonym: DBZP
MF: C₁₈H₂₂N₂ • 2HCl
FW: 339.3
Purity: ≥98%
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

Benzylpiperazines (BZP) are stimulant drugs which inhibit monoamine uptake.^{1,2} BZP and substituted derivatives, such as 1-(3,4-methylenedioxybenzyl)piperazine, have been identified as designer drugs or drugs of abuse.^{3,4} 1,4-Dibenzylpiperazine (hydrochloride) is a common impurity produced as a reaction byproduct during the synthesis of BZP. The biological and toxicological properties of this compound are not known. This product is intended for research and forensic applications.

References

1. Baumann, M.H., Clark, R.D., Budzynski, A.G., *et al.* N-substituted piperazines abused by humans mimic the molecular mechanism of 3,4-methylenedioxymethamphetamine (MDMA, or 'Ecstasy'). *Neuropsychopharmacology* **30(3)**, 550-560 (2005).
2. Hill, S.L. and Thomas, S.H. Clinical toxicology of newer recreational drugs. *Clin. Toxicol. (Phila)* **49(8)**, 705-719 (2011).
3. Kikura-Hanajiri, R., Uchiyama, N., and Goda, Y. Survey of current trends in the abuse of psychotropic substances and plants in Japan. *Leg. Med. (Tokyo)* **13(3)**, 109-115 (2011).
4. Arbo, M.D., Bastos, M.L., and Carmo, H.F. Piperazine compounds as drugs of abuse. *Drug Alcohol Depend.* **122(3)**, 174-185 (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, 07/07/2023

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