

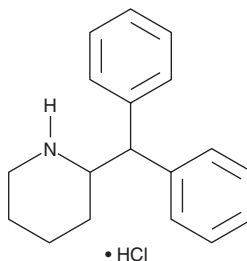
PRODUCT INFORMATION



2-DPMP (hydrochloride)

Item No. 11481

CAS Registry No.: 5807-81-8
Formal Name: 2-(diphenylmethyl)-piperidine, monohydrochloride
Synonyms: Desoxypipradrol, 2-Diphenylmethylpiperidine
MF: C₁₈H₂₁N • HCl
FW: 287.8
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

2-DPMP (hydrochloride) is structurally related to pipradrol and methylphenidate (Ritalin), which are psychostimulatory piperadines that inhibit monoamine transporters, preventing the reuptake of dopamine and norepinephrine.^{1,2} This compound has recently been identified in recreational drugs and considered the likely source of toxicity associated with prolonged psychiatric symptoms.³⁻⁵ This product is intended for research and forensic applications.

References

1. Berends, H.I., Nijlant, J.M.M., Movig, K.L.L., et al. *Eur. J. Phys. Rehabil. Med.* **45(4)**, 621-630 (2009).
2. Heron, C., Costentin, J., and Bonnet, J.-J. *Eur. J. Pharmacol.* **264**, 391-398 (1994).
3. Hill, S.L. and Thomas, S.H. *Clin. Toxicol. (Phila)* **49**, 705-719 (2011).
4. Murray, D.B., Potts, S., Haxton, C., et al. *Clin. Toxicol. (Phila)* **50**, 108-113 (2012).
5. Gibbons, S. *Clin. Toxicol. (Phila)* **50**, 15-24 (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, 07/17/2023

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