

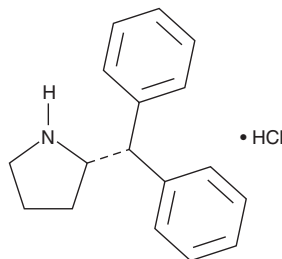
PRODUCT INFORMATION



(S)-Desoxy-D2PM (hydrochloride)

Item No. 90010945

CAS Registry No.: 188398-87-0
Formal Name: 2-(diphenylmethyl)-pyrrolidine, monohydrochloride
Synonym: (S)-2-Diphenylmethylpyrrolidine
MF: C₁₇H₁₉N • HCl
FW: 404.4
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

(S)-Desoxy-D2PM (hydrochloride) is most commonly used as a chiral solvating agent for NMR analysis of chiral compounds.¹ Interestingly, it is also, structurally, closely related to 2-DPMP (Item No. 11481), a dopamine transporter inhibitor and psychoactive stimulant.² The physiological and toxicological properties of (S)-desoxy-D2PM have not been explored.

Reference

1. Bailey, D.J., O'Hagan, D., and Tavasli, M. A short synthesis of (S)-2-(diphenylmethyl)pyrrolidine, a chiral solvating agent for NMR analysis. *Tetrahedron Asymmetry* **8**(1), 149-153 (1997).
2. Schmitt, K.C., Zhen, J., Kharkar, P., *et al.* Interaction of cocaine-, benztropine-, and GBR12909-like compounds with wildtype and mutant human dopamine transporters: Molecular features that differentially determine antagonist binding properties. *J. Neurochem.* **107**(4), 928-940 (2008).

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, 06/23/2023

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