

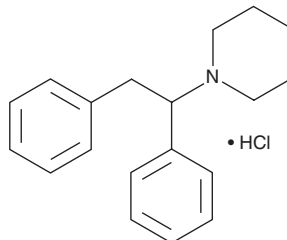
PRODUCT INFORMATION



Diphenidine (hydrochloride)

Item No. 15518

CAS Registry No.: 28383-15-5
Formal Name: 1-(1,2-diphenylethyl)-piperidine, monohydrochloride
Synonyms: DEP, 1-(1,2-diphenylethyl) piperidine
MF: C₁₉H₂₃N • HCl
FW: 301.9
Purity: ≥98%
UV/Vis.: λ_{max}: 258 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

Diphenidine (hydrochloride) (Item No. 15518) is an analytical reference standard categorized as a dissociative anesthetic.¹ Diphenidine has been found in samples seized by law enforcement and has a high potential for abuse.^{1,2} This product is intended for research and forensic applications.

References

1. Gerace, E., Bovetto, E., and DiCorcia, D. A case of nonfatal intoxication associated with the recreational use of diphenidine. *J. Forensic Sci.* **62(4)**, 1107-1111 (2017).
2. Jones, L.E., Stewart, A., Peters, K.L., et al. Infrared and raman screening of seized novel psychoactive substances: A large scale study of >200 samples. *Analyst* **141(3)**, 902-909 (2016).

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