

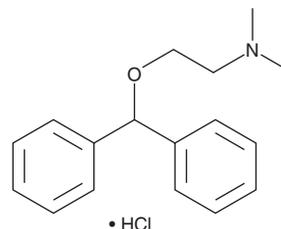
PRODUCT INFORMATION



Diphenhydramine (hydrochloride)

Item No. 22489

CAS Registry No.: 147-24-0
Formal Name: 2-(diphenylmethoxy)-N,N-dimethylethanamine, monohydrochloride
Synonyms: Diphenhydramine, DPH
MF: C₁₇H₂₁NO • HCl
FW: 291.8
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

Diphenhydramine (hydrochloride) (Item No. 22489) is an analytical reference standard that is categorized as an antihistamine.¹ It increases hallucinations and manic behavior associated with use of products marketed as bath salts.² This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 11158).

References

1. Thurmond, R.L., Gelfand, E.W., and Dunford, P.J. The role of histamine H₁ and H₄ receptors in allergic inflammation: The search for new antihistamines. *Nat. Rev. Drug Discov.* **7**(1), 41-53 (2008).
2. Gunderson, E.W., Kirkpatrick, M.G., Willing, L.M., et al. Intranasal substituted cathinone "bath salts" psychosis potentially exacerbated by diphenhydramine. *J. Addict. Med.* **7**(3), 163-168 (2013).

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, 06/30/2023

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