

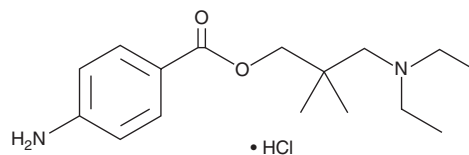
PRODUCT INFORMATION



Dimethocaine (hydrochloride)

Item No. 11159

CAS Registry No.: 553-63-9
Formal Name: 3-(diethylamino)-2,2-dimethyl-1-(4-aminobenzoate)-1-propanol, monohydrochloride
MF: C₁₆H₂₆N₂O₂ • HCl
FW: 314.9
Purity: ≥98%
UV/Vis.: λ_{max}: 222, 297 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

Dimethocaine is a local anesthetic that, because of its similarity in action to cocaine, has potential for abuse. This compound completely inhibits dopamine uptake in rat striatal synaptosomes with an IC₅₀ value of 1.2 μM comparable to that of cocaine (IC₅₀ = 0.7 μM).¹ As a result, dimethocaine dose-dependently substitutes for cocaine in drug discrimination tests in rats and rhesus monkeys.^{2,3} This product is intended for forensic and research purposes.

References

1. Woodward, J.J., Compton, D.M., Balster, R.L., *et al.* In vitro and in vivo effects of cocaine and selected local anesthetics on the dopamine transporter. *Eur. J. Pharmacol.* **277(1)**, 7-13 (1995).
2. Graham, J.H. and Balster, R.L. Cocaine-like discriminative stimulus effects of procaine, dimethocaine and lidocaine in rats. *Psychopharmacology (Berl)* **110(3)**, 287-294 (1993).
3. Wilcox, K.M., Kimmel, H.L., Lindsey, K.P., *et al.* In vivo comparison of the reinforcing and dopamine transporter effects of local anesthetics in rhesus monkeys. *Synapse* **58(4)**, 220-228 (2005).

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, 07/13/2023

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