

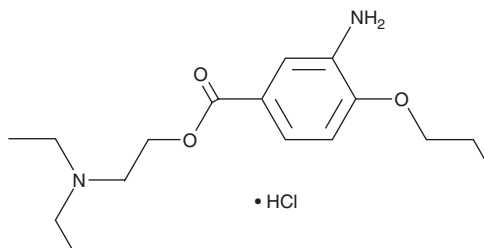
PRODUCT INFORMATION



Proxymetacaine (hydrochloride)

Item No. 39840

CAS Registry No.: 5875-06-9
Formal Name: 3-amino-4-propoxy-benzoic acid, 2-(diethylamino)ethyl ester, monohydrochloride
Synonym: Proparacaine
MF: $C_{16}H_{26}N_2O_3 \cdot HCl$
FW: 330.9
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Proxymetacaine (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the proxymetacaine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Proxymetacaine (hydrochloride) is slightly soluble (0.1-1 mg/ml) in chloroform.

Description

Proxymetacaine is an anesthetic.¹ Intrathecal administration of proxymetacaine reduces motor function, decreases righting behaviors in the placement test, and increases latency to paw withdrawal in the hot plate test in rats (ED_{50} s = 0.24, 0.3, and 0.16 μ mol/kg, respectively). Topical application of proxymetacaine (0.5% w/v) reduces corneal sensitivity in the corneal touch test in dogs.² Formulations containing proxymetacaine have been used to induce anesthesia prior to ocular surgery.

References

1. Hung, C.-H., Wang, J.-J., Chen, Y.-C., *et al.* Intrathecal oxybuprocaine and proxymetacaine produced potent and long-lasting spinal anesthesia in rats. *Neurosci. Lett.* **454**(3), (2009).
2. Herring, I.P., Bobofchak, M.A., Landry, M.P., *et al.* Duration of effect and effect of multiple doses of topical ophthalmic 0.5% proparacaine hydrochloride in clinically normal dogs. *Am. J. Vet. Res.* **66**(1), 77-80 (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, 07/18/2024

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