

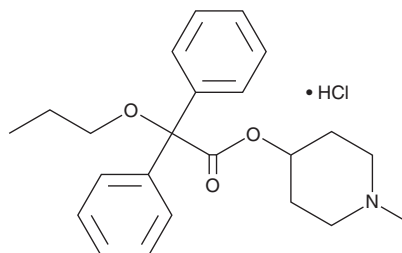
PRODUCT INFORMATION



Propiverine (hydrochloride)

Item No. 36049

CAS Registry No.: 54556-98-8
Formal Name: α -phenyl- α -propoxy-benzeneacetic acid, 1-methyl-4-piperidinyl ester, monohydrochloride
MF: $C_{23}H_{29}NO_3 \cdot HCl$
FW: 403.9
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

Propiverine (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the propiverine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Propiverine (hydrochloride) is slightly soluble in ethanol, DMSO, and dimethyl formamide.

Description

Propiverine is an antagonist of muscarinic acetylcholine receptors (mAChRs).¹ It binds to M_1 and M_2 mAChRs in rat bladder and submaxillary gland, respectively ($K_{iS} = 0.3$ and $0.19 \mu M$, respectively), and inhibits carbachol-induced contraction of bladder strips isolated from patients with normal bladder activity and detrusor overactivity associated with benign prostatic hyperplasia ($pA_{2S} = 7.7$ and 7.6 , respectively).² Formulations containing propiverine have been used in the treatment of overactive bladder.

References

1. Yamada, S., Ito, Y., Taki, Y., *et al.* The N-oxide metabolite contributes to bladder selectivity resulting from oral propiverine: Muscarinic receptor binding and pharmacokinetics. *Drug Metab. Dispos.* **38(8)**, 1314-1321 (2010).
2. Yamanishi, T., Kaga, K., Fuse, M., *et al.* The role of muscarinic receptor subtypes on carbachol-induced contraction of normal human detrusor and overactive detrusor associated with benign prostatic hyperplasia. *J. Pharmacol. Sci.* **128(2)**, 65-70 (2015).

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, 12/19/2022

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