

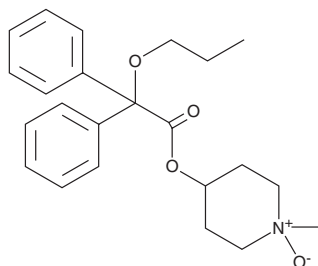
PRODUCT INFORMATION



Propiverine N-oxide

Item No. 41013

CAS Registry No.: 111071-96-6
Formal Name: α -phenyl- α -propoxy-benzeneacetic acid, 1-methyl-1-oxido-4-piperidinyl ester
MF: $C_{23}H_{29}NO_4$
FW: 383.5
Purity: $\geq 95\%$
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

Propiverine N-oxide is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Propiverine N-oxide is soluble in methanol and acetonitrile.

Description

Propiverine N-oxide is an active metabolite of the muscarinic acetylcholine receptor antagonist propiverine (Item No. 36049).¹ *In vivo*, propiverine N-oxide (0.422 mg/kg) reduces sacral anterior root stimulation-induced bladder contractions and salivary flow in mini pigs.

Reference

1. Scheepe, J.R., Braun, P.M., Jünemann, K.P., *et al.* Effects of propiverine and its metabolite propiverine-N-oxide on bladder contraction and salivation in mini pigs. *Urol. Int.* **81**(4), 468-473 (2008).

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, 11/18/2024

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