

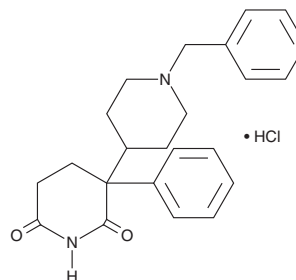
PRODUCT INFORMATION



Benzetimide (hydrochloride)

Item No. 34527

CAS Registry No.: 5633-14-7
Formal Name: 3-phenyl-1'-(phenylmethyl)-[3,4'-bipiperidine]-2,6-dione, monohydrochloride
Synonyms: Dioxatrine, R 4929, Spasmentral
MF: C₂₃H₂₆N₂O₂ • HCl
FW: 398.9
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Benzetimide (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the benzetimide (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Benzetimide (hydrochloride) is slightly soluble in chloroform (sonicated).

Benzetimide (hydrochloride) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Benzetimide is a muscarinic acetylcholine receptor (mAChR) antagonist.¹ It induces mydriasis and inhibits pilocarpine-induced salivation and lacrimation in rats (ED₅₀s = 0.06 and 0.04 mg/kg, respectively).²

References

1. Gray, J.A., Lüllmann, H., Mitchelson, F.J., *et al.* Stereoselective binding in cardiac tissue of the enantiomers of benzetimide, and antimuscarinic drug. *Br. J. Pharmacol.* **56(4)**, 485-490 (1976).
2. Janssen, P.A.J. and Niemegeers, C.J.E. The peripheral and central anticholinergic properties of benzetimide (R 4929) and other atropine-like drugs as measured in a new anti-pilocarpine test in rats. *Psychopharmacologia* **11(3)**, 231-254 (1967).

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, 11/16/2022

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