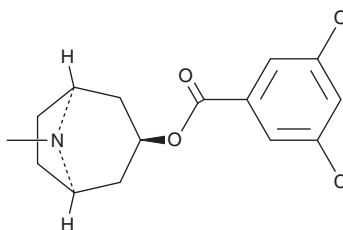


PRODUCT INFORMATION



MDL 72222
Item No. 43507

CAS Registry No.: 40796-97-2
Formal Name: 3,5-dichloro-benzoic acid, (3-endo)-8-methyl-8-azabicyclo[3.2.1]oct-3-yl ester
Synonym: Bemisetron
MF: C₁₅H₁₇Cl₂NO₂
FW: 314.2
Purity: ≥98%
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

MDL 72222 is supplied as a solid. A stock solution may be made by dissolving the MDL 72222 in the solvent of choice, which should be purged with an inert gas. MDL 72222 is slightly soluble (0.1-1 mg/ml) in ethanol.

Description

MDL 72222 is an antagonist of the serotonin (5-HT) receptor subtype 5-HT₃ (IC₅₀ = 0.82 nM).¹ It is selective for 5-HT₃ over 5-HT₁ and 5-HT₂ receptors and α₁-, α₂-, and β-adrenergic receptors (IC₅₀s = >10,000 nM for all), as well as the histamine H₁ receptor (IC₅₀ = 3,800 nM). Intraplantar administration of MDL 72222 (10 and 100 μg/animal) inhibits formalin-induced paw licking and complete Freund's adjuvant-induced decreases in paw withdrawal latency in mice.² It inhibits emesis induced by cisplatin (Item No. 13119) in dogs when administered at a dose of 0.5 mg/kg.³

References

1. Fozard, J.R. MDL 72222: A potent and highly selective antagonist at neuronal 5-hydroxytryptamine receptors. *Naunyn Schmiedebergs Arch. Pharmacol.* **326(1)**, 36-44 (1984).
2. Giordano, J. and Rogers, L.V. Peripherally administered serotonin 5-HT₃ receptor antagonists reduce inflammatory pain in rats. *Eur. J. Pharmacol.* **170(1-2)**, 83-86 (1989).
3. Fukui, H., Yamamoto, M., and Sato, S. Vagal afferent fibers and peripheral 5-HT₃ receptors mediate cisplatin-induced emesis in dogs. *Jpn. J. Pharmacol.* **59(2)**, 221-226 (1992).

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, 06/17/2025

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