

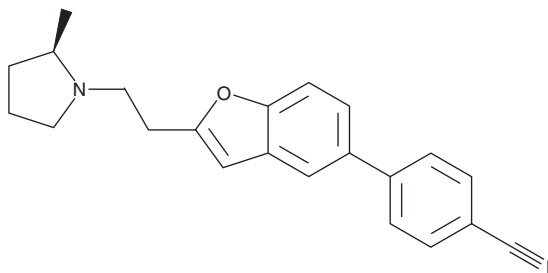
PRODUCT INFORMATION



ABT-239

Item No. 45575

CAS Registry No.: 460746-46-7
Formal Name: 4-[2-[2-[(2R)-2-methyl-1-pyrrolidinyl]ethyl]-5-benzofuran]-benzonitrile
MF: C₂₂H₂₂N₂O
FW: 330.4
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

ABT-239 is supplied as a solid. A stock solution may be made by dissolving the ABT-239 in the solvent of choice, which should be purged with an inert gas. ABT-239 is slightly soluble (0.1-1 mg/ml) in ethanol and DMSO.

Description

ABT-239 is a histamine H₃ receptor antagonist (K_s = 0.316 and 1.258 nM, respectively, for the human and rat receptors).¹ *In vivo*, ABT-239 (0.1-1 mg/kg) inhibits the dipsogenic response induced by (R)-(-)-α-methylhistamine (Item No. 25601) in mice. It improves social memory in adult and aged rats and attenuates methamphetamine-induced hyperactivity in mice. ABT-239 (3 mg/kg) decreases escape latency in the Morris water maze in a rat model of chronic stress.²

References

1. Fox, G.B., Esbenshade, T.A., Pan, J.B., *et al.* Pharmacological properties of ABT-239 [4-(2-[2-[(2R)-2-Methylpyrrolidinyl]ethyl]-benzofuran-5-yl)benzonitrile]: II. Neurophysiological characterization and broad preclinical efficacy in cognition and schizophrenia of a potent and selective histamine H₃ receptor antagonist. *J. Pharmacol. Exp. Ther.* **313**(1), 176-190 (2005).
2. Trofimiuk, E., Wielgat, P., and Car, H. Selective H₃ antagonist (ABT-239) differentially modifies cognitive function under the impact of restraint stress. *Front. Syst. Neurosci.* **14**, 614810 (2021).

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, 04/28/2026

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