

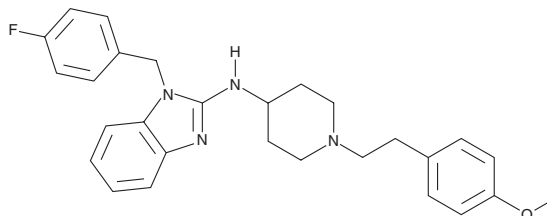
PRODUCT INFORMATION



Astemizole

Item No. 16967

CAS Registry No.: 68844-77-9
Formal Name: 1-[(4-fluorophenyl)methyl]-N-[1-[2-(4-methoxyphenyl)ethyl]-4-piperidinyl]-1H-benzimidazol-2-amine
Synonym: NSC 329963
MF: C₂₈H₃₁FN₄O
FW: 458.6
Purity: ≥98%
UV/Vis.: λ_{max}: 215, 250, 285 nm
Supplied as: A crystalline 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

Astemizole is supplied as a crystalline solid. A stock solution may be made by dissolving the astemizole in the solvent of choice, which should be purged with an inert gas. Astemizole is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of astemizole in these solvents is approximately 5, 20, and 10 mg/ml, respectively.

Astemizole is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, astemizole should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Astemizole has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Astemizole is a non-sedating antihistamine that antagonizes the histamine H1 receptor with a K_i value of 4.4 nM.¹ It is considerably less potent at muscarinic acetylcholine receptors (K_i = 2.4 μM).² Astemizole is also a potent blocker of ether-a-go-go-related gene (ERG) potassium channels (IC₅₀ = ~1 nM) that has been used to study long QT syndrome type 2.^{3,4} Furthermore, because it can target oncogenic ERG1 and ERG potassium channels, astemizole has been explored as a potential antineoplastic agent for decreasing proliferation of various cancer cells.⁵

References

1. Ahn, H.S. and Barnett, A. Selective displacement of [³H]mepyramine from peripheral vs. central nervous system receptors by loratadine, a non-sedating antihistamine. *Eur. J. Pharmacol.* **127(1-2)**, 153-155 (1986).
2. Kubo, N., Shirakawa, S., Kuno, T., *et al.* Antimuscarinic effects of antihistamines: Quantitative evaluation by receptor-binding assay. *Jpn. J. Pharmacol.* **43(3)**, 277-282 (1987).
3. Du, L.P., Tsai, K.C., Li, M.Y., *et al.* The pharmacophore hypotheses of I_{Kr} potassium channel blockers: Novel class III antiarrhythmic agents. *Bioorg. Med. Chem. Lett.* **14(18)**, 4771-4777 (2004).
4. Ficker, E., Obejero-Paz, C.A., Zhao, S., *et al.* The binding site for channel blockers that rescue misprocessed human long QT syndrome type 2 ether-a-gogo-related gene (HERG) mutations. *J. Biol. Chem.* **277(7)**, 4989-4998 (2002).
5. García-Quiroz, J., García-Becerra, R., Santos-Martínez, N., *et al.* In vivo dual targeting of the oncogenic Ether-à-go-go-1 potassium channel by calcitriol and astemizole results in enhanced antineoplastic effects in breast tumors. *BMC Cancer* **14**, 2-10 (2014).

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, 11/02/2022

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