

PRODUCT INFORMATION

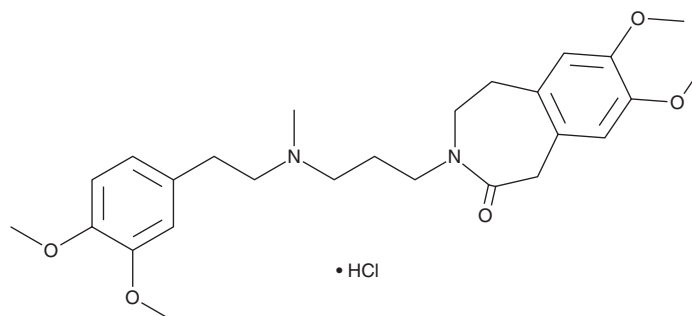


Zatebradine (hydrochloride)

Item No. 18608

CAS Registry No.: 91940-87-3
Formal Name: 3-[3-[[2-(3,4-dimethoxyphenyl)ethyl]methylamino]propyl]-1,3,4,5-tetrahydro-7,8-dimethoxy-2H-3-benzazepin-2-one, monohydrochloride

Synonym: UL-FS 49
MF: C₂₆H₃₆N₂O₅ • HCl
FW: 493.0
Purity: ≥95%
UV/Vis.: λ_{max}: 280 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

Zatebradine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the zatebradine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Zatebradine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of zatebradine (hydrochloride) in these solvents is approximately 1, 20, and 25 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of zatebradine (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of zatebradine (hydrochloride) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Zatebradine is a bradycardic compound that blocks hyperpolarization-activated inward current (I_h) through cyclic nucleotide-gated cation (HCN) channels in sinoatrial node cells (IC_{50} s = 1.83, 2.21, 1.9, and 1.88 μ M, respectively for HCN1-4 *in vitro*).^{1,2} Additionally, it can block voltage-gated outward K^+ (I_K) currents and related neuronal hyperpolarization-activated inward current ($I_{h,n}$) channels, but exhibits little or no activity for L-type Ca^{2+} (I_{Ca}) currents.^{3,4} When assessed through telemetric ECG recording in mice, zatebradine reduced heart rate from 600 to 200 beats per minute with an ED_{50} value of 1.8 mg/kg and induced increasing arrhythmia at concentrations >10 mg/kg.¹

References

1. Stieber, J., Wieland, K., Stöckl, G., *et al.* *Mol. Pharmacol.* **69**(4), 1328-1337 (2006).
2. Melchiorre, M., Del Lungo, M., Guandalini, L., *et al.* *J. Med. Chem.* **53**(18), 6773-6777 (2010).
3. Satoh, T.-O. and Yamada, M. *Pflugers. Arch.* **443**(4), 532-540 (2002).
4. BoSmith, R.E., Briggs, I., and Sturgess, N.C. *Br. J. Pharmacol.* **110**(1), 343-349 (1993).

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, 11/07/2022

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