

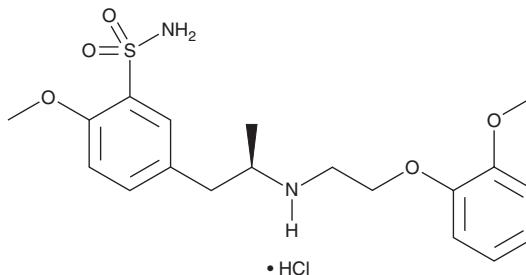
PRODUCT INFORMATION



Tamsulosin (hydrochloride)

Item No. 24020

CAS Registry No.: 106463-17-6
Formal Name: 5-[(2R)-2-[[2-(2-ethoxyphenoxy)ethyl]amino]propyl]-2-methoxy-benzenesulfonamide, monohydrochloride
Synonyms: LY253351, YM-12617-1
MF: C₂₀H₂₈N₂O₅S • HCl
FW: 445.0
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

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

Description

Tamsulosin is a potent antagonist of α_1 -adrenergic receptors (α_1 -ARs; K_i = 0.229 nM in a radioligand binding assay).¹ It is 3,800-fold selective for α_1 -ARs over α_2 -ARs (K_i = 871 nM). Tamsulosin antagonizes norepinephrine-induced contraction of isolated rabbit aorta (pA_2 = 10.11) but has no effect on contraction stimulated by histamine, serotonin (Item No. 14332), angiotensin II (Item No. 17150), or prostaglandin F_{2 α} (Item No. 16010). *In vivo*, tamsulosin reverses the pressor effect of phenylephrine (Item Nos. 17205 | 18619) in pithed rats. It reversibly reduces fertility in male rats when administered at a dose of 0.15 mg/kg.² Tamsulosin (1-100 μ g/kg) also reduces prostatic pressure in a dose-dependent manner with minimal hypotensive effects in anesthetized dogs.³ Formulations containing tamsulosin have been used for the treatment of benign prostatic hyperplasia.

References

1. Honda, K., Takenaka, T., Miyata-Osawa, A., *et al.* Studies on YM-12617: A selective and potent antagonist of postsynaptic α_1 -adrenoceptors. *Naunyn Schmiedeberg's Arch. Pharmacol.* **328**(3), 264-272 (1985).
2. Ratnasooriya, W.D. and Wadsworth, R.M. Tamsulosin, a selective alpha 1-adrenoceptor antagonist, inhibits fertility of male rats. *Andrologia* **26**(2), 107-110 (1994).
3. Sudoh, K., Tanaka, H., Inagaki, O., *et al.* Effect of tamsulosin, a novel alpha 1-adrenoceptor antagonist, on urethral pressure profile in anesthetized dogs. *J. Auton. Pharmacol.* **16**(3), 147-154 (1996).

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, 12/19/2022

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