

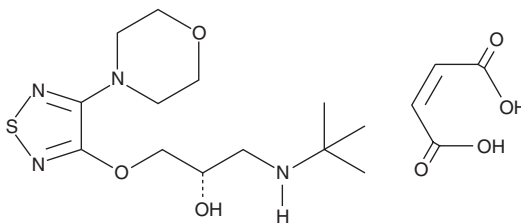
PRODUCT INFORMATION



Timolol (maleate)

Item No. 13974

CAS Registry No.: 26921-17-5
Formal Name: (2S)-1-[(1,1-dimethylethyl)amino]-3-[[4-(4-morpholinyl)-1,2,5-thiadiazol-3-yl]oxy]-2-propanol, 2Z-butenedioate
Synonyms: MK-950, (S)-Timolol (maleate), WP934
MF: C₁₃H₂₄N₄O₃S • C₄H₄O₄
FW: 432.5
Purity: ≥95%
UV/Vis.: λ_{max}: 209, 296 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

Timolol (maleate) is supplied as a crystalline solid. A stock solution may be made by dissolving the timolol (maleate) in the solvent of choice, which should be purged with an inert gas. Timolol (maleate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of timolol (maleate) in these solvents is approximately 0.2, 16, and 20 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 timolol (maleate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of timolol (maleate) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Timolol is a non-selective β -adrenergic receptor antagonist with log K_d values of -8.27, -9.86, and -6.8 for binding to human β_1 -, β_2 -, and β_3 -adrenoceptors, respectively.¹ It has been reported that only the (S) enantiomer contributes to the β -blocking effects of racemic timolol, but the weakly active (R) isomer maintains a beneficial effect on intraocular pressure without the undesirable side-effect of bronchial constriction caused by non-selective action of (S)-timolol on β_1 and β_2 receptors.^{2,3} Timolol has been used alone and in fixed combinations with either prostaglandin analogs or carbonic anhydrase inhibitors to reduce intraocular pressure in research models of ocular hypertension and glaucoma.^{4,5}

References

1. Baker, J.G. *Br. J. Pharmacol.* **144**(3), 317-322 (2005).
2. Tosi, G., Zironi, F., Caselli, E., et al. *Synthesis* **10**, 1625-1628 (2004).
3. Mehvar, R. and Brocks, D.R. *J. Pharm. Pharm. Sci.* **4**(2), 185-200 (2001).
4. Cheng, J.-W., Cheng, S.-W., Gao, L.-D., et al. *PLoS One* **7**(9), 45079 (2012).
5. Li, N., Chen, X., Zhou, Y., et al. *Clin. Exp. Ophthalmol.* **34**(8), 755-764 (2006).

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, 10/18/2022

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