

PRODUCT INFORMATION

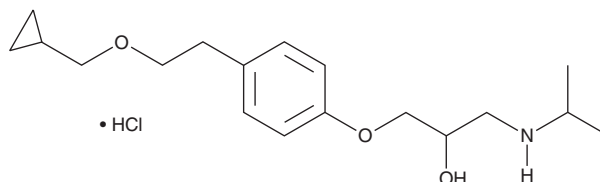


Betaxolol (hydrochloride)

Item No. 26285

CAS Registry No.: 63659-19-8
Formal Name: 1-[4-[2-(cyclopropylmethoxy)ethyl]phenoxy]-3-[(1-methylethyl)amino]-2-propanol, monohydrochloride

Synonym: SL 75212
MF: $C_{18}H_{29}NO_3 \cdot HCl$
FW: 343.9
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 224, 276 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 5 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Betaxolol (hydrochloride) (Item No. 26285) is an analytical reference standard categorized as a β_1 -adrenergic receptor antagonist.¹ Formulations containing β -adrenergic receptor antagonists have been used to enhance physical performance in athletes.^{2,3} This product is intended for use in analytical forensic applications. This product is also available as a general research tool (Item No. 18625).

References

1. Baker, J.G. The selectivity of β -adrenoceptor antagonists at the human β_1 , β_2 and β_3 adrenoceptors. *Br. J. Pharmacol.* **144**(3), 317-322 (2005).
2. Myung, S.-W. and Jo, C.-H. Gas chromatograph-mass spectrometric method for the determination of carvedilol and its metabolites in human urine. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **822**(1-2), 70-77 (2005).
3. Gupta, P. and Goyal, R.N. Amino functionalized graphene oxide and polymer nanocomposite based electrochemical platform for sensitive assay of anti-doping drug atenolol in biological fluids. *J. Electrochem. Soc.* **163**(13), B601-B608 (2016).

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, 07/19/2023

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