

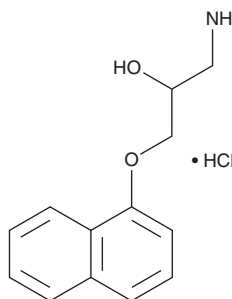
PRODUCT INFORMATION



Norpropranolol (hydrochloride)

Item No. 36011

CAS Registry No.: 62618-09-1
Formal Name: 1-amino-3-(1-naphthalenyloxy)-2-propanol, monohydrochloride
Synonyms: (±)-Desisopropylpropranolol, N-Desisopropylpropranolol
MF: C₁₃H₁₅NO₂ • HCl
FW: 253.7
Purity: ≥98%
UV/Vis.: λ_{max}: 214, 229 nm
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

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

Description

Norpropranolol is an active metabolite of the β-adrenergic receptor (β-AR) antagonist propranolol (Item No. 23349).^{1,2} It is formed from propranolol by the cytochrome P450 (CYP) isoform CYP1A2.² Norpropranolol binds to β-ARs (K_d = 42 nM in turkey erythrocytes) and inhibits (-)-isoproterenol-induced activation of adenylate cyclase with a K_i value of 175 nM.³

References

1. Bilezikian, J.P., Dornfeld, A.M., and Gammon, D.E. Structure-binding-activity analysis of beta-adrenergic amines—I. Binding to the beta receptor and activation of adenylate cyclase. *Biochem. Pharmacol.* **27(10)**, 1445-1454 (1978).
2. Masubuchi, Y., Hosokawa, S., Horie, T., *et al.* Cytochrome P450 isozymes involved in propranolol metabolism in human liver microsomes. The role of CYP2D6 as ring-hydroxylase and CYP1A2 as N-desisopropylase. *Drug Metab. Dispos.* **22(6)**, 909-915 (1994).
3. Bilezikian, J.P., Dornfeld, A.M., and Gammon, D.E. Structure-binding-activity analysis of beta-adrenergic amines—I. Binding to the beta receptor and activation of adenylate cyclase. *Biochem. Pharmacol.* **27(10)**, 1445-1454 (1978).

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, 12/15/2022

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