

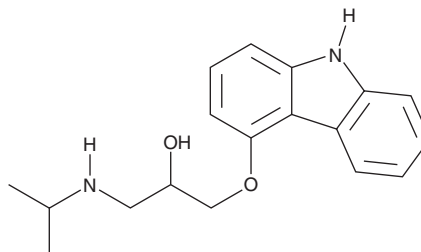
PRODUCT INFORMATION



Carazolol

Item No. 18628

CAS Registry No.: 57775-29-8
Formal Name: 1-(9H-carbazol-4-yloxy)-3-[(1-methylethyl)amino]-2-propanol
Synonyms: BM 51052, (±)-Carazolol
MF: C₁₈H₂₂N₂O₂
FW: 298.4
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 243, 286, 319, 332 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

Carazolol is supplied as a crystalline solid. A stock solution may be made by dissolving the carazolol in the solvent of choice, which should be purged with an inert gas. Carazolol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of carazolol in ethanol is approximately 5 mg/ml and approximately 15 mg/ml in DMSO and DMF.

Carazolol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, carazolol should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Carazolol has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Carazolol is a β -adrenergic receptor (β -AR) ligand (K_d s = 0.2, 0.03, and 4.47 nM for the human β_1 -, β_2 -, and β_3 -ARs, respectively).¹ It reduces epinephrine-induced increases in mean arterial blood pressure and heart rate in conscious pigs when administered at a dose of 10 μ g/kg.² Carazolol has been found in wastewater effluent.³ Formulations containing carazolol have been used to reduce stress in livestock during transport.

References

1. Baker, J.G. The selectivity of β -adrenoceptor agonists at human β_1 -, β_2 - and β_3 -adrenoceptors. *Br. J. Pharmacol.* **160**(5), 1048-1061 (2010).
2. Gregory, N.G. and Wilkins, L.J. The effect of carazolol on the cardiovascular responses to adrenaline in stress sensitive pigs. *Vet. Res. Commun.* **5**(3), 277-283 (1982).
3. Ternes, T.A. Occurrence of drugs in German sewage treatment plants and rivers. *Water Res.* **32**(11), 3245-3260 (1998).

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, 12/08/2022

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