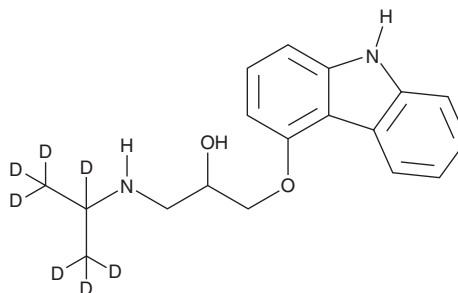


PRODUCT INFORMATION



Carazolol-d₇ Item No. 35560

CAS Registry No.: 1173021-02-7
Formal Name: 1-((9H-carbazol-4-yl)oxy)-3-((propan-2-yl-d₇)amino)propan-2-ol
Synonym: (±)-Carazolol-d₇
MF: C₁₈H₁₅D₇N₂O₂
FW: 305.4
Chemical Purity: ≥98% (Carazolol)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₇); ≤1% d₀
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

Carazolol-d₇ is intended for use as an internal standard for the quantification of carazolol (Item No. 18628) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Carazolol-d₇ is supplied as a solid. A stock solution may be made by dissolving the carazolol-d₇ in the solvent of choice, which should be purged with an inert gas. Carazolol-d₇ is slightly soluble in DMSO and methanol.

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, 12/12/2022

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