

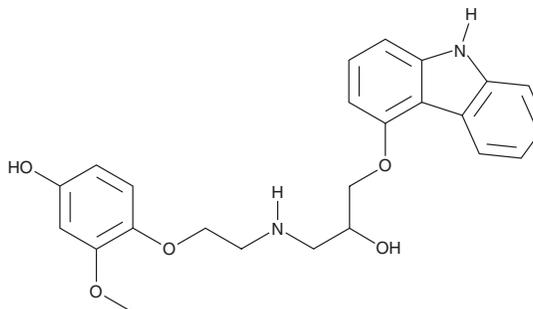
PRODUCT INFORMATION



4'-hydroxyphenyl Carvedilol

Item No. 27172

CAS Registry No.: 142227-49-4
Formal Name: 4-[2-[[3-(9H-carbazol-4-yloxy)-2-hydroxypropyl]amino]ethoxy]-3-methoxy-phenol
Synonym: BM 140686
MF: C₂₄H₂₆N₂O₅
FW: 422.5
Purity: ≥95%
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

4'-hydroxyphenyl Carvedilol is supplied as a solid. A stock solution may be made by dissolving the 4'-hydroxyphenyl carvedilol in the solvent of choice, which should be purged with an inert gas. 4'-hydroxyphenyl Carvedilol is slightly soluble in acetonitrile and methanol.

Description

4'-hydroxyphenyl Carvedilol is a metabolite of carvedilol (Item No. 15418) that is a more potent β -adrenergic receptor antagonist than carvedilol but has weaker vasodilatory activity.¹ It is formed through oxidation primarily by the cytochrome P450 (CYP) isoform CYP2D6.² 4'-hydroxyphenyl Carvedilol inhibits the formation of DPPH radicals in a cell-free assay.³

References

1. Patel, D.P., Sharma, P., Sanyal, M., *et al.* UPLC-MS/MS assay for the simultaneous quantification of carvedilol and its active metabolite 4'-hydroxyphenyl carvedilol in human plasma to support a bioequivalence study in healthy volunteers. *Biomed. Chromatogr.* **27(8)**, 974-986 (2013).
2. Oldham, H.G. and Clarke, S.E. *In vitro* identification of the human cytochrome P450 enzymes involved in the metabolism of R(+)- and S(-)-carvedilol. *Drug Metab. Dispos.* **25(8)**, 970-977 (1997).
3. Malig, T.C., Ashkin, M.R., Burman, A.L., *et al.* Comparison of free-radical inhibiting antioxidant properties of carvedilol and its phenolic metabolites. *Medchemcomm.* **8(3)**, 606-615 (2017).

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, 11/29/2022

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