

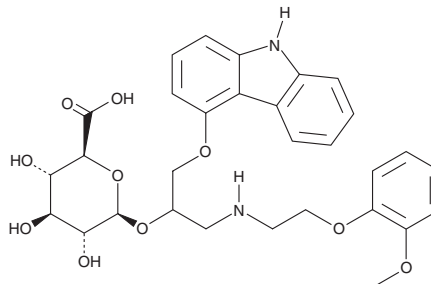
PRODUCT INFORMATION



(R,S)-Carvedilol Glucuronide

Item No. 27173

CAS Registry No.: 114869-83-9
Formal Name: 2-(9H-carbazol-4-yloxy)-1-[[[2-(2-methoxyphenoxy)ethyl]amino]methyl]ethyl β-D-glucopyranosiduronic acid
MF: C₃₀H₃₄N₂O₁₀
FW: 582.6
Purity: ≥95% (mixture of diastereomers)
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

(R,S)-Carvedilol glucuronide is supplied as a solid. A stock solution may be made by dissolving the (R,S)-carvedilol glucuronide in the solvent of choice, which should be purged with an inert gas. (R,S)-Carvedilol glucuronide is slightly soluble in methanol.

Description

(R,S)-Carvedilol glucuronide is a racemic mixture of the carvedilol (Item No. 15418) metabolites (R)-carvedilol glucuronide and (S)-carvedilol glucuronide.¹ (R)-Carvedilol glucuronide is formed *via* glucuronidation of (R)-carvedilol by the UDP-glucuronosyltransferase (UGT) isoforms UGT1A1 and UGT2B4. (S)-Carvedilol glucuronide is formed *via* glucuronidation of (S)-carvedilol by UGT2B4 and UGT2B7.

Reference

1. Hanioka, N., Tanaka, S., Moriguchi, Y., *et al.* Stereoselective glucuronidation of carvedilol in human liver and intestinal microsomes. *Pharmacology* **90**(3-4), 117-124 (2012).

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, 12/19/2022

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