

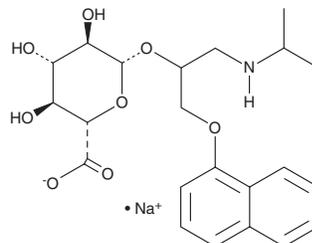
PRODUCT INFORMATION



Propranolol β -D-Glucuronide (sodium salt)

Item No. 22382

Formal Name: β -D-glucopyranosiduronic acid, 1-[[[(1-methylethyl)amino]methyl]-2-(1-naphthalenyloxy)ethyl], monosodium salt
MF: $C_{22}H_{28}NO_8 \cdot Na$
FW: 457.5
Purity: $\geq 95\%$ (mixture of diastereomers)
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Propranolol β -D-glucuronide (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the propranolol β -D-glucuronide (sodium salt) in the solvent of choice. Propranolol β -D-glucuronide (sodium salt) is soluble in the organic solvent methanol, which should be purged with an inert gas.

Description

Propranolol β -D-glucuronide is a metabolite of the β -adrenergic receptor antagonist propranolol (Item No. 23349).¹ It is formed from propranolol by the UDP-glucuronosyltransferase (UGT) isoforms UGT1A7, UGT1A9, UGT1A10, and UGT2A1.

Reference

1. Yang, F., Liu, S., Wolber, G., *et al.* Complete reaction phenotyping of propranolol and 4-hydroxypropranolol with the 19 enzymes of the human UGT1 and UGT2 families. *Int. J. Mol. Sci.* **23(13)**, 7476 (2022).

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, 03/09/2023

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