

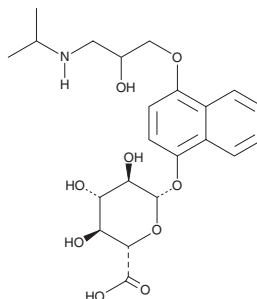
# PRODUCT INFORMATION



## 4-hydroxy Propranolol $\beta$ -D-Glucuronide

Item No. 22383

**CAS Registry No.:** 94731-13-2  
**Formal Name:** 4-[2-hydroxy-3-[(1-methylethyl)amino]propoxy]-1-naphthalenyl  $\beta$ -D-glucopyranosiduronic acid  
**MF:** C<sub>22</sub>H<sub>29</sub>NO<sub>9</sub>  
**FW:** 451.5  
**Purity:**  $\geq 95\%$  (mixture of diastereomers)  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

4-hydroxy Propranolol  $\beta$ -D-glucuronide is supplied as a solid. A stock solution may be made by dissolving the 4-hydroxy propranolol  $\beta$ -D-glucuronide in water. We do not recommend storing the aqueous solution for more than one day.

### Description

4-hydroxy Propranolol  $\beta$ -D-glucuronide is a metabolite of the  $\beta$ -adrenergic receptor antagonist 4-hydroxy propranolol (Item No. 18630).<sup>1</sup> It is formed from propranolol (Item No. 23349) via a 4-hydroxy propranolol intermediate by the UDP-glucuronosyltransferase (UGT) isoforms UGT1A9 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, 03/31/2026

#### 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