

# PRODUCT INFORMATION



## Nebivolol O-β-D-Glucuronide

Item No. 26657

**Formal Name:** (2S,3S,4S,5R,6R)-6-(1-(6-fluorochroman-2-yl)-2-((2-(6-fluorochroman-2-yl)-2-hydroxyethyl)amino)ethoxy)-3,4,5-trihydroxytetrahydro-2H-pyran-2-carboxylic acid

**MF:**  $C_{28}H_{33}F_2NO_{10}$

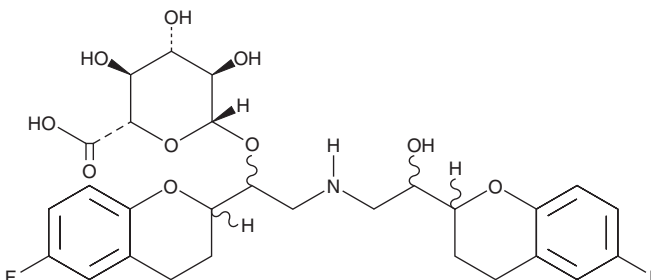
**FW:** 581.6

**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

Nebivolol O-β-D-glucuronide is supplied as a solid. A stock solution may be made by dissolving the nebivolol O-β-D-glucuronide in the solvent of choice, which should be purged with an inert gas. Nebivolol O-β-D-glucuronide is slightly soluble in methanol and chloroform (warmed).

### Description

Nebivolol O-β-D-glucuronide is an active metabolite of the  $\beta_1$ -adrenergic receptor antagonist nebivolol (Item No. 23660).<sup>1,2</sup> This product is a mixture of D-nebivolol-11-O-β-D-glucuronide, L-nebivolol-11-O-β-D-glucuronide, D-nebivolol-11'-O-β-D-glucuronide, and L-nebivolol-11'-O-β-D-glucuronide.

### References

1. Vieira, C.P., Neves, D.V., Cesarino, E.J., *et al.* An indirect stereoselective analysis of nebivolol glucuronides in plasma by LC-MS/MS: Application to clinical pharmacokinetics. *J. Pharm. Biomed. Anal.* **144**, 25-30 (2017).
2. O'Donnell, J.P., Owens, W., Duncan, J., *et al.* Glucuronidated Nebivolol Metabolites. *Forest Laboratories Holdings Ltd.* **US 2007/0014734 A1** (2007).

#### 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](http://WWW.CAYMANCHEM.COM)