

PRODUCT INFORMATION

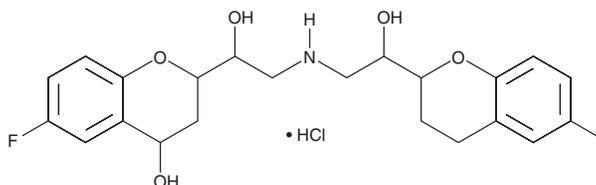


4-hydroxy Nebivolol (hydrochloride)

Item No. 26658

CAS Registry No.: 2717115-85-8
Formal Name: 6-fluoro-α-[[[2-(6-fluoro-3,4-dihydro-2H-1-benzopyran-2-yl)-2-hydroxyethyl]amino methyl]-3,4-dihydro-4-hydroxy-2H-1-benzopyran-2-methanol, monohydrochloride

MF: C₂₂H₂₅F₂NO₅ • HCl
FW: 457.9
Purity: ≥98% (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

4-hydroxy Nebivolol (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the 4-hydroxy nebivolol (hydrochloride) in the solvent of choice, which should be purged with an inert gas. 4-hydroxy Nebivolol (hydrochloride) is soluble in the organic solvent DMSO.

Description

4-hydroxy Nebivolol is a major metabolite of nebivolol (Item No. 23660).¹ It is formed by the hydroxylation of nebivolol by the cytochrome P450 (CYP) isoform CYP2D6.

Reference

1. Hu, X., Lan, T., Dai, D., *et al.* Evaluation of 24 CYP2D6 variants on the metabolism of nebivolol in vitro. *Drug Metab. Dispos.* **44(11)**, 1828-1831 (2016).

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, 01/27/2023

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