

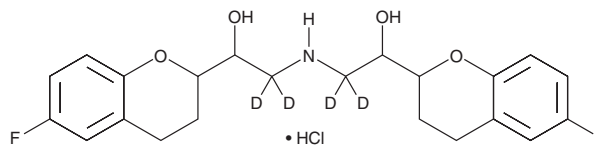
# PRODUCT INFORMATION



## (±)-Nebivolol-d<sub>4</sub> (hydrochloride)

Item No. 26659

**CAS Registry No.:** 2701283-32-9  
**Formal Name:** 2,2'-azanediyibis(1-(6-fluorochroman-2-yl)ethan-2,2-d<sub>2</sub>-1-ol), monohydrochloride  
**MF:** C<sub>22</sub>H<sub>21</sub>D<sub>4</sub>F<sub>2</sub>NO<sub>4</sub> • HCl  
**FW:** 445.9  
**Chemical Purity:** ≥98% (mixture of diastereomers; Nebivolol)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>4</sub>); ≤1% d<sub>0</sub>  
**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-d<sub>4</sub> (hydrochloride) is intended for use as an internal standard for the quantification of nebivolol (Item No. 23660) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

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

### Description

(±)-Nebivolol-d<sub>4</sub> is intended for use as an internal standard for the quantification of nebivolol (Item No. 23660) by GC- or LC-MS. Nebivolol is an antagonist of the β<sub>1</sub>-adrenergic receptor (β<sub>1</sub>-AR; IC<sub>50</sub> = 7.41 nM).<sup>1</sup> It is selective for β<sub>1</sub>- over β<sub>2</sub>-ARs (IC<sub>50</sub> = 251 nM), as well as the serotonin (5-HT) receptor subtypes 5-HT<sub>1A</sub> and 5-HT<sub>2</sub> and the α<sub>1</sub>- and α<sub>2</sub>-adrenergic, histamine H<sub>1</sub>, and dopamine D<sub>2</sub> receptors (IC<sub>50</sub>s = 27.5, 2,239, 3,162, >10,000, 5,623, and 10,000 nM, respectively). Nebivolol induces vasodilation in isolated mouse renal arteries (EC<sub>50</sub> = 11.36 μM) and decreases contraction of isolated human left ventricular trabeculae induced by isoproterenol (Item No. 15592; IC<sub>50</sub> = 7 μM).<sup>2,3</sup> Nebivolol inhibits proliferation of primary human coronary artery smooth muscle cells (HCASMCs) in the presence and absence of growth factors (IC<sub>50</sub>s = 6.1, 6.8, 6.4, and 7.7 μM for HCASMCs grown in media containing no growth factor, PDGF-BB, basic FGF, and TGF-β1, respectively).<sup>4</sup> It is also an inhibitor of the severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) main protease (M<sup>Pro</sup>), also known as 3C-like protease (3CL<sup>Pro</sup>; IC<sub>50</sub> = 60.2 μg/ml), and inhibits SARS-CoV-2 pathogenicity *in vitro* (IC<sub>50</sub> = 0.03 μg/ml).<sup>5</sup> Formulations containing nebivolol have been used in the treatment of hypertension.

### References

1. Pauwels, P.J., Gommeren, W., Van Lommen, G., *et al.* *Mol. Pharmacol.* **34**(6), 843-851 (1988).
2. Georgescu, A., Pluteanu, F., Flonta, M.L., *et al.* *Pharmacology* **81**(2), 110-117 (2008).
3. Brixius, K., Bundkirchen, A., Böck, B., *et al.* *Br. J. Pharmacol.* **133**(8), 1330-1338 (2001).
4. Brehm, B.R., Wolf, S.C., Bertsch, D., *et al.* *Cardiovasc. Res.* **49**(2), 430-439 (2001).
5. Hamed, M.I.A., Darwish, K.M., Soltane, R., *et al.* *RSC Adv.* **11**(56), 35536-35558 (2021).

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