

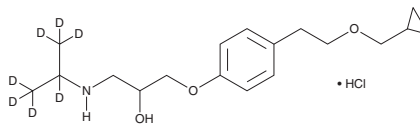
# PRODUCT INFORMATION



## Betaxolol-d<sub>7</sub> (hydrochloride)

Item No. 39815

**CAS Registry No.:** 1219802-92-2  
**Formal Name:** 1-(4-(2-(cyclopropylmethoxy)ethyl)phenoxy)-3-((propan-2-yl-d<sub>7</sub>)amino)propan-2-ol, monohydrochloride  
**MF:** C<sub>18</sub>H<sub>22</sub>D<sub>7</sub>NO<sub>3</sub> • HCl  
**FW:** 350.9  
**Chemical Purity:** ≥98% (Betaxolol)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>7</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

Betaxolol-d<sub>7</sub> (hydrochloride) is intended for use as an internal standard for the quantification of betaxolol (Item No. 18625) 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).

Betaxolol-d<sub>7</sub> (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the betaxolol-d<sub>7</sub> (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Betaxolol-d<sub>7</sub> (hydrochloride) is slightly soluble in chloroform and methanol.

### Description

Betaxolol is an antagonist of the β<sub>1</sub>-adrenergic receptor (β<sub>1</sub>-AR; K<sub>i</sub> = 23 nM).<sup>1</sup> It is selective for the β<sub>1</sub>-AR over the β<sub>2</sub>-AR (K<sub>i</sub> = 790 nM). Betaxolol is cytotoxic to 10.014 pRSV-T human corneal epithelial cells.<sup>2</sup> It reduces intraocular pressure in cynomolgus monkeys with ocular hypertension induced by laser trabeculoplasty when administered at a dose of 150 μg/animal.<sup>3</sup> Betaxolol also inhibits sodium influx induced by the neurotoxin veratridine in rat cortical synaptosomes (IC<sub>50</sub> = 28.3 μM).<sup>4</sup>

### References

1. Wellstein, A., Palm, D., and Belz, G.G. Affinity and selectivity of β-adrenoceptor antagonists *in vitro*. *J. Cardiovasc. Pharmacol.* **8(Suppl. 11)**, S36-S40 (1986).
2. Pozarowska, D., Pozarowski, P., and Darzynkiewicz, Z. Cytometric assessment of cytostatic and cytotoxic effects of topical glaucoma medications on human epithelial corneal line cells. *Cytometry B Clin. Cytom.* **78(2)**, 130-137 (2010).
3. Sharif, N.A., Xu, S.X., Crider, J.Y., *et al.* Levobetaxolol (Betaxon™) and other β-adrenergic antagonists: Preclinical pharmacology, IOP-lowering activity and sites of action in human eyes. *J. Ocul. Pharmacol. Ther.* **17(4)**, 305-317 (2001).
4. Chidlow, G., Melena, J., and Osborne, N.N. Betaxolol, a β<sub>1</sub>-adrenoceptor antagonist, reduces Na<sup>+</sup> influx into cortical synaptosomes by direct interaction with Na<sup>+</sup> channels: Comparison with other β-adrenoceptor antagonists. *Br. J. Pharmacol.* **130(4)**, 759-766 (2000).

#### 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, 10/11/2023

#### 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