

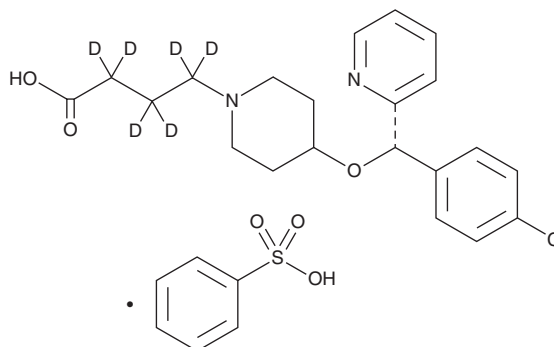
# PRODUCT INFORMATION



## Bepotastine-d<sub>6</sub> (besylate)

Item No. 40233

**CAS Registry No.:** 2469626-71-7  
**Formal Name:** 4-[(S)-(4-chlorophenyl)-2-pyridinylmethoxy]-1-piperidinebutanoic-d<sub>6</sub> acid, monobenzenesulfonate  
**Synonym:** Bepotastine besilate-d<sub>6</sub>  
**MF:** C<sub>21</sub>H<sub>19</sub>ClD<sub>6</sub>N<sub>2</sub>O<sub>3</sub> • C<sub>6</sub>H<sub>6</sub>O<sub>3</sub>S  
**FW:** 553.1  
**Chemical Purity:** ≥95% (Bepotastine)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>6</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

Bepotastine-d<sub>6</sub> (besylate) is intended for use as an internal standard for the quantification of bepotastine (Item No. 23721) 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).

Bepotastine-d<sub>6</sub> (besylate) is supplied as a solid. A stock solution may be made by dissolving the bepotastine-d<sub>6</sub> (besylate) in the solvent of choice, which should be purged with an inert gas. Bepotastine-d<sub>6</sub> (besylate) is slightly soluble in DMSO (warmed) and methanol.

### Description

Bepotastine is an antagonist of the histamine H<sub>1</sub> receptor that is selective over H<sub>3</sub>, α<sub>1</sub>-, α<sub>2</sub>-, and β-adrenergic, dopamine D<sub>2long</sub>, serotonin 5-HT<sub>2</sub>, muscarinic acetylcholine, and benzodiazepine receptors.<sup>1</sup> It reduces dye leakage from the nasal passages of rats acutely sensitized to an antigen (ED<sub>50</sub> = 0.03 mg/kg) and inhibits histamine-induced bronchoconstriction in the anesthetized dog (ED<sub>50</sub> = 3.2 μg/kg).<sup>2,3</sup> Bepotastine prevents conjunctival vascular hyperpermeability in a guinea pig model of conjunctivitis in a dose-dependent manner.<sup>4</sup> Formulations containing bepotastine have been used in the treatment of itching associated with allergic conjunctivitis.

### References

1. Kato, M., Nishida, A., Aga, Y., *et al.* Pharmacokinetic and pharmacodynamic evaluation of central effect of the novel antiallergic agent betotastine besilate. *Arzneimittelforschung*. **47(10)**, 1116-1124 (1997).
2. Murata, T., Matsumoto, Y., Suzuki, T., *et al.* Effect of betotastine besilate (TAU-284), a novel anti-allergic agent, on experimental allergic rhinitis. *Arerugi* **46(7)**, 576-584 (1997).
3. Matsubara, S., Ono, C., Yamazaki, N., *et al.* Inhibitory effects of bepotastine, a novel anti-allergic drug on histamine-induced bronchoconstriction in anesthetized dogs. *Yakuri to Chiryō* **25(4)**, 895-900 (1997).
4. Kida, T., Fuji, A., Sakai, O., *et al.* Bepotastine besilate, a highly selective histamine H<sub>1</sub> receptor antagonist, suppresses vascular hyperpermeability and eosinophil recruitment in in vitro and in vivo experimental allergic conjunctivitis models. *Exp. Eye Res.* **91(1)**, 85-91 (2010).

#### 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, 11/28/2023

#### 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