

PRODUCT INFORMATION



Trospium-d₈ (chloride)

Item No. 26686

Formal Name: (1R,3s,5S)-3-(2-hydroxy-2,2-diphenylacetoxy)spiro[bicyclo[3.2.1]octane-8,1'-pyrrolidin]-8-ium-2',2',3',3',4',4',5',5'-d₈, monochloride

Synonym: Relaspium-d₈

MF: C₂₅H₂₂D₈NO₃ • Cl

FW: 436.0

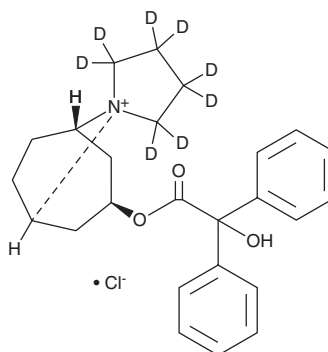
Chemical Purity: ≥98% (Trospium)

Deuterium Incorporation: ≥99% deuterated forms (d₁-d₈); ≤1% d₀

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

Trospium-d₈ (chloride) is intended for use as an internal standard for the quantification of trospium (Item No. 21279) 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).

Trospium-d₈ (chloride) is supplied as a solid. A stock solution may be made by dissolving the trospium-d₈ (chloride) in the solvent of choice. Trospium-d₈ (chloride) is soluble in the organic solvent DMSO, which should be purged with an inert gas.

Description

Trospium is an antagonist of muscarinic (M) receptors (K_is = 0.50-2.3 nM for human M₁₋₅ recombinant receptors, respectively).¹ Formulations containing antimuscarinics, including trospium, are used to manage overactive bladder by relieving urgency, frequency, and incontinence.¹⁻³

References

1. Hegde, S.S. Muscarinic receptors in the bladder: From basic research to therapeutics. *Br. J. Pharmacol.* **147**(2), S80-S87 (2006).
2. Maggiore, U.L.R., Salvatore, S., Alessandri, F., *et al.* Pharmacokinetics and toxicity of antimuscarinic drugs for overactive bladder treatment in females. *Expert Opin. Drug Metab. Toxicol.* **8**(11), 1387-1408 (2012).
3. Michel, M.C. and de la Rosette, J.J.M.C.H. Role of muscarinic receptor antagonists in urgency and nocturia. *BJU Int.* **96**(Suppl 1), 37-42 (2005).

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, 12/20/2022

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