

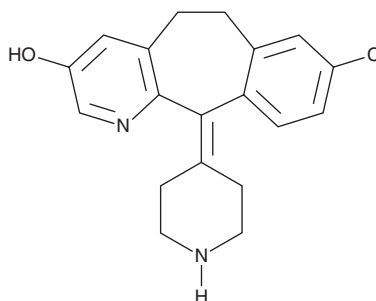
PRODUCT INFORMATION



3-hydroxy Desloratadine

Item No. 21847

CAS Registry No.: 119410-08-1
Formal Name: 8-chloro-6,11-dihydro-11-(4-piperidinylidene)-5H-benzo[5,6]cyclohepta[1,2-b]pyridin-3-ol
Synonym: SCH 45581
MF: C₁₉H₁₉ClN₂O
FW: 326.8
Purity: ≥98%
UV/Vis.: λ_{max}: 245, 291 nm
Supplied as: A crystalline 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

3-hydroxy Desloratadine is supplied as a crystalline solid. A stock solution may be made by dissolving the 3-hydroxy desloratadine in the solvent of choice, which should be purged with an inert gas. 3-hydroxy Desloratadine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 3-hydroxy desloratadine in ethanol is approximately 5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

3-hydroxy Desloratadine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 3-hydroxy desloratadine should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 3-hydroxy Desloratadine has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

3-hydroxy Desloratadine is a major metabolite of desloratadine (Item No. 16931), a tricyclic antagonist of the histamine H₁ receptor.¹

Reference

1. Afrime, M., Gupta, S., Banfield, C., *et al.* A pharmacokinetic profile of desloratadine in healthy adults, including elderly. *Clin. Pharmacokinet.* **41(Suppl 1)**, 13-19 (2002).

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, 01/16/2023

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