

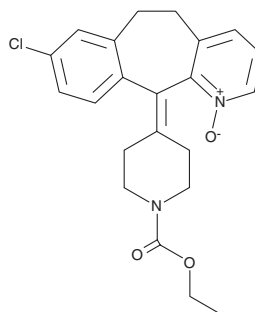
PRODUCT INFORMATION



Loratadine N-oxide

Item No. 35244

CAS Registry No.: 165739-62-8
Formal Name: 4-(8-chloro-5,6-dihydro-1-oxido-11H-benzo[5,6]cyclohepta[1,2-b]pyridin-11-ylidene)-1-piperidinecarboxylic acid, ethyl ester
Synonym: SCH 38554
MF: C₂₂H₂₃ClN₂O₃
FW: 398.9
Purity: ≥95%
UV/Vis.: λ_{max}: 259 nm
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

Loratadine N-oxide is supplied as a solid. A stock solution may be made by dissolving the loratadine N-oxide in the solvent of choice, which should be purged with an inert gas. Loratadine N-oxide is slightly soluble in chloroform and methanol.

Description

Loratadine N-oxide is a metabolite of the histamine H₁ receptor inverse agonist loratadine (Item No. 15625).¹

Reference

1. Habicht, S.C., Duan, P., Vinueza, N.R., *et al.* Liquid chromatography/tandem mass spectrometry utilizing ion-molecule reactions and collision-activated dissociation for the identification of N-oxide drug metabolites. *J. Pharm. Biomed. Anal.* **51(4)**, 805-811 (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, 11/19/2024

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