

PRODUCT INFORMATION



Rupatadine (fumarate)

Item No. 29478

CAS Registry No.: 182349-12-8
Formal Name: 8-chloro-6,11-dihydro-11-[1-[(5-methyl-3-pyridinyl)methyl]-4-piperidinylidene]-5H-benzo[5,6]cyclohepta[1,2-b]pyridine, (2E)-2-butenedioate

MF: C₂₆H₂₆ClN₃ • C₄H₄O₄

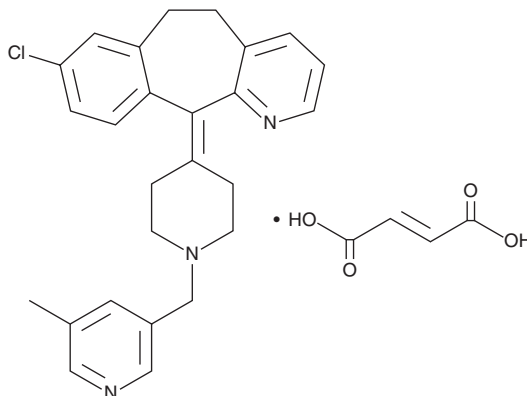
FW: 532.0

Purity: ≥98%

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

Rupatadine (fumarate) is supplied as a crystalline solid. A stock solution may be made by dissolving the rupatadine (fumarate) in the solvent of choice, which should be purged with an inert gas. Rupatadine (fumarate) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of rupatadine (fumarate) in these solvents is approximately 1 mg/ml.

Rupatadine (fumarate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, rupatadine (fumarate) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Rupatadine (fumarate) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Rupatadine is an antagonist of histamine H₁ and platelet-activating factor (PAF) receptors.¹ It binds to rabbit platelet membranes and guinea pig cerebellum membranes with apparent K_i values of 0.55 and 0.1 μM, respectively, in radioligand binding assays. It inhibits contraction of isolated guinea pig ileum induced by histamine (pA₂ = 9.29) and platelet aggregation induced by PAF in washed rabbit platelets (pA₂ = 6.68). Rupatadine prevents histamine- or PAF-induced hypotension in rats (ID₅₀s = 1.4 and 0.44 mg/kg, respectively) and reverses histamine- or PAF-induced bronchoconstriction in guinea pigs (ID₅₀s = 113 and 9.6 mg/kg, respectively). It does not reduce locomotor activity or potentiate sleep induced by barbiturates in mice, indicating a lack of sedative effects.

Reference

1. Merlos, M., Giral, M., Balsa, D., *et al.* Rupatadine, a new potent, orally active dual antagonist of histamine and platelet-activating factor (PAF). *J. Pharmacol. Exp. Ther.* **280**(1), 114-121 (1997).

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, 12/09/2022

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