

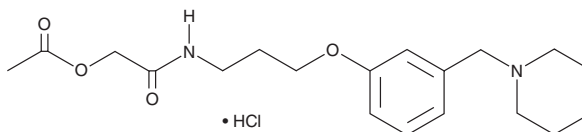
PRODUCT INFORMATION



Roxatidine Acetate (hydrochloride)

Item No. 21248

CAS Registry No.: 93793-83-0
Formal Name: 2-(acetyloxy)-N-[3-[3-(1-piperidinylmethyl)phenoxy]propyl]-acetamide, monohydrochloride
Synonyms: HOE 760, TZU-0460
MF: C₁₉H₂₈N₂O₄ • HCl
FW: 384.9
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 277 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

Roxatidine acetate (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the roxatidine acetate (hydrochloride) in the solvent of choice. Roxatidine acetate (hydrochloride) is soluble in organic solvents such as ethanol and DMSO. It is also soluble in water. The solubility of roxatidine acetate (hydrochloride) in ethanol, DMSO, and water is approximately 12, 78, and 77 mg/ml, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

Roxatidine acetate is a competitive histamine H₂ receptor antagonist that upon oral absorption is rapidly converted to its active metabolite roxatidine.¹ By inhibiting the binding of histamine to H₂ receptors, roxatidine reduces both intracellular cAMP concentrations and gastric acid secretion by parietal cells.¹

Reference

1. Murdoch, D., and McTavish, D. Roxatidine acetate. A review of its pharmacodynamic and pharmacokinetic properties, and its therapeutic potential in peptic ulcer disease and related disorders. *Drugs* **42(2)**, 240-260 (1991).

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, 10/24/2022

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