

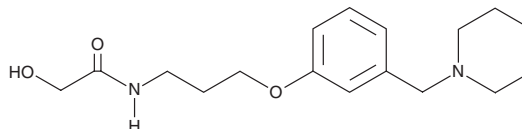
PRODUCT INFORMATION



Roxatidine

Item No. 36111

CAS Registry No.: 78273-80-0
Formal Name: 2-hydroxy-N-[3-[3-(1-piperidinylmethyl)phenoxy]propyl]-acetamide
MF: C₁₇H₂₆N₂O₃
FW: 306.4
Purity: ≥95%
Supplied as: A neat oil
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 is supplied as a neat oil. A stock solution may be made by dissolving the roxatidine in the solvent of choice, which should be purged with an inert gas. Roxatidine is soluble in the organic solvent DMSO at a concentration of up to 25 mg/ml.

Description

Roxatidine is a histamine H₂ receptor antagonist and major active metabolite of roxatidine acetate (Item No. 21248).¹ Roxatidine reduces histamine-induced adenylate cyclase production in guinea pig parietal cells (IC₅₀ = 0.8 μM). It inhibits histamine-induced hydrogen ion accumulation in the same cells (pA₂ = 7.03). Roxatidine (200 mg/kg) reduces small intestinal lesion area in a rat model of gastric mucosal injury induced by indomethacin (Item No. 70270).²

References

1. Sewing, K.-F., Beil, W., and Hannemann, H. Comparative pharmacology of histamine H₂-receptor antagonists. *Drugs* **35(Suppl. 3)**, 25-29 (1988).
2. Umegaki, E., Yoda, Y., Tokioka, S., et al. Protective effect of roxatidine against indomethacin induced small intestinal mucosal injury in rats. *J. Gastroenterol. Hepatol.* **25(Suppl. 1)**, S35-S40 (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, 11/15/2022

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