

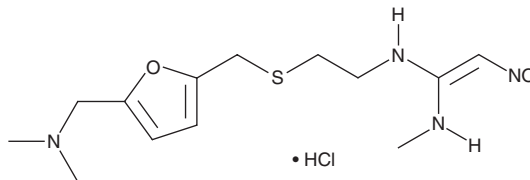
PRODUCT INFORMATION



Ranitidine (hydrochloride)

Item No. 16939

CAS Registry No.: 66357-59-3
Formal Name: N'-[2-[[[5-[(dimethylamino)methyl]-2-furanyl]methyl]thio]ethyl]-N-methyl-2-nitro-1,1-ethenediamine, monohydrochloride
MF: C₁₃H₂₂N₄O₃S • HCl
FW: 350.9
Purity: ≥98%
UV/Vis.: λ_{max}: 230, 325 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

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

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of ranitidine (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of ranitidine (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Ranitidine is a histamine H₂ receptor antagonist.¹ It reverses histamine-induced relaxation of isolated rat uterine horn (pA₂ = 6.9) as well as histamine-induced increases in contraction frequency in isolated guinea pig right atrium (pA₂ = 7.2). Ranitidine (0.03-3 mg/kg, i.v.) inhibits histamine- and pentagastrin-induced gastric acid secretion in a dose-dependent manner in anesthetized rats. Formulations containing ranitidine have been used in the treatment and prevention of heartburn and gastroesophageal reflux disease (GERD).

Reference

1. Daly, M.J., Humphray, J.M., and Stables, R. Some *in vitro* and *in vivo* actions of the new histamine H₂-receptor antagonist, ranitidine. Br. J. Pharmacol. 72(1), 49-54 (1981).

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, 11/16/2022

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