

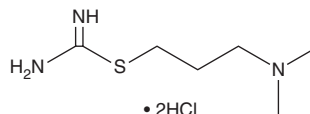
PRODUCT INFORMATION



Dimaprit (hydrochloride)

Item No. 29418

CAS Registry No.: 23256-33-9
Formal Name: carbamimidithioic acid,
3-(dimethylamino)propyl ester,
dihydrochloride
MF: C₆H₁₅N₃S • 2HCl
FW: 234.2
Purity: ≥95%
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

Dimaprit (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the dimaprit (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Dimaprit (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of dimaprit (hydrochloride) in these solvents is approximately 30 mg/ml.

Dimaprit (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, dimaprit (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Dimaprit (hydrochloride) 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

Dimaprit is a histamine H₂ receptor agonist with a K_i value of 44 μM in guinea pig right atrium.^{1,2} It is selective for histamine H₂ over H₁ and H₃ receptors with relative potencies of 71, <0.0001, and <0.008, respectively, compared to histamine. Dimaprit (6 μM) inhibits histamine release from isolated peritoneal mast cells in a rat model of anaphylaxis induced by *A. suum* extract, which is not affected by pretreatment with mepyramine (Item No. 20978), an H₁ receptor inverse agonist, or cimetidine (Item No. 18743), an H₂ receptor antagonist.³ It stimulates gastric acid secretion *in vivo* in anaesthetized rats and cats with relative potencies of 19.5 and approximately 400 to 500, respectively, compared to histamine.⁴ Dimaprit inhibits 2-(2-pyridyl)ethylamine-induced bronchoconstriction in guinea pigs in a dose-dependent manner.⁵

References

1. Hill, S.J. Distribution, properties, and functional characteristics of three classes of histamine receptor. *Pharmacol. Rev.* **42(1)**, 45-83 (1990).
2. Rising, T.J. and Steward, A. The determination of receptor constants for histamine H₂-agonists in the guinea-pig isolated right atrium using an irreversible H₂-antagonist. *Br. J. Pharmacol.* **87(1)**, 211-216 (1986).
3. Kohno, S., Ogawa, K., Nabe, T., *et al.* Dimaprit, a histamine H₂-agonist, inhibits anaphylactic histamine release from mast cells and the decreased release is restored by thioperamide (H₃-antagonist), but not by cimetidine (H₂-antagonist). *Jpn. J. Pharmacol.* **62(1)**, 75-79 (1993).
4. Parsons, M.E., Owen, D.A.A., Ganellin, C.R., *et al.* Dimaprit-[S-[3-(N,N-dimethylamino)propyl]isothioure] – A highly specific histamine H₂-receptor agonist. Part 1. Pharmacology. *Agents Actions* **7(1)**, 31-37 (1977).
5. Drazen, J.M., Schneider, M.W., and Venugopalan, C.S. Bronchodilator activity of dimaprit in the guinea pig *in vitro* and *in vivo*. *Eur. J. Pharmacol.* **55(3)**, 233-239 (1979).

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, 12/07/2022

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