

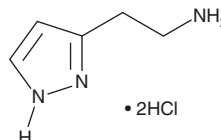
PRODUCT INFORMATION



Betazole (hydrochloride)

Item No. 41712

CAS Registry No.: 138-92-1
Formal Name: 1H-pyrazole-3-ethanamine, dihydrochloride
Synonyms: Ametazole, Gastramine
MF: $C_5H_9N_3 \cdot 2HCl$
FW: 184.1
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Betazole (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the betazole (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Betazole (hydrochloride) is soluble (≥ 10 mg/ml) in DMSO.

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 betazole (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. Betazole (hydrochloride) is soluble (≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Betazole is a histamine H₂ receptor agonist.¹ It induces acid secretion in isolated human gastric glands ($EC_{50} = 120 \mu M$), an effect mediated by histamine H₂ receptors. Betazole (3-10 mg/kg) increases pepsin and acid levels in gastric juice in dogs.²

References

1. Leth, R., Elander, B., Haglund, U., *et al.* Histamine H₂-receptor of human and rabbit parietal cells. *Am. J. Physiol.* **253**(Pt. 1), G497-G501 (1987).
2. Ishimori, A., Sakurada, H., Kawamura, T., *et al.* Comparison of dose-response curves between acid and pepsin to betazole hydrochloride in the dog with gastric fistula. *Gastroenterol. Jpn.* **10**(3), 221-224 (1975).

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, 10/03/2024

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