

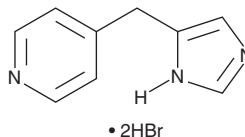
PRODUCT INFORMATION



Immethridine (hydrobromide)

Item No. 29502

CAS Registry No.: 699020-93-4
Formal Name: 4-(1H-imidazol-5-ylmethyl)-
pyridine, dihydrobromide
MF: $C_9H_9N_3 \cdot 2HBr$
FW: 321.0
Purity: $\geq 98\%$
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

Immethridine (hydrobromide) is supplied as a solid. A stock solution may be made by dissolving the immethridine (hydrobromide) in the solvent of choice, which should be purged with an inert gas. Immethridine (hydrobromide) is soluble in organic solvents such as DMSO. Immethridine (hydrobromide) is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Immethridine is a potent histamine H_3 receptor agonist ($K_i = 0.85$ nM).¹ It is selective for histamine H_3 over histamine H_4 receptors ($K_i = 245$ nM). Immethridine inhibits cAMP-stimulated β -galactosidase transcription in SK-N-MC cells expressing human H_3 receptors ($EC_{50} = 0.18$ nM) and induces a concentration-dependent decrease of the electrically induced twitch contraction in isolated guinea pig ileum. *In vivo*, immethridine (30 mg/kg, s.c.) inhibits hydrochloric acid-induced formation of gastric lesions in rats.²

References

1. Kitbunnadaj, R., Zuiderveld, O.P., Christophe, B., et al. Identification of 4-(1H-imidazol-4(5)-ylmethyl)pyridine (immethridine) as a novel, potent, and highly selective histamine H_3 receptor agonist. *J. Med. Chem.* **47(10)**, 2414-2417 (2004).
2. Adami, M., Pozzoli, C., Leurs, R., et al. Histamine H_3 receptors are involved in the protective effect of ghrelin against HCl-induced gastric damage in rats. *Pharmacology* **86(5-6)**, 259-266 (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, 12/09/2022

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