

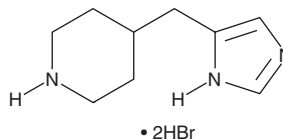
PRODUCT INFORMATION



Immepip (hydrobromide)

Item No. 29518

CAS Registry No.: 164391-47-3
Formal Name: 4-(1H-imidazol-5-ylmethyl)-
piperidine, dihydrobromide
MF: $C_9H_{15}N_3 \cdot 2HBr$
FW: 327.1
Purity: $\geq 95\%$
Supplied as: A crystalline 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

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

Description

Immepip is a histamine H_3 receptor agonist ($K_i = 0.4$ nM in SK-N-MC cell membranes expressing the human receptor).¹ It is selective for histamine H_3 over H_1 and H_2 receptors in CHO cell membranes expressing the guinea pig and human receptors, respectively (K_i s = >16 μM for both) but also agonizes H_4 receptors ($K_i = 9$ nM in SK-N-MC cell membranes expressing the human receptor).^{1,2} Immepip (5 mg/kg, s.c.) decreases hypothalamic histamine release in anesthetized rats.³ It decreases flinching in the formalin test in rats when administered at doses of 5 and 30 mg/kg and inhibits formalin-induced paw edema at 30 mg/kg.⁴

References

1. Liu, C., Ma, X., Jiang, X., *et al.* Cloning and pharmacological characterization of a fourth histamine receptor (H_4) expressed in bone marrow. *Mol. Pharmacol.* **59**(3), 420-426 (2001).
2. Vollinga, R.C., de Koning, J.P., Jansen, F.P., *et al.* A new potent and selective histamine H_3 receptor agonist, 4-(1H-imidazol-4-ylmethyl)piperidine. *J. Med. Chem.* **37**(3), 332-333 (1994).
3. Jansen, F.P., Mochizuki, T., Yamamoto, Y., *et al.* In vivo modulation of rat hypothalamic histamine release by the histamine H_3 receptor ligands, immepip and clobenpropit. Effects of intrahypothalamic and peripheral application. *Eur. J. Pharmacol.* **362**(2-3), 149-155 (1998).
4. Cannon, K.E., Leurs, R., and Hough, L.B. Activation of peripheral and spinal histamine H_3 receptors inhibits formalin-induced inflammation and nociception, respectively. *Pharmacol. Biochem. Behav.* **88**(1), 122-129 (2007).

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