

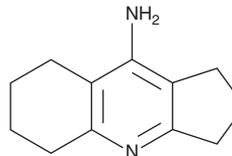
PRODUCT INFORMATION



Ipidacrine

Item No. 34529

CAS Registry No.: 62732-44-9
Formal Name: 2,3,5,6,7,8-hexahydro-1H-cyclopenta[b]quinolin-9-amine
Synonym: NIK-247
MF: C₁₂H₁₆N₂
FW: 188.3
Purity: ≥98%
UV/Vis.: λ_{max}: 215, 251 nm
Supplied as: A 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

Ipidacrine is supplied as a solid. A stock solution may be made by dissolving the ipidacrine in the solvent of choice, which should be purged with an inert gas. Ipidacrine is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ipidacrine in these solvents is approximately 25 mg/ml.

Ipidacrine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ipidacrine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Ipidacrine 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

Ipidacrine is an inhibitor of acetylcholinesterase (AChE; IC₅₀ = 270 nM).¹ It induces long-term potentiation (LTP) in rat hippocampal slices, an effect that can be blocked by the muscarinic antagonists atropine (Item No. 12008) or pirenzepine (Item No. 29527), when used at a concentration of 10 μM.² Ipidacrine (0.03-1 mg/kg) prevents memory deficits induced by the anticholinergic agent scopolamine in the passive avoidance test in rats.²

References

1. Ogura, H., Kosasa, T., Kuriya, Y., *et al.* Comparison of inhibitory activities of donepezil and other cholinesterase inhibitors on acetylcholinesterase and butyrylcholinesterase in vitro. *Methods Find. Exp. Clin. Pharmacol.* **22(8)**, 609-613 (2000).
2. Yoshida, S. and Suzuki, N. Antiamnesic and cholinomimetic side-effects of the cholinesterase inhibitors, physostigmine, tacrine and NIK-247 in rats. *Eur. J. Pharmacol.* **250(1)**, 117-124 (1993).

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, 11/22/2022

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