

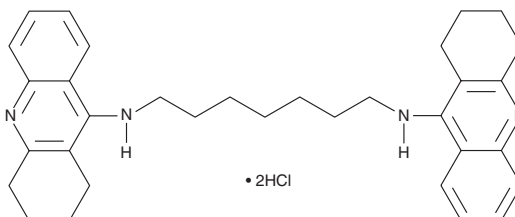
Product Information



bis(7)-Tacrine

Catalog No. 10005836

CAS Registry No.: 224445-12-9
Formal Name: N,N'-*bis*(1,2,3,4-tetrahydro-9-acridinyl)-1,7-heptanediamine, dihydrochloride
Synonym: 1,7-N-heptylene-*bis*-9,9'-amino-1,2,3,4-tetrahydro-acridine
MF: C₃₃H₄₀N₄ • 2HCl
FW: 565.6
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that *bis*(7)-tacrine be stored as supplied at -20°C. It should be stable for at least two years.

bis(7)-Tacrine is supplied as a crystalline solid. A stock solution may be made by dissolving the *bis*(7)-tacrine in an organic solvent purged with an inert gas. *bis*(7)-Tacrine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of *bis*(7)-tacrine in these solvents is approximately 20 mg/ml.

bis(7)-Tacrine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, *bis*(7)-tacrine should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. *bis*(7)-Tacrine has a solubility of approximately 0.5 mg/ml in a 1:6 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Tacrine is an amino acridine compound that inhibits acetylcholinesterase (AChE), and has been proposed as a clinical treatment for Alzheimer's disease.¹ *bis*(7)-Tacrine is a tacrine dimer, linked *via* a 7-carbon alkyl spacer. It inhibits AChE with an IC₅₀ of 0.40 nM, making it more than 1,000 times more potent than tacrine.² *bis*(7)-Tacrine also protects against hydrogen peroxide induced apoptosis in rat pheochromocytoma cells.³

References

1. Giacobini, E. Cholinesterase inhibitors for Alzheimer's disease therapy: From tacrine to future applications. *Neurochem. Int.* **32**, 413-419 (1998).
2. Carlier, P.R., Han, Y.F., Chow, E.S.H., *et al.* Evaluation of short-tether *Bis*-THA AChE inhibitors. A further test of the dual binding site hypothesis. *Bioorg. Med. Chem.* **7**, 351-357 (1999).
3. Hong, W., Carlier, P.R., Wing-Lok, H., *et al.* Attenuation of scopolamine-induced deficits in navigational memory performance in rats by *bis*(7)-tacrine, a novel dimeric AChE inhibitor. *Acta Pharmacol. Sin.* **20**(3), 211-217 (1999).

Related Product

Tacrine (hydrochloride) - Cat. No. 70240

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 02/09/2010

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com