

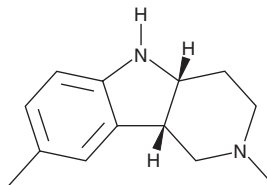
PRODUCT INFORMATION



Stobadine

Item No. 36082

CAS Registry No.: 85202-17-1
Formal Name: 2,3,4,4aR,5,9bS-hexahydro-2,8-dimethyl-1H-pyrido[4,3-b]indole
Synonym: Stobadin
MF: C₁₃H₁₈N₂
FW: 202.3
Purity: ≥98%
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

Stobadine is supplied as a solid. A stock solution may be made by dissolving the stobadine in the solvent of choice, which should be purged with an inert gas. Stobadine is soluble in the organic solvent DMSO at a concentration of approximately 100 mM.

Description

Stobadine is a pyridoindole antioxidant.¹ It scavenges hydroxyl, peroxy, and alkoxy radicals and quenches singlet oxygen in cell-free assays. *Ex vivo*, stobadine (2 mg/kg) reduces lipid peroxidation induced by hypoxia and reoxygenation in rat forebrain.² It improves survival in a rat model of carotid artery ligation-induced cerebral ischemia. Stobadine (1 mg/kg, i.v.) increases myocardial blood flow and decreases infarct area in a dog model of myocardial infarction induced by acute occlusion of the left coronary artery.¹ Dietary administration of stobadine (0.05% w/w) delays the development of cataracts and reduces eye lens protein oxidation and plasma levels of malondialdehyde (MDA) in a rat model of diabetes induced by streptozotocin (STZ; Item No. 13104).³

References

1. Horáková, L. and Štolc, S. Antioxidant and pharmacodynamic effects of pyridoindole stobadine. *Gen. Pharmacol.* **30**(5), 627-638 (1998).
2. Štolc, S., Vlkolinský, R., and Pavlásek, J. Neuroprotection by the pyridoindole stobadine: A minireview. *Brain Res. Bull.* **42**(5), 335-340 (1997).
3. Kyselova, Z., Gajdosik, A., Gajdosikova, A., *et al.* Effect of the pyridoindole antioxidant stobadine on development of experimental diabetic cataract and on lens protein oxidation in rats: Comparison with vitamin E and BHT. *Mol. Vis.* **11**, 56-65 (2005).

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, 12/19/2022

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