

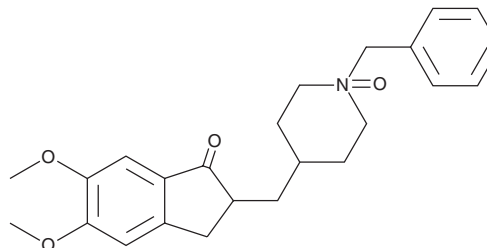
PRODUCT INFORMATION



Donepezil N-oxide

Item No. 27174

CAS Registry No.: 120013-84-5
Formal Name: 2,3-dihydro-5,6-dimethoxy-2-[[[1-oxido-1-(phenylmethyl)-4-piperidiny]methyl]-1H-inden-1-one
MF: C₂₄H₂₉NO₄
FW: 395.5
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

Donepezil N-oxide is supplied as a solid. A stock solution may be made by dissolving the donepezil N-oxide in the solvent of choice, which should be purged with an inert gas. Donepezil N-oxide is slightly soluble in chloroform and methanol.

Description

Donepezil N-oxide is an active metabolite of the acetylcholinesterase inhibitor donepezil (Item No. 13245).¹ Donepezil N-oxide inhibits cholinesterase (ChE) activity in human erythrocytes in a concentration-dependent manner, exhibiting 45.5% inhibition when used at a concentration of 20 μM.²

References

1. Groppa, F., Coin, A., De Rosa, G., *et al.* Monitoring plasma levels of donepezil, 5-O-desmethyl-donepezil, 6-O-desmethyl-donepezil, and donepezil-N-oxide by a novel HPLC method in patients with Alzheimer disease. *Ther. Drug Monit.* **38(1)**, 108-113 (2016).
2. Atto, R.A., Behnan, A.B., and Abachi, F.T. In vitro kinetic study of donepezil N-oxide metabolites *Iraq. J. Pharm.* **11(2)**, 93-101 (2011).

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, 11/28/2022

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