

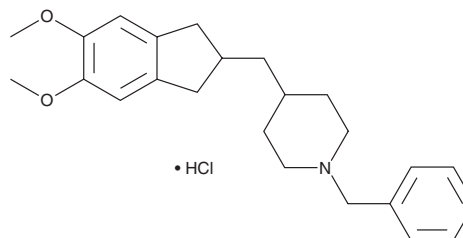
PRODUCT INFORMATION



Deoxy Donepezil (hydrochloride)

Item No. 30045

CAS Registry No.: 1034439-57-0
Formal Name: 4-[(2,3-dihydro-5,6-dimethoxy-1H-inden-2-yl)methyl]-1-(phenylmethyl)piperidine, monohydrochloride
MF: C₂₄H₃₁NO₂ • HCl
FW: 402.0
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Deoxy donepezil (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the deoxy donepezil (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Deoxy donepezil (hydrochloride) is slightly soluble in methanol.

Description

Deoxy donepezil is a potential impurity found in bulk preparations of the acetylcholinesterase inhibitor donepezil.¹ It is produced as a by-product during donepezil synthesis.²

References

1. Krishna Reddy, K.V.S.R., Moses Babu, J., Anil Kumar, P., et al. Identification and characterization of potential impurities of donepezil. *J. Pharm. Biomed. Anal.* **35(5)**, 1047-1058 (2004).
2. Krizmanic, I., Lerman, L., Samardzic, Z., et al. Impurities of donepezil. *PLIVA Hrvatska d.o.o., Zagreb (HR) US 2009/0298879 A1* (2009).

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, 02/01/2022

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