

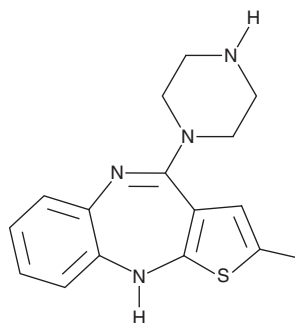
PRODUCT INFORMATION



N-desmethyl Olanzapine

Item No. 33422

CAS Registry No.: 161696-76-0
Formal Name: 2-methyl-4-(1-piperazinyl)-10H-thieno[2,3-b][1,5]benzodiazepine
Synonyms: LY170055, 4'-N-desmethyl Olanzapine
MF: C₁₆H₁₈N₄S
FW: 298.4
Purity: ≥95%
UV/Vis.: λ_{max}: 227, 272 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

N-desmethyl Olanzapine is supplied as a solid. A stock solution may be made by dissolving the N-desmethyl olanzapine in the solvent of choice, which should be purged with an inert gas. N-desmethyl Olanzapine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of N-desmethyl olanzapine in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

N-desmethyl Olanzapine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, N-desmethyl olanzapine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. N-desmethyl Olanzapine has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

N-desmethyl Olanzapine is a major metabolite of the atypical antipsychotic agent olanzapine (Item No. 11937).^{1,2} N-desmethyl Olanzapine is formed *via* oxidative metabolism of olanzapine by the cytochrome 450 (CYP) isoform CYP1A2.¹

References

1. Ring, B.J., Catlow, J., Lindsay, T.J., *et al.* Identification of the human cytochromes P450 responsible for the in vitro formation of the major oxidative metabolites of the antipsychotic agent olanzapine. *J. Pharmacol. Exp. Ther.* **276**(2), 658-666 (1996).
2. Chiu, J.A. and Franklin, R.B. Analysis and pharmacokinetics of olanzapine (LY170053) and two metabolites in rat plasma using reversed-phase HPLC with electrochemical detection. *J. Pharm. Biomed. Anal.* **14**(5), 609-615 (1996).

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, 01/03/2023

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