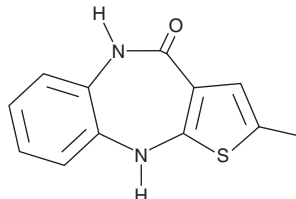


PRODUCT INFORMATION

Olanzapine-related Compound B

Item No. 33786

CAS Registry No.: 221176-49-4
Formal Name: 5,10-dihydro-2-methyl-4H-thieno[2,3-b][1,5]benzodiazepin-4-one
Synonym: LY301664
MF: C₁₂H₁₀N₂OS
FW: 230.3
Purity: ≥98%
UV/Vis.: λ_{max}: 231, 262 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

Olanzapine-related compound B is supplied as a solid. A stock solution may be made by dissolving the olanzapine-related compound B in the solvent of choice, which should be purged with an inert gas. Olanzapine-related compound B is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of olanzapine-related compound B in ethanol is approximately 15 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Olanzapine-related compound B is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, olanzapine-related compound B should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Olanzapine-related compound B has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Olanzapine-related compound B is a potential impurity found in commercial preparations of olanzapine.¹ It is a degradation product formed by exposure to oxidative stress.

Reference

1. Hiriyan, S.G., Basavaiah, K., Goud, P.S.K., *et al.* Identification and characterization of olanzapine degradation products under oxidative stress conditions. *Acta Chromatogr.* **20**(1), 81–93 (2008).

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, 11/15/2022

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