

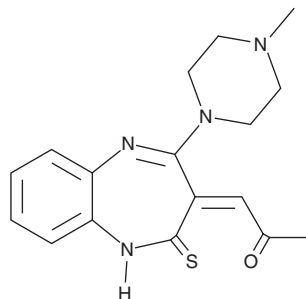
PRODUCT INFORMATION



Olanzapine Thiolactam Impurity

Item No. 27784

CAS Registry No.: 1017241-36-9
Formal Name: 1Z-[4,5-dihydro-2-(4-methyl-1-piperazinyl)-4-thioxo-3H-1,5-benzodiazepin-3-ylidene]-2-propanone
Synonym: Olanapine Ketothiolactam
MF: C₁₇H₂₀N₄OS
FW: 328.4
Purity: ≥95%
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 thiolactam impurity is supplied as a solid. A stock solution may be made by dissolving the olanzapine thiolactam impurity in the solvent of choice, which should be purged with an inert gas. Olanzapine thiolactam impurity is slightly soluble in DMSO (warmed) and methanol (warmed).

Description

Olanzapine thiolactam impurity is a potential impurity found in commercial preparations of olanzapine.¹ It is a degradation product formed during storage or exposure to thermal stress.

Reference

1. Baertschi, S.W., Brunner, H., Bunnell, C.A., *et al.* Isolation, identification, and synthesis of two oxidative degradation products of olanzapine (LY170053) in solid oral formulations. *J. Pharm. Sci.* **97**(2), 883-892 (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, 10/04/2022

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