

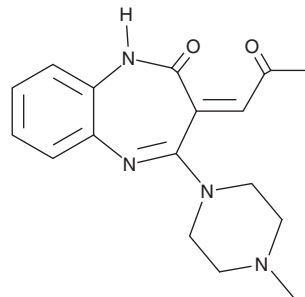
PRODUCT INFORMATION



Olanzapine Lactam Impurity

Item No. 33792

CAS Registry No.: 1017241-34-7
Formal Name: (3Z)-1,3-dihydro-4-(4-methyl-1-piperazinyl)-3-(2-oxopropylidene)-2H-1,5-benzodiazepin-2-one
MF: C₁₇H₂₀N₄O₂
FW: 312.4
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 lactam impurity is supplied as a solid. A stock solution may be made by dissolving the olanzapine lactam impurity in the solvent of choice, which should be purged with an inert gas. Olanzapine lactam impurity is slightly soluble in methanol and DMSO.

Description

Olanzapine lactam impurity is a potential impurity found in commercial preparations of olanzapine.¹ It is a degradation product formed by exposure to thermal or oxidative 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, 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