

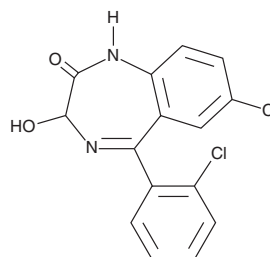
PRODUCT INFORMATION



Lorazepam

Item No. 15891

CAS Registry No.: 846-49-1
Formal Name: 7-chloro-5-(2-chlorophenyl)-1,3-dihydro-3-hydroxy-2H-1,4-benzodiazepin-2-one
Synonyms: Apo-Lorazepam, NSC 279758, Wy 4036
MF: C₁₅H₁₀Cl₂N₂O₂
FW: 321.2
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥2 years



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

Description

Lorazepam (Item No. 15891) is an analytical reference material categorized as a benzodiazepine.^{1,2} Lorazepam has anxiolytic properties and is a metabolite of diazepam (Item Nos. 28623 | 15554).^{2,3} Formulations containing lorazepam have been used in the treatment of anxiety. Lorazepam is regulated as a Schedule IV compound in the United States. This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards for reference materials.

References

1. Miller, L.G., Greenblatt, D.J., Barnhill, J.G., *et al.* Chronic benzodiazepine administration. I. Tolerance is associated with benzodiazepine receptor downregulation and decreased γ -aminobutyric acid_A receptor function. *J. Pharmacol. Exp. Ther.* **246(1)**, 170-176 (1988).
2. Baldwin, D., Woods, R., Lawson, R., *et al.* Efficacy of drug treatments for generalised anxiety disorder: Systematic review and meta-analysis. *BMJ* **342**, d1199 (2011).
3. Moosmann, B., Bisel, P., and Auwärter, V. Characterization of the designer benzodiazepine diazepam and preliminary data on its metabolism and pharmacokinetics. *Drug Test Anal.* **6(7-8)**, 757-763 (2014).

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, 03/13/2024

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