

PRODUCT INFORMATION



Lormetazepam

Item No. 21989

CAS Registry No.: 848-75-9

Formal Name: 7-chloro-5-(2-chlorophenyl)-1,3-dihydro-3-hydroxy-1-methyl-2H-1,4-benzodiazepin-2-one

Synonyms: N-methyl Lorazepam, Ro 5-5516, Wy 4082

MF: $C_{16}H_{12}Cl_2N_2O_2$

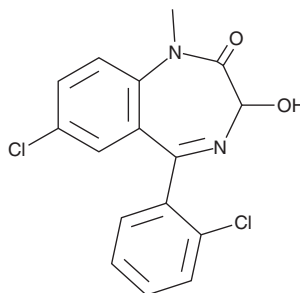
FW: 335.2

Purity: $\geq 98\%$

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥ 5 years



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

Description

Lormetazepam (Item No. 21989) is an analytical reference standard categorized as a benzodiazepine.¹ Lormetazepam is an active metabolite of diclazepam (Item Nos. 28623 | 15554) and is associated with misuse.^{2,3} Formulations containing lormetazepam have been used in the treatment of insomnia. Lormetazepam is regulated as a Schedule IV compound in the United States. This product is intended for research and forensic applications.

References

1. Hümpel, M., Illi, V., Milius, W., *et al.* The pharmacokinetics and biotransformation of the new benzodiazepine lormetazepam in humans. I. Absorption, distribution, elimination and metabolism of lormetazepam-5-¹⁴C. *Eur. J. Drug Metab. Pharmacokinet.* **4(4)**, 237-243 (1979).
2. Moosmann, B., Bisel, P., and Auwärter, V. Characterization of the designer benzodiazepine diclazepam and preliminary data on its metabolism and pharmacokinetics. *Drug Test Anal.* **6(7-8)**, 757-763 (2014).
3. Magro, L., Faccini, M., and Leone, R. Lormetazepam addiction. *Neuropathology of drug addictions and substance misuse*. Preedy, V., editor, 1st edition, Academic Press (2016).

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, 06/29/2023

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