

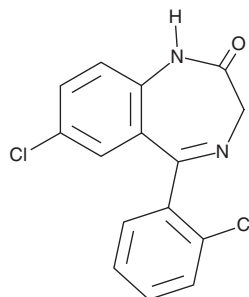
PRODUCT INFORMATION



Delorazepam

Item No. 15938

CAS Registry No.: 2894-67-9
Formal Name: 7-chloro-5-(2-chlorophenyl)-1,3-dihydro-2H-1,4-benzodiazepin-2-one
Synonyms: Chlordesmethyl diazepam, NSC 169895, Ro 5-3027
MF: C₁₅H₁₀Cl₂N₂O
FW: 305.1
Purity: ≥98%
UV/Vis.: λ_{max}: 227, 319 nm
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

Delorazepam is a benzodiazepine with anxiolytic effects.¹ It is a metabolite of the designer drug diclazepam that can be detected in either the serum or urine.² Delorazepam can also be produced by enzymatic treatment of lorazepam with β-glucuronidase.³ This compound is intended for forensic and research applications.

References

1. Paradiso, L. and Rapisarda, V. The psychosedative effect of chlordesmethyl diazepam on anxiety states in neurotic and psychotic inpatients. *Clin. Ther.* **3(3)**, 194-199 (1980).
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. Fu, S., Molnar, A., Bowron, P., et al. Reduction of temazepam to diazepam and lorazepam to delorazepam during enzymatic hydrolysis. *Anal. Bioanal. Chem.* **400(1)**, 153-164 (2011).

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