

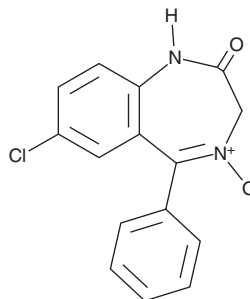
PRODUCT INFORMATION



Demoxepam

Item No. 21937

CAS Registry No.: 963-39-3
Formal Name: 7-chloro-1,3-dihydro-5-phenyl-2H-1,4-benzodiazepin-2-one, 4-oxide
Synonyms: Chlordiazepoxide Lactam, NSC 46007, NSC 46077, NSC 169898
MF: C₁₅H₁₁ClN₂O₂
FW: 286.7
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥3 years



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

Description

Demoxepam (Item No. 21937) is an analytical reference standard categorized as a benzodiazepine. Demoxepam is an active metabolite of chlordiazepoxide (Item No. 14261).^{1,2} It is also a potential impurity in commercial preparations of chlordiazepoxide.³ This product is intended for research and forensic applications.

References

1. Sidransky, H., Verney, E., Cosgrove, J.W., *et al.* Effect of benzodiazepines on tryptophan binding to rat hepatic nuclei. *Toxicol. Pathol.* **20(3 Pt 1)**, 350-356 (1992).
2. Koechlin, B.A., and D'Arconte, L. Determination of chlordiazepoxide (librium) and of a metabolite of lactam character in plasma of humans, dogs, and rats by a specific spectrofluorometric micro method. *Anal. Biochem.* **5**, 195-207 (1963).
3. French, W.N., Matsui, F.F., and Smith, S.J. Determination of major impurity in chlordiazepoxide formulations and drug substance. *J. Pharm. Sci.* **64(9)**, 1545-1547 (1975).

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, 02/01/2024

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