

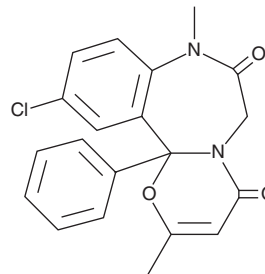
PRODUCT INFORMATION



Ketazolam

Item No. 11449

CAS Registry No.: 27223-35-4
Formal Name: 11-chloro-8,12b-dihydro-2,8-dimethyl-12b-phenyl-4H-[1,3]oxazino[3,2-d][1,4]benzodiazepine-4,7(6H)-dione
Synonyms: NSC 338158, U 28774
MF: C₂₀H₁₇ClN₂O₃
FW: 368.8
Purity: ≥95%
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

Ketazolam is a benzodiazepine derivative that was used as an anxiolytic in the 1980s.^{1,2} Tolerance within fifteen days, physical dependence, and benzodiazepine withdrawal syndrome are associated with ketazolam use.³ This product is intended for forensic purposes.

References

1. Rickels, K., Csanalosi, I., Greisman, P., *et al.* Ketazolam and diazepam in anxiety: A controlled study. *J. Clin. Pharmacol.* **20(10)**, 581-589 (1980).
2. Kim, K.K., Sirman, A., Trainor, F.S., *et al.* Anxiolytic efficacy and safety of ketazolam compared with diazepam and placebo. *Clin. Ther.* **3(1)**, 9-14 (1980).
3. Higgitt, A., Fonagy, P., and Lader, M. The natural history of tolerance to the benzodiazepines. *Psychol. Med. Monogr. Suppl.* **13**, 1-55 (1988).

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, 07/13/2023

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