

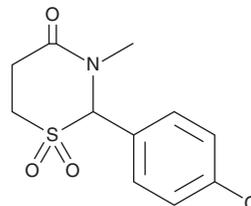
PRODUCT INFORMATION



Chlormezanone

Item No. 25716

CAS Registry No.: 80-77-3
Formal Name: 2-(4-chlorophenyl)tetrahydro-3-methyl-4H-1,3-thiazin-4-one, 1,1-dioxide
Synonyms: (±)-Chlormezanone, NSC 169108
MF: C₁₁H₁₂ClNO₃S
FW: 273.7
Purity: ≥98%
UV/Vis.: λ_{max}: 228 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Chlormezanone is supplied as a crystalline solid. A stock solution may be made by dissolving the chlormezanone in the solvent of choice, which should be purged with an inert gas. Chlormezanone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of chlormezanone in these solvents is approximately 2, 20, and 10 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of chlormezanone can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of chlormezanone in PBS, pH 7.2, is approximately 0.1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Chlormezanone is a centrally acting muscle relaxant.¹ It induces paralysis in mice, guinea pigs, dogs, cats, and monkeys (EC₅₀s = 133, <200, 100-250, 50, and 50 mg/kg, respectively). Chlormezanone increases ataxia induced by pentobarbital and alcohol in mice and blocks the crossed extensor reflex of the spine without affecting the knee-jerk reflex in cats. It also induces head drop in rabbits.

Reference

1. Gesler, R.M. and Coulston, F. Pharmacologic and toxicologic studies of two new centrally acting skeletal muscle relaxants, chlormethazanone and dichlormethazanone. *Toxicol. Appl. Pharmacol.* **1(2)**, 168-184 (1959).

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, 10/06/2022

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