

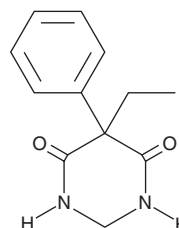
PRODUCT INFORMATION



Primidone

Item No. 19277

CAS Registry No.: 125-33-7
Formal Name: 5-ethylidihydro-5-phenyl-4,6(1H,5H)-pyrimidinedione
Synonym: NSC 41701
MF: C₁₂H₁₄N₂O₂
FW: 218.3
Purity: ≥98%
UV/Vis.: λ_{max}: 206 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

Primidone (Item No. 19277) is an analytical reference standard categorized as a barbiturate. Primidone partially substitutes for the barbiturate barbitone and reduces barbitone withdrawal-induced convulsions in rats in a model of barbitone dependence.¹ Formulations containing primidone have been used in the treatment of various seizure types. This product is intended for research and forensic applications.

Reference

1. Norton, P.R. The effects of drugs on barbiturate withdrawal convulsions in the rat. *J. Pharm. Pharmacol.* **22(10)**, 763-766 (1970).

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, 01/11/2024

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