

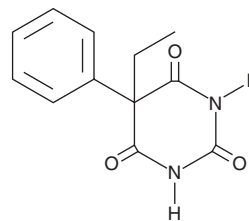
PRODUCT INFORMATION



Phenobarbital

Item No. 9001494

CAS Registry No.: 50-06-6
Formal Name: 5-ethyl-5-phenyl-2,4,6(1H,3H,5H)-pyrimidinetrione
Synonyms: NSC 9848, NSC 128143, Phenobarbitone
MF: C₁₂H₁₂N₂O₃
FW: 232.2
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥2 years



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

Description

Phenobarbital (Item No. 9001494) is an analytical reference material categorized as a barbiturate. Phenobarbital is an active metabolite of primidone (Item Nos. 40161 | 19277).^{1,2} It has anticonvulsant activity in rats.² Phenobarbital has been found as an adulterant in cocaine.³ Phenobarbital is regulated as a Schedule IV compound in the United States. This product is intended for research and forensic applications.

This product qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

References

1. Nagaki, S., Ratnaraj, N., and Patsalos, P.N. Blood and cerebrospinal fluid pharmacokinetics of primidone and its primary pharmacologically active metabolites, phenobarbital and phenylethylmalonamide in the rat. *Eur. J. Drug Metab. Pharmacokinet.* **24(3)**, 255-264 (1999).
2. Chiu, P. and Burnham, W.M. The effect of anticonvulsant drugs on convulsions triggered by direct stimulation of the brainstem. *Neuropharmacology* **21(4)**, 355-359 (1982).
3. Fucci, N. and De Giovanni, N. Adulterants encountered in the illicit cocaine market. *Forensic Sci. Int.* **95(3)**, 247-252 (1998).

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, 02/29/2024

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