

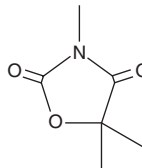
# PRODUCT INFORMATION



## Trimethadione

Item No. 23893

CAS Registry No.: 127-48-0  
Formal Name: 3,5,5-trimethyl-2,4-oxazolidinedione  
Synonyms: NSC 15799, NSC 169503  
MF:  $C_6H_9NO_3$   
FW: 143.1  
Purity:  $\geq 98\%$   
Supplied as: A solid  
Storage:  $-20^\circ\text{C}$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

Trimethadione is supplied as a solid. A stock solution may be made by dissolving the trimethadione in the solvent of choice, which should be purged with an inert gas. Trimethadione is slightly soluble in chloroform and methanol.

### Description

Trimethadione is an anticonvulsant.<sup>1,2</sup> It decreases the levels of  $\alpha$ -oxoglutarate, malate, and fumarate in the brain as well as plasma glucose levels in rats.<sup>1</sup> Trimethadione increases latency to first pentetrazole-induced clonic seizure in mice ( $ED_{50} = 10 \text{ mg/kg}$ ).<sup>2</sup> It also inhibits absence-like seizures but not tonic convulsions in spontaneously epileptic rats.<sup>3</sup>

### References

1. Nahorski, S.R. Biochemical effects of the anticonvulsants trimethadione, ethosuximide and chlordiazepoxide in rat brain. *J. Neurochem.* **19(8)**, 1937-1946 (1972).
2. Frey, H.H. Determination of the anticonvulsant potency of unmetabolized trimethadione. *Acta. Pharmacol. Toxicol. (Copenh)* **27(4)**, 295-300 (1969).
3. Sasa, M., Ohno, Y., Ujihara, H., et al. Effects of antiepileptic drugs on absence-like and tonic seizures in the spontaneously epileptic rat, a double mutant rat. *Epilepsia* **29(5)**, 505-513 (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, 10/17/2022

#### 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