

PRODUCT INFORMATION



Cambendazol

Item No. 20841

CAS Registry No.: 26097-80-3
Formal Name: N-[2-(4-thiazolyl)-1H-benzimidazol-6-yl]-carbamic acid, 1-methylethyl ester

Synonyms: Isopropyl 2-(4-thiazolyl)-5-benzimidazolecarbamate, MK-905, NSC 377071

MF: C₁₄H₁₄N₄O₂S

FW: 302.4

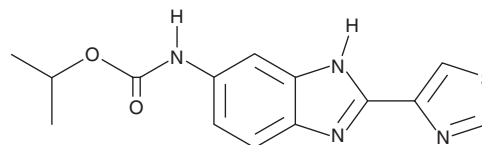
Purity: ≥98%

UV/Vis.: λ_{max}: 235, 324 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

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

Description

Cambendazol is an antihelmintic that, at a dose of 50 mg/kg/day in mice, eradicates *S. ratti* adult worms from intestine and *S. stercoralis* larva from muscle.¹ It inhibits polymerization of microtubules isolated from bovine brain with an IC₅₀ value of 64.2 μM.² Formulations containing cambendazol have been used to treat strongyloidiasis in humans and have been studied as potential chemotherapeutics.^{3,4}

References

1. Grove, D. I. *Strongyloides ratti* and *S. stercoralis*: The effects of thiabendazole, mebendazole, and cambendazole in infected mice. *Am. J. Trop. Med. Hyg.* **31(3 Pt 1)**, 469-476 (1982).
2. Friedman, P. A. and Platzer, E.G. Interaction of anthelmintic benzimidazoles and benzimidazole derivatives with bovine brain tubulin. *Biochim. Biophys. Acta.* **544(3)**, 605-614 (1978).
3. Bichalho, S.A., Leão, O.J., and Pena, Q. Jr. Cambendazole in the treatment of human strongyloidiasis. *Am. J. Trop. Med. Hyg.* **32(5)**, 1181-1183 (1983).
4. Dunn, W.A. Jr., Akin, D.E., Progulske-Fox, A., et al. Methods and compositions for the treatment of cancers and pathogenic infections. *Patent Application Publication*. US20100285012. 1-7 (2010).

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, 10/20/2022

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