

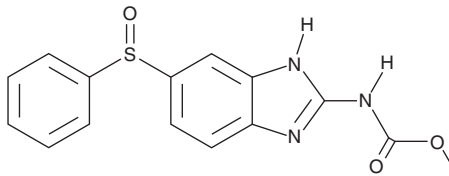
PRODUCT INFORMATION



Oxfendazole

Item No. 29742

CAS Registry No.: 53716-50-0
Formal Name: N-[6-(phenylsulfinyl)-1H-benzimidazol-2-yl]-carbamic acid, methyl ester
Synonym: Oxfenbendazole, Fenbendazole sulfoxide, HOE 8051
MF: C₁₅H₁₃N₃O₃S
FW: 315.3
Purity: ≥98%
UV/Vis.: λ_{max}: 229, 297 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

Oxfendazole is supplied as a crystalline solid. A stock solution may be made by dissolving the oxfendazole in the solvent of choice, which should be purged with an inert gas. Oxfendazole is soluble in organic solvents DMSO and dimethyl formamide (DMF). The solubility of oxfendazole in DMSO is approximately 1 mg/ml. Oxfendazole is slightly soluble in DMF.

Oxfendazole is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, oxfendazole should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Oxfendazole has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Oxfendazole is a benzimidazole anthelmintic.^{1,2} *In vivo*, oxfendazole is curative in pig models of *H. rubidus* or *A. suum* infection when administered at doses of 3 and 6 mg/kg, respectively.¹ It is lethal to *H. contortus* and *T. circumcincta* when administered to infected cattle (LD₅₀s = 4.28 and 3.61 mg/kg of host weight, respectively).² Formulations containing oxfendazole have been used in the treatment and prevention of helminth infections in livestock.

References

1. Kingsbury, P.A., Rowlands, D.T., Reid, J.F., *et al.* Anthelmintic activity of oxfendazole in pigs. *Vet. Rec.* **108(1)**, 10-11 (1981).
2. Le Jambre, L.F., Prichard, P.K., Hennessy, D.R., *et al.* Efficiency of oxfendazole administered as a single dose or in a controlled release capsule against benzimidazole-resistant *Haemonchus contortus*, *Ostertagia circumcincta* and *Trichostrongylus colubriformis*. *Res. Vet. Sci.* **31(3)**, 289-294 (1981).

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, 11/03/2022

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