

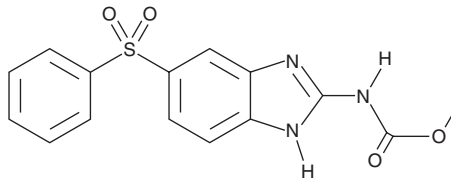
PRODUCT INFORMATION



Fenbendazole sulfone

Item No. 20921

CAS Registry No.: 54029-20-8
Formal Name: N-[6-(phenylsulfonyl)-1H-benzimidazol-2-yl]-
carbamic acid, methyl ester
Synonyms: HOE 5151, Oxfendazole
MF: C₁₅H₁₃N₃O₄S
FW: 331.4
Purity: ≥98%
UV/Vis.: λ_{max}: 290 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

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

Description

Fenbendazole sulfone is a minor metabolite of the anthelmintic fenbendazole (Item No. 19687).¹ Fenbendazole undergoes oxidation to form oxfendazole, which is further oxidized to form fenbendazole sulfone.

Reference

1. Short, C.R., Flory, W., Hsieh, L.C., *et al.* The oxidative metabolism of fenbendazole: A comparative study. *J. Vet. Pharmacol. Ther.* **11(1)**, 50-55 (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, 10/18/2022

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