

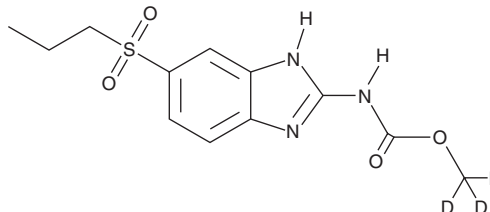
PRODUCT INFORMATION



Albendazole sulfone-d₃

Item No. 35129

CAS Registry No.: 1448345-60-5
Formal Name: N-[6-(propylsulfonyl)-1H-benzimidazol-2-yl]-carbamic acid, methyl-d₃ ester
Synonyms: ABZ-SO₂-d₃, ABZ-SOO-d₃
MF: C₁₂H₁₂D₃N₃O₄S
FW: 300.3
Chemical Purity: ≥90% (Albendazole sulfone)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
UV/Vis.: λ_{max}: 223 nm
Supplied as: A 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

Albendazole sulfone-d₃ is intended for use as an internal standard for the quantification of albendazole sulfone (Item No. 35445) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Albendazole sulfone-d₃ is supplied as a solid. A stock solution may be made by dissolving the albendazole sulfone-d₃ in the solvent of choice, which should be purged with an inert gas. Albendazole sulfone-d₃ is soluble in the organic solvent dimethyl formamide at a concentration of approximately 1 mg/ml. It is also slightly soluble in DMSO.

Description

Albendazole sulfone is an active metabolite of the benzimidazole anthelmintic albendazole (Item No. 23705).^{1,2} It inhibits spore formation of *E. hellem*, *E. cuniculi*, and *E. intestinalis* *in vitro* (IC₉₀S = <0.0001, <0.001, and <0.1 µg/ml, respectively).²

References

1. Baliharová, V., Velík, J., Fimanová, K., *et al.* Inhibitory effect of albendazole and its metabolites on cytochromes P450 activities in rat and mouflon *in vitro*. *Pharmacol. Rep.* **57**(1), 97-106 (2005).
2. Ridoux, O. and Drancourt, M. In vitro susceptibilities of the microsporidia *Encephalitozoon cuniculi*, *Encephalitozoon hellem*, and *Encephalitozoon intestinalis* to albendazole and its sulfoxide and sulfone metabolites. *Antimicrob. Agents and Chemother.* **42**(12), 3301-3303 (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, 12/12/2022

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