

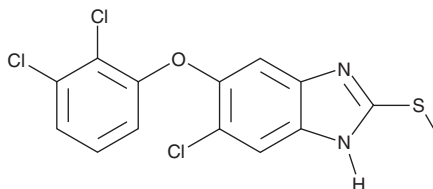
PRODUCT INFORMATION



Triclabendazole

Item No. 18883

CAS Registry No.: 68786-66-3
Formal Name: 6-chloro-5-(2,3-dichlorophenoxy)-2-(methylthio)-1H-benzimidazole
Synonym: TCBZ
MF: C₁₄H₉Cl₃N₂OS
FW: 359.7
Purity: ≥98%
UV/Vis.: λ_{max}: 222, 254, 306 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

Triclabendazole is supplied as a crystalline solid. A stock solution may be made by dissolving the triclabendazole in the solvent of choice, which should be purged with an inert gas. Triclabendazole is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

Description

Triclabendazole is a benzimidazole anthelmintic.¹ It eliminates both immature and adult flukes in a sheep model of *F. hepatica* infection when administered at a dose of 10 mg/kg.² Triclabendazole (5 μM) also protects against glucose- or α-synuclein-induced apoptosis in *S. cerevisiae* via inhibition of adenylyl cyclase in the Ras-adenylyl cyclase-protein kinase A (PKA) nutrient sensing pathway.³

References

1. Shrifi, Y., Farahnak, A., Golestani, A., *et al.* Triclabendazole effect on protease enzyme activity in the excretory- secretory products of *Fasciola hepatica* *in vitro*. *Iran. J. parasitol.* **9(1)**, 107-113 (2014).
2. Smeal, M.G. and Hall, C.A. The activity of triclabendazole against immature and adult *Fasciola hepatica* infections in sheep. *Aust. Vet. J.* **60(11)**, 329-331 (1983).
3. Lee, Y.J., Burlet, E., Wang, S., *et al.* Triclabendazole protects yeast and mammalian cells from oxidative stress: Identification of a potential neuroprotective compound. *Biochem. Biophys. Res. Commun.* **414(1)**, 205-208 (2011).

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, 10/05/2022

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