

PRODUCT INFORMATION

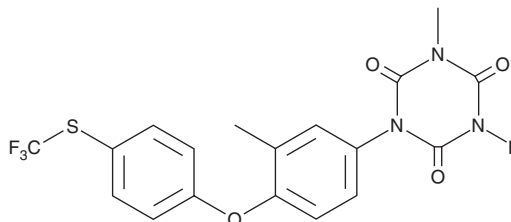


Toltrazuril

Item No. 20767

CAS Registry No.: 69004-03-1
Formal Name: 1-methyl-3-[3-methyl-4-[4-[(trifluoromethyl)thio]phenoxy]phenyl]-1,3,5-triazine-2,4,6(1H,3H,5H)-trione

Synonyms: BAY-Vi-9142
MF: $C_{18}H_{14}F_3N_3O_4S$
FW: 425.4
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 242 nm
Supplied as: A crystalline solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Toltrazuril is supplied as a crystalline solid. A stock solution may be made by dissolving the toltrazuril in the solvent of choice, which should be purged with an inert gas. Toltrazuril is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of toltrazuril in ethanol is approximately 1 mg/ml and approximately 25 mg/ml in DMSO and DMF.

Toltrazuril is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, toltrazuril should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Toltrazuril 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

Toltrazuril is an antiprotozoal agent that is widely used to prevent *coccidia* parasite infection in veterinary applications.¹⁻⁴ The effectiveness of treatment appears to depend on the coccidian species, development stage of the parasite, and level of infection.^{1,4}

References

1. Alnassan, A.A., Shehata, A.A., Kotsch, M., *et al.* Efficacy of early treatment with toltrazuril in prevention of coccidiosis and necrotic enteritis in chickens. *Avian Pathol.* **42**(5), 482-490 (2013).
2. Bosco, A., Rinaldi, L., Cappelli, G., *et al.* Metaphylactic treatment strategies with toltrazuril and diclazuril and growth performance of buffalo calves exposed to a natural eimeria infection. *Vet. Parasitol.* **212**(3-4), 408-410 (2015).
3. Zechner, G., Bauer, C., Jacobs, J.R., *et al.* Efficacy of diclazuril and toltrazuril in the prevention of coccidiosis in dairy calves under field conditions. *Vet. Rec.* **176**(5), 126 (2015).
4. Sokół, R., Gesek, M., Raś-Noryńska, M., *et al.* Toltrazuril (Baycox) treatment against coccidiosis caused by *Eimeria* sp. in Japanese quails (*Coturnix coturnix japonica*). *Pol. J. Vet. Sci.* **17**(3), 465-468 (2014).

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