

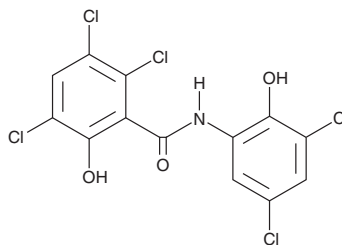
PRODUCT INFORMATION



Oxyclozanide

Item No. 46177

CAS Registry No.: 2277-92-1
Formal Name: 2,3,5-trichloro-N-(3,5-dichloro-2-hydroxyphenyl)-6-benzamide
MF: $C_{13}H_6Cl_5NO_3$
FW: 401.5
Purity: $\geq 98\%$
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

Oxyclozanide is supplied as a solid. A stock solution may be made by dissolving the oxyclozanide in the solvent of choice, which should be purged with an inert gas. Oxyclozanide is soluble (≥ 10 mg/ml) in ethanol and DMSO.

Description

Oxyclozanide is a salicylanilide anthelmintic.^{1,2} *In vitro*, it uncouples oxidative phosphorylation in rat liver mitochondria and increases oxygen uptake and death in sheep liver flukes (*F. hepatica*).¹ Oxyclozanide (22.5 mg/kg) decreases the survival of rumen fluke (*C. daubneyi*) larvae and adult worms by 82 and 95.9%, respectively, in goats.² Formulations containing oxyclozanide have been used in the treatment of fascioliasis in cattle, sheep, and goats.

References

1. Corbett, J.R. and Goose, A.J. The biochemical mode of action of the fasciolicides nitroxylin, hexachlorophane and oxyclozanide. *Biochem. J.* **121**(3), (1971).
2. Paraud, C., Gaudin, C., Pors, I., *et al.* Efficacy of oxyclozanide against the rumen fluke *Calicophoron daubneyi* in experimentally infected goats. *Vet. J.* **180**(2), 265-267 (2009).

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, 07/24/2026

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