

PRODUCT INFORMATION



Closantel

Item No. 20531

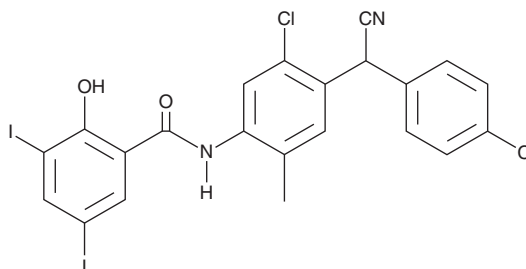
CAS Registry No.: 57808-65-8
Formal Name: N-[5-chloro-4-[(4-chlorophenyl)cyanomethyl]-2-methylphenyl]-2-hydroxy-3,5-diiodo-benzamide

Synonym: NSC 335306
MF: C₂₂H₁₄Cl₂N₂O₂
FW: 663.1
Purity: ≥98%
UV/Vis.: λ_{max}: 222, 333 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



Laboratory Procedures

Closantel is supplied as a crystalline solid. A stock solution may be made by dissolving the closantel in the solvent of choice. Closantel is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of closantel in these solvents is approximately 20 mg/ml.

Closantel is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, closantel should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Closantel has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Closantel is a salicylanilide compound with narrow-spectrum anthelmintic actions. It is effective against certain roundworms, flukes, and myiasis but is not effective against tapeworms or most external parasites.¹

Reference

1. Vieira, J.N., Filho, F.S.M., Ferreira, G.F., *et al.* In vitro susceptibility of nematophagous fungi to antiparasitic drugs: Interactions and implications for biological control. *Braz. J. Biol.* (2016).

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, 11/28/2016

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