

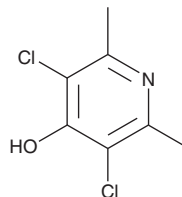
PRODUCT INFORMATION



Clopidol

Item No. 26800

CAS Registry No.: 2971-90-6
Formal Name: 3,5-dichloro-2,6-dimethyl-4-pyridinol
Synonyms: Farmcoccid, Meticlorpindol, NSC 253479, WR 61112
MF: C₇H₇Cl₂NO
FW: 192.0
Purity: ≥98%
UV/Vis.: λ_{max}: 210, 270 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

Clopidol is supplied as a crystalline solid. A stock solution may be made by dissolving the clopidol in the solvent of choice, which should be purged with an inert gas. Clopidol is sparingly soluble in 10 mM NaOH.

Description

Clopidol is a coccidiostat.¹ It prevents mortality of chickens infected with the parasites *E. tenella*, *E. necatrix*, *E. acervuline*, *E. mivati*, *E. brunetti*, and *E. maxima*. It has antimalarial activity against *P. berghei* in mice and *P. cynomologi* in rhesus macaques.² Clopidol is a potential contaminant in food poultry.³ Formulations containing clopidol have been used in the prevention of coccidiosis in chickens.

References

1. Reid, W.M. and Brewer, R.N. Efficacy studies on meticlorpindol as a coccidiostat. *Poultry Sci.* **46(3)**, 638-642 (1967).
2. Markley, L.D., Van Heertum, J.C., and Doorenbos, H.E. Antimalarial activity of clopidol, 3,5-dichloro-2,6-dimethyl-4-pyridinol, and its esters, carbonates, and sulfonates. *J. Med. Chem.* **15(11)**, 1188-1189 (1972).
3. MacNeil, J.D. Committee on drugs and related topics. Drug residues in animal tissues. *J. AOAC Int.* **86(1)**, 116-127 (2003).

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, 10/31/2022

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