

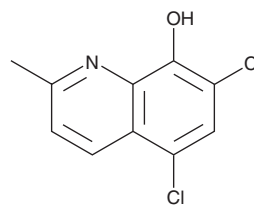
PRODUCT INFORMATION



Chlorquinaldol

Item No. 29396

CAS Registry No.: 72-80-0
Formal Name: 5,7-dichloro-2-methyl-8-quinolinol
MF: C₁₀H₇Cl₂NO
FW: 228.1
Purity: ≥98%
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

Chlorquinaldol is supplied as a crystalline solid. A stock solution may be made by dissolving the chlorquinaldol in the solvent of choice, which should be purged with an inert gas. Chlorquinaldol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of chlorquinaldol in these solvents is approximately 5, 20, and 30 mg/ml, respectively.

Description

Chlorquinaldol is a monohydroxyquinoline antimicrobial agent.¹ It is active against a variety of fungi (MICs = 1-10 µg/ml), as well as clinical isolates of *S. aureus*, *S. epidermidis*, *E. faecalis*, *S. pyogenes*, and *P. acnes* Gram-positive (MICs = 0.016-32 mg/L) and *E. coli*, *P. mirabilis*, *P. aeruginosa*, and *Enterobacter* Gram-negative bacteria (MICs = 8-512 mg/L).^{1,2} Chlorquinaldol (100 µg/ml) is also active against clinical isolates of *Staphylococcus* from bovine mastitis.³

References

1. Littman, M.L. Antimycotic effect of chlorquinaldol. *Trans. N. Y. Acad. Sci.* **18(2)**, 161-174 (1955).
2. Bortolin, M., Bidossi, A., De Vecchi, E., *et al.* *In vitro* antimicrobial activity of chlorquinaldol against microorganisms responsible for skin and soft tissue infections: Comparative evaluation with gentamicin and fusidic acid. *Front. Microbiol.* **8**, 1039 (2017).
3. Fratta, I., Mann, P.H., and Sigg, E.B. Bacteriophage typing and sensitivity testing to chlorquinaldol of Staphylococci isolated from bovine mastitis. *Can. J. Comp. Med. Vet. Sci.* **25(4)**, 96-97 (1961).

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, 11/15/2022

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