

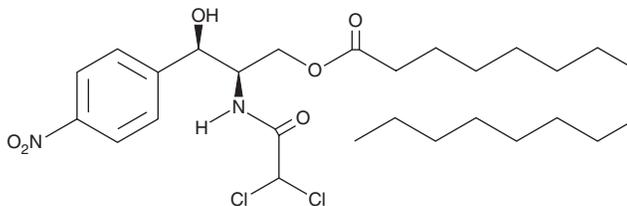
PRODUCT INFORMATION



Chloramphenicol Palmitate

Item No. 26663

CAS Registry No.: 530-43-8
Formal Name: hexadecanoic acid, (2R,3R)-2-[(2,2-dichloroacetyl)amino]-3-hydroxy-3-(4-nitrophenyl)propyl ester
MF: C₂₇H₄₂Cl₂N₂O₆
FW: 561.5
Purity: ≥99%
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

Chloramphenicol palmitate is supplied as a solid. A stock solution may be made by dissolving the chloramphenicol palmitate in the solvent of choice, which should be purged with an inert gas. Chloramphenicol palmitate is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Chloramphenicol palmitate is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Chloramphenicol palmitate is an orally bioavailable ester prodrug form of the antibiotic chloramphenicol.¹ It is hydrolyzed in the small intestine to release chloramphenicol. Formulations containing chloramphenicol palmitate were previously used in the treatment of severe bacterial infections.

Reference

1. Ambrose, P.J. Clinical pharmacokinetics of chloramphenicol and chloramphenicol succinate. *Clin. Pharmacokinet.* **9(3)**, 222-238 (1984).

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, 12/21/2022

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