

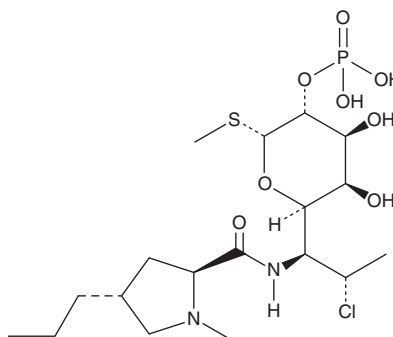
PRODUCT INFORMATION



Clindamycin Phosphate

Item No. 35611

CAS Registry No.: 24729-96-2
Formal Name: methyl 7-chloro-6,7,8-trideoxy-6-[[[(2S,4R)-1-methyl-4-propyl-2-pyrrolidinyl]carbonyl]amino]-1-thio-L-threo- α -D-galacto-octopyranoside, 2-(dihydrogen phosphate)
Synonyms: Clindamycin 2-Phosphate, DARE-BV1, NSC 618653, U-28508
MF: $C_{18}H_{34}ClN_2O_8PS$
FW: 505.0
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

Clindamycin phosphate is supplied as a solid. A stock solution may be made by dissolving the clindamycin phosphate in the solvent of choice, which should be purged with an inert gas. Clindamycin phosphate is soluble in the organic solvent DMSO at a concentration of approximately 2 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of clindamycin phosphate can be prepared by directly dissolving the solid in aqueous buffers. The solubility of clindamycin phosphate in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Clindamycin phosphate is a lincosamide antibiotic.¹ It is active against *P. acnes* *in vitro* (MIC = 0.02 $\mu\text{g/ml}$). Topical administration of clindamycin phosphate (1% w/w) reduces the number of bacteria at the burn site in a mouse model of *P. acnes* burn infection. Formulations containing clindamycin phosphate have been used in the treatment of acne.

Reference

1. Komagata, Y., Komiyama, K., and Nomura, S. Fundamental studies on antibacterial activity of clindamycin against *Propionibacterium acnes*. *Jpn. J. Antibiot.* **51**(2), 130-136 (1998).

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, 12/12/2022

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