

PRODUCT INFORMATION



Clindamycin 3-phosphate

Item No. 46097

CAS Registry No.: 28708-34-1

Formal Name: (2S-*trans*)-methyl 7-chloro-6,7,8 trideoxy-6-[[[(1-methyl-4-propyl-2-pyrrolidinyl) carbonyl]amino]-1-thio-L-*threo*- α -D-galacto-octopyranoside, 3-(dihydrogen phosphate)

MF: $C_{18}H_{34}ClN_2O_8PS$

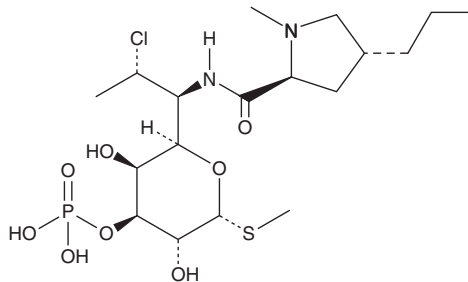
FW: 505.0

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Clindamycin 3-phosphate is supplied as a solid. A stock solution may be made by dissolving the clindamycin 3-phosphate in the solvent of choice, which should be purged with an inert gas. Clindamycin 3-phosphate is slightly soluble (0.1-1 mg/ml) in DMSO and a 1:1 solution of acetonitrile:water.

Description

Clindamycin 3-phosphate is a microbial biotransformation product and prodrug form of the lincosamide antibiotic clindamycin (Item Nos. 15006 | 35581).^{1,2} It is formed from clindamycin by *S. coelicolor*, resulting in inactivation of the antibiotic activity of clindamycin.² Clindamycin 3-phosphate increases survival in a mouse model of *S. aureus* infection.¹

References

1. Brodasky, T.F. and Lewis, C. An *in vivo-in vitro* comparison of the 2- and 3-phosphate esters of clindamycin. *J. Antibiot. (Tokyo)* **25(4)**, 230-238 (1972).
2. Coats, J.H. and Argoudelis, A.D. Microbial transformation of antibiotics: Phosphorylation of clindamycin by *Streptomyces coelicolor* Müller. *J. Bacteriol.* **108(1)**, 459-464 (1971).

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, 08/17/2026

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