

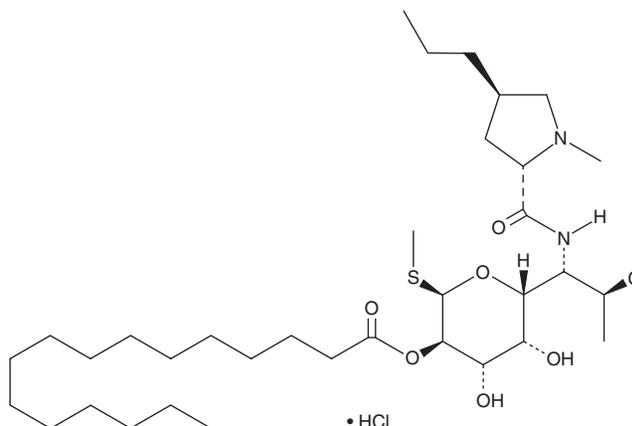
# PRODUCT INFORMATION



## Clindamycin Palmitate (hydrochloride)

Item No. 29102

**CAS Registry No.:** 25507-04-4  
**Formal Name:** 2-hexadecanoate methyl 7-chloro-6,7,8-trideoxy-6-[[[(2S,4R)-1-methyl-4-propyl-2-pyrrolidiny]carbonyl]amino]-1-thio-L-threo- $\alpha$ -D-galacto-octopyranoside, monohydrochloride  
**Synonym:** Clindamycin 2-Palmitate  
**MF:** C<sub>34</sub>H<sub>63</sub>ClN<sub>2</sub>O<sub>6</sub>S • HCl  
**FW:** 699.9  
**Purity:** ≥95%  
**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

Clindamycin palmitate (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the clindamycin palmitate (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Clindamycin palmitate (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of clindamycin palmitate (hydrochloride) in ethanol is approximately 30 mg/ml and approximately 25 mg/ml in DMSO and DMF.

### Description

Clindamycin palmitate is a water-soluble prodrug form of the antibiotic clindamycin (Item No. 15006).<sup>1</sup> It undergoes rapid hydrolysis *in vivo* to release active clindamycin. Formulations containing clindamycin palmitate have been used in the treatment of various bacterial infections.

### Reference

1. Bharathi, C., Jayaram, P., Sunder Raj, J., *et al.* Identification, isolation and characterization of impurities of clindamycin palmitate hydrochloride. *J. Pharm. Biomed. Anal.* **48(4)**, 1211-1218 (2008).

#### 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, 12/05/2022

#### 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