

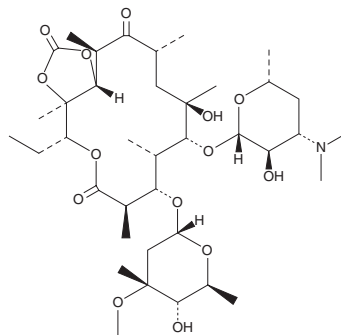
PRODUCT INFORMATION



Davercin

Item No. 21348

CAS Registry No.: 55224-05-0
Formal Name: cyclic 11,12-carbonate erythromycin
Synonym: Erythromycin cyclocarbonate
MF: C₃₈H₆₅NO₁₄
FW: 759.9
Purity: ≥98%
UV/Vis.: λ_{max}: 205 nm
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

Davercin is supplied as a solid. A stock solution may be made by dissolving the davercin in the solvent of choice, which should be purged with an inert gas. Davercin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of davercin in ethanol is approximately 30 mg/ml and approximately 15 mg/ml in DMSO and DMF.

Davercin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, davercin should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Davercin has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Davercin is a macrolide antibiotic that inhibits bacterial protein synthesis by targeting the 50S ribosomal subunit, blocking the progression of nascent polypeptide chains.¹ It has broad-spectrum activity against various Gram-positive and Gram-negative bacteria (MICs = 0.02-50 µg/ml). Davercin is non-toxic to mice and rats (LD₅₀s = 4.05 and 5.9 g/kg, respectively). Formulations containing davercin have been used to treat a variety of urogenital bacterial infections.

Reference

1. Bojarska-Dahlig, H., Slawinski, W., Roślik-Kamińska, D., *et al.* L-aspartate of erythromycin A cyclic 11,12-carbonate, a new semisynthetic erythromycin derivative. *J. Antibiot. (Tokyo)* **29(9)**, 907-914 (1976).

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, 11/14/2022

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