

PRODUCT INFORMATION



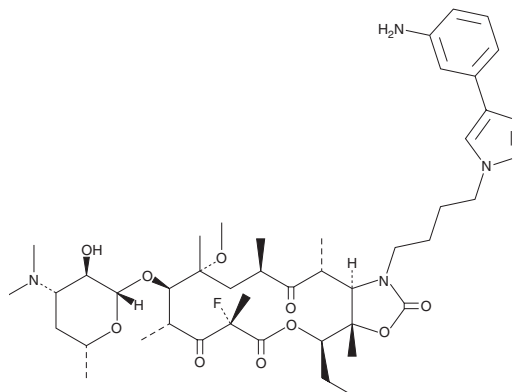
Solithromycin

Item No. 21681

CAS Registry No.: 760981-83-7

Formal Name: (3aS,4R,7S,9R,10R,11R,13R,15R,15aR)-1-[4-[4-(3-aminophenyl)-1H-1,2,3-triazol-1-yl]butyl]-4-ethyl-7-fluorooctahydro-11-methoxy-3a,7,9,11,13,15-hexamethyl-10-[[3,4,6-trideoxy-3-(dimethylamino)-β-D-xyllo-hexopyranosyl]oxy]-2H-oxacyclotetradecino[4,3-d]oxazole-2,6,8,14(1H,7H,9H)-tetrone

MF: $C_{43}H_{65}FN_6O_{10}$
FW: 845.0
Purity: ≥98%
UV/Vis.: λ_{max} : 233, 311 nm
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

Solithromycin is supplied as a crystalline solid. A stock solution may be made by dissolving the solithromycin in the solvent of choice, which should be purged with an inert gas. Solithromycin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of solithromycin in these solvents is approximately 2.5 and 5 mg/ml, respectively.

Description

Solithromycin is a macrolide antibiotic with activity against Gram-positive and Gram-negative bacteria, including *C. trachomatis* and *C. pneumoniae* (MICs = 0.125-0.5 and 0.25-1 µg/ml, respectively).¹ It also inhibits growth of multidrug resistant *P. falciparum* *in vitro* (IC₅₀s = 0.5 nM for both Dd2 and 7G8 strains).² Solithromycin (100 mg/kg) administered in a four-dose series eradicates *P. berghei* infection in mice.

References

1. Roblin, P.M., Kohlhoff, S.A., Parker, C., *et al.* *In vitro* activity of CEM-101, a new fluoroketolide antibiotic, against *Chlamydia trachomatis* and *Chlamydia (Chlamydophila) pneumoniae*. *Antimicrob. Agents Chemother.* **54**(3), 1358-1359 (2010).
2. Wittlin, S., Eklund, E.H., Craft, J.C., *et al.* *In vitro* and *in vivo* activity of solithromycin (CEM-101) against *Plasmodium* species. *Antimicrob. Agents Chemother.* **56**(2), 703-707 (2012).

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, 11/29/2022

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