

PRODUCT INFORMATION



Tulathromycin A

Item No. 19202

CAS Registry No.: 217500-96-4

Formal Name: (2R,3S,4R,5R,8R,10R,11R,12S,13S,14R)-13-[[[2,6-dideoxy-3-C-methyl-3-O-methyl-4-C-[(propylamino)methyl]- α -L-ribo-hexopyranosyl]oxy]-2-ethyl-3,4,10-trihydroxy-3,5,8,10,12,14-hexamethyl-11-[[[3,4,6-trideoxy-3-(dimethylamino)- β -D-xylo-hexopyranosyl]oxy]-1-oxa-6-azacyclopentadecan-15-one

Synonym:

MF: $C_{41}H_{79}N_3O_{12}$

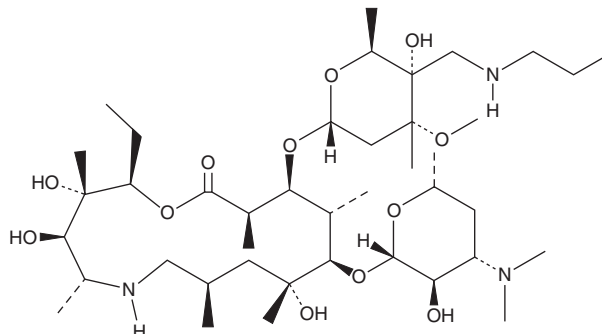
FW: 806.1

Purity: $\geq 95\%$

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

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

Tulathromycin A is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, tulathromycin A should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Tulathromycin A 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

Tulathromycin A is a macrolide antibiotic that inhibits bacterial protein synthesis ($IC_{50} = 0.26 \mu\text{M}$) by targeting the 50S ribosomal subunit, blocking the progression of nascent polypeptide chains.¹ It is effective against *M. haemolytica* and *P. multocida* (MICs = 2 and 0.5 $\mu\text{g/ml}$, respectively), which cause respiratory tract infections in cattle and swine.¹

Reference

1. Andersen, N.M., Poehlsgaard, J., Warrass, R., *et al.* Inhibition of protein synthesis on the ribosome by tildipirosin compared with other veterinary macrolides. *Antimicrob. Agents Chemother.* **56**(11), 6033-6036 (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, 11/23/2022

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