

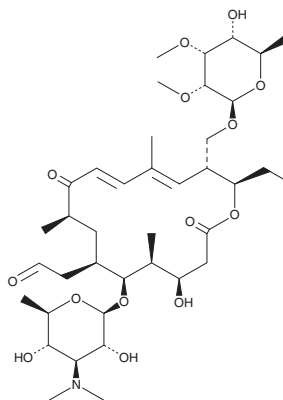
PRODUCT INFORMATION



Desmycosin

Item No. 41972

CAS Registry No.: 11032-98-7
Formal Name: 4^A-O-de(2,6-dideoxy-3-C-methyl- α -L-ribo-hexopyranosyl)-tylosin
Synonyms: Demycarosyltylosin, Oxacyclohexadecane, Tylosin B
MF: C₃₉H₆₅NO₁₄
FW: 771.9
Purity: $\geq 90\%$
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

Desmycosin is supplied as a solid. A stock solution may be made by dissolving the desmycosin in the solvent of choice, which should be purged with an inert gas. Desmycosin is soluble (≥ 10 mg/ml) in ethanol. Desmycosin is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Desmycosin is a macrolide antibiotic.¹ It is also an acidic degradation product of tylosin (Item No. 22599).² Desmycosin is active against *H. influenzae*, *S. aureus*, *S. pneumoniae*, and *S. pyogenes* (MICs = 4, 1, <0.125 , and <0.125 $\mu\text{g/ml}$, respectively).¹

References

1. Iveković, D., Lopotar, N., Brajsa, K., *et al.* Electrochemical reduction of desmycosin, structure investigation and antibacterial evaluation of the resulting products. *Eur. J. Pharm. Sci.* **18**(5), 323-328 (2003).
2. Thompson, T.S., van den Heever, J.P., and Komarnicki, J.A.F. Tylosin A and desmycosin in honey by salting-out assisted liquid-liquid extraction and aqueous normal phase ultraperformance liquid chromatography-tandem mass spectrometry. *Anal. Bioanal. Chem.* **411**(24), 6509-6518 (2019).

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, 10/07/2025

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