

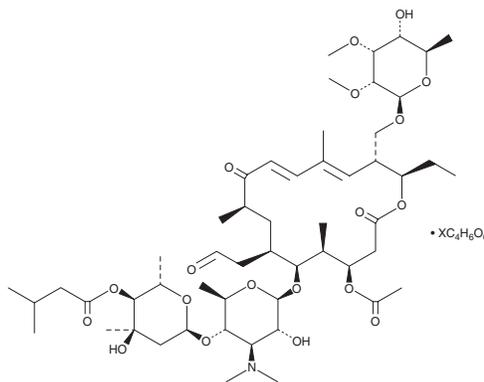
# PRODUCT INFORMATION



## Tylvalosin (tartrate)

Item No. 33849

**CAS Registry No.:** 63428-13-7  
**Formal Name:** 3-acetate 4<sup>B</sup>-(3-methylbutanoate) tylosin, 2R,3R-dihydroxybutanedioate  
**Synonym:** Acetylisovaleryltylosin tartrate  
**MF:** C<sub>53</sub>H<sub>87</sub>NO<sub>19</sub> • XC<sub>4</sub>H<sub>6</sub>O<sub>6</sub>  
**FW:** 1,042.3  
**Purity:** ≥80%  
**UV/Vis.:** λ<sub>max</sub>: 290 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

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

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of tylvalosin (tartrate) can be prepared by directly dissolving the solid in aqueous buffers. Tylvalosin (tartrate) is slightly soluble in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

Tylvalosin is a macrolide antibiotic.<sup>1</sup> It is active against *S. aureus*, *E. coli*, and *P. multocida* (MICs = 2, 128, and 128 µg/ml, respectively). Tylvalosin (5 and 10 µg/ml) reduces LPS-induced production of pro-inflammatory cytokines, prostaglandin E<sub>2</sub> (PGE<sub>2</sub>; Item No. 14010), and nitric oxide in RAW 264.7 cells.<sup>2</sup> It reduces the severity of lung lesions in pigs infected with *M. hyopneumoniae* and *P. multocida* when administered at a dose of 10 mg/kg per day in the drinking water.<sup>3</sup> Formulations containing tylvalosin have been used in the treatment and metaphylaxis of enzootic pneumonia caused by *M. hyopneumoniae* in pigs.

### References

1. Yang, B., Lei, Z., Zhao, Y., et al. Combination susceptibility testing of common antimicrobials *in vitro* and the effects of sub-MIC of antimicrobials on *Staphylococcus aureus* biofilm formation. *Front. Microbiol.* **8**, 2125 (2017).
2. Zhao, Z., Tang, X., Zhao, X., et al. Tylvalosin exhibits anti-inflammatory property and attenuates acute lung injury in different models possibly through suppression of NF-κB activation. *Biochem. Pharmacol.* **90(1)**, 73-87 (2014).
3. Lopez Rodriguez, A., Berge, A.C., Ramage, C., et al. Evaluation of the clinical efficacy of a water soluble formulation of tylvalosin in the control of enzootic pneumonia associated with *Mycoplasma hyopneumoniae* and *Pasteurella multocida* in pigs. *Porcine Health Manag.* **6(1)**, 39 (2020).

#### 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, 11/07/2022

#### 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