

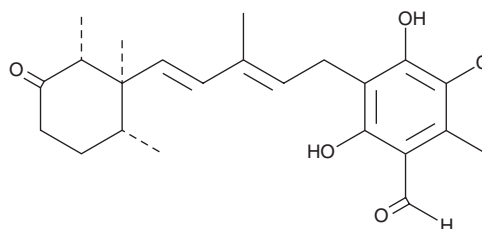
# PRODUCT INFORMATION



## Ascochlorin

Item No. 19437

**CAS Registry No.:** 26166-39-2  
**Formal Name:** 3-chloro-4,6-dihydroxy-2-methyl-5-[(2E,4E)-3-methyl-5-[(1R,2R,6R)-1,2,6-trimethyl-3-oxocyclohexyl]-2,4-pentadien-1-yl]-benzaldehyde  
**Synonyms:** Antibiotic LL-Z1272y, Ilicicolin D, NSC 287492  
**MF:** C<sub>23</sub>H<sub>29</sub>ClO<sub>4</sub>  
**FW:** 404.9  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Fungus/*Acremonium* sp.



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Ascochlorin is supplied as a solid. A stock solution may be made by dissolving the ascochlorin in the solvent of choice, which should be purged with an inert gas. Ascochlorin is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

### Description

Ascochlorin is an isoprenoid antibiotic and antiviral that has diverse effects on mammalian cells.<sup>1,2</sup> It suppresses PMA-induced invasion in renal carcinoma cells (IC<sub>50</sub> = ~10 μM) by inhibiting the expression of matrix metalloproteinase-9.<sup>3</sup> At 2 μM, ascochlorin profoundly increases the expression of p53 by increasing protein stability in cancer cells, and, at 50 μM, it inhibits signaling through STAT3.<sup>4,5</sup> Ascochlorin also binds and inhibits the mitochondrial cytochrome bc<sub>1</sub> complex, blocking reduction of cytochrome *b* by ubiquinone.<sup>6</sup>

### References

1. Tamura, G., Suzuki, S., Takatsuki, A., et al. *J. Antibiot. (Tokyo)* **21(9)**, 539-544 (1968).
2. Takatsuki, A., Tamura, G., and Arima, K. *App. Microbiol.* **17(6)**, 825-829 (1969).
3. Hong, S., Park, K.-K., Magae, J., et al. *J. Biol. Chem.* **280(26)**, 25202-20209 (2005).
4. Jeong, J.-H., Nakajima, H., Magae, J., et al. *Int. J. Cancer* **124(12)**, 2797-2803 (2009).
5. Dai, X., Ahn, K.S., Kim, C., et al. *Mol. Oncol.* **9(4)**, 818-833 (2015).
6. Berry, E.A., Huang, L., Lee, D.-W., et al. *Biochim. Biophys. Acta.* **1797(3)**, 360-370 (2016).

#### 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, 03/13/2026

#### 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