

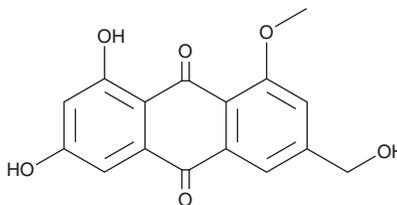
PRODUCT INFORMATION



Carviolin

Item No. 29634

CAS Registry No.: 478-35-3
Formal Name: 1,3-dihydroxy-6-(hydroxymethyl)-8-methoxy-9,10-anthracenedione
MF: C₁₆H₁₂O₆
FW: 300.3
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Penicillium* sp.



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

Laboratory Procedures

Carviolin is supplied as a solid. A stock solution may be made by dissolving the carviolin in the solvent of choice, which should be purged with an inert gas. Carviolin is soluble in ethanol, methanol, dichloromethane, and DMSO.

Description

Carviolin is an anthraquinone fungal metabolite that has been found in *Z. longicaudata* with immunosuppressive and antitrypanosomal activities.^{1,2} It inhibits LPS- or concanavalin A-induced proliferation of mouse splenocytes (IC₅₀s = 4 and 4.5 μg/ml, respectively).¹ Carviolin is active against *T. b. brucei* (MIC = 41.66 μM).²

References

1. Fujimoto, H., Nakamura, E., Okuyama, E., *et al.* Six immunosuppressive features from an ascomycete, *Zopfiella longicaudata*, found in a screening study monitored by immunomodulatory activity. *Chem. Pharm. Bull. (Tokyo)* **52(8)**, 1005-1008 (2004).
2. Aly, A.H., Debbab, A., Clements, C.M., *et al.* NF kappa B inhibitors and antitrypanosomal metabolites from endophytic fungus *Penicillium* sp. isolated from *Limonium tubiflorum*. *Bioorg. Med. Chem.* **19(1)**, 414-421 (2011).

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, 12/08/2022

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