

PRODUCT INFORMATION



Remisporine B

Item No. 31669

CAS Registry No.: 571194-06-4
Formal Name: (5bS,7S,7aR,7bR,14bS)-7,7a,7b,8,14,14b-hexahydro-4,7,13-trihydroxy-2,11-dimethyl-5,14-dioxo-furo[2',3',4':3,4]indeno[1,2-b:6,7-b']bis[1]benzopyran-5b,7(5H)-dicarboxylic acid, dimethyl ester

MF: C₃₀H₂₄O₁₂

FW: 576.5

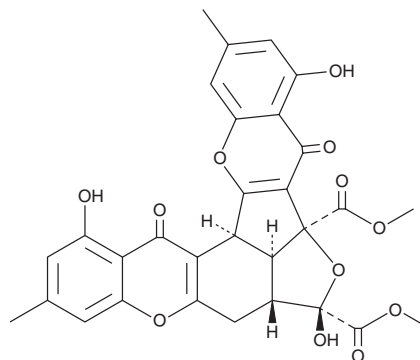
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Remispora maritima*



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

Laboratory Procedures

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

Description

Remisporine B is a polyketide and derivative of remisporine A that has been found in *Penicillium* and has immunosuppressant activity.^{1,2} It inhibits LPS- or concanavalin A-induced proliferation of isolated mouse splenic lymphocytes (IC₅₀s = 30.1 and 32.4 μg/ml, respectively).¹

References

1. Liu, H., Chen, S., Liu, W., *et al.* Polyketides with immunosuppressive activities from mangrove endophytic fungus *Penicillium* sp. ZJ-SY₂. *Mar. Drugs* **14**(12), 217 (2016).
2. Kong, F., and Carter, G.T. Remisporine B, a novel dimeric chromenone derived from spontaneous Diels-Alder reaction of remisporine A. *Tetrahedron Lett.* **44**(15), 3119-3122 (2003).

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, 08/11/2023

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