

PRODUCT INFORMATION



Echinulin

Item No. 22711

CAS Registry No.: 1859-87-6
Formal Name: (3S,6S)-3-[[2-(1,1-dimethyl-2-propen-1-yl)-5,7-bis(3-methyl-2-buten-1-yl)-1H-indol-3-yl]methyl]-6-methyl-2,5-piperazinedione

MF: C₂₉H₃₉N₃O₂

FW: 461.6

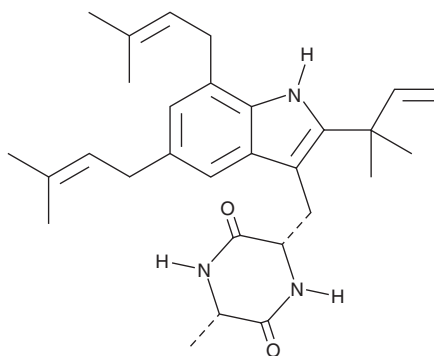
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Aspergillus* sp.



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

Laboratory Procedures

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

Description

Echinulin is a diketopiperazine metabolite found in fungi that, when administered to rabbits, at a dose of 10 mg/kg, leads to liver and lung damage.¹ It has no activity at the δ , κ , or μ -opioid receptors or the cannabinoid CB₁ and CB₂ receptors.²

References

1. Ali, M., Mohammed, N., Alnaqeeb, M.A., *et al.* Toxicity of echinulin from *Aspergillus chevalieri* in rabbits. *Toxicol. Lett.* **48(3)**, 235-241 (1989).
2. Gao, J., Leon, F., Radwan, M.M., *et al.* Benzyl derivatives with *in vitro* binding affinity for human opioid and cannabinoid receptors from the fungus *Eurotium repens*. *J. Nat. Prod.* **74(7)**, 1636-1639 (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, 02/19/2024

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