

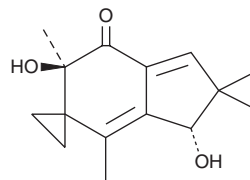
PRODUCT INFORMATION



Illudin M

Item No. 22062

CAS Registry No.: 1146-04-9
Formal Name: (3'S,6'R)-2',3'-dihydro-3',6'-dihydroxy-2',2',4',6'-tetramethyl-spiro[cyclopropane-1,5'-[5H]inden]-7'(6'H)-one
Synonyms: DR-15977, (-)-Illudin M, NSC 400978, NSC 626370
MF: C₁₅H₂₀O₃
FW: 248.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

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

Description

Illudin M is a cytotoxic sesquiterpene from the fungus *O. illudens* that alkylates DNA.¹ It is cytotoxic to HL-60 human leukemia cells at 6-100 nM.² Illudin M was used as a base structure for the development of potential anticancer agents with less toxicity and improved efficacy.^{1,3} It is closely related to illudin S (Item No. 17451), which has been the focus of additional studies.

References

1. Jacobson, J.M., Feinman, L., Liebes, L., *et al.* Pharmacokinetics, safety, and antiviral effects of hypericin, a derivative of St. John's wort plant, in patients with chronic hepatitis C virus infection. *Antimicrob. Agents and Chemother.* **45(2)**, 517-524 (2001).
2. Kelner, M.J., McMorris, T.C., Beck, W.T., *et al.* Preclinical evaluation of illudins as anticancer agents. *Cancer Research* **47(12)**, 3186-3189 (1987).
3. Schobert, R., Knauer, S., Seibt, S., *et al.* Anticancer active illudins: Recent developments of a potent alkylating compound class. *Current Medicinal Chemistry* **18**, 790-807 (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, 05/25/2017

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