

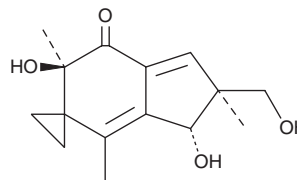
# PRODUCT INFORMATION



## Illudin S

Item No. 17451

**CAS Registry No.:** 1149-99-1  
**Formal Name:** (2'S,3'R,6'R)-2',3'-dihydro-3',6'-dihydroxy-2'-(hydroxymethyl)-2',4',6'-trimethyl-spiro[cyclopropane-1,5'-[5H]inden]-7'(6'H)-one  
**Synonyms:** Lampterol, NSC 400979, NSC 626369  
**MF:** C<sub>15</sub>H<sub>20</sub>O<sub>4</sub>  
**FW:** 264.3  
**Purity:** ≥98%  
**Supplied as:** A light yellow solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Fungus/*Pleurotus lampas*



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

### Laboratory Procedures

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

Illudin S is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

Illudins are fungal sesquiterpenes that, through their unique DNA alkylating actions, have anticancer potential.<sup>1,2</sup> Illudin S is a cytotoxic illudin that is converted, intracellularly, to metabolites that cause a complete block of cell cycling at the G<sub>1</sub>-S phase interface, particularly in myeloid and T-lymphocyte leukemia cells (IC<sub>50</sub> = 6-11 nM).<sup>1</sup> T-lymphocyte leukemia CEM cells that are resistant to doxorubicin (Item No. 15007), epipodophyllotoxins, and 1-β-D-arabinofuranosylcytosine display only 2-fold increased resistance to illudin S.<sup>1</sup> Illudin S metabolites induce DNA damage that is not repaired by the processes that counter conventional DNA alkylating agents.<sup>2-4</sup>

### References

1. Kelner, M.J., McMorris, T.C., Beck, W.T., *et al. Cancer Res.* **47(12)**, 3186-3189 (1987).
2. Schobert, R., Knauer, S., Seibt, S., *et al. Curr. Med. Chem.* **18**, 790-807 (2011).
3. Kelner, M.J., McMorris, T.C., Estes, L., *et al. Biochem. Pharmacol.* **48**, 403-409 (1994).
4. Jaspers, N.G.J., Raams, A., Kelner, M.J., *et al. DNA Repair (Amst)* **1(12)**, 1027-1038 (2002).

#### 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, 08/08/2023

#### 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