

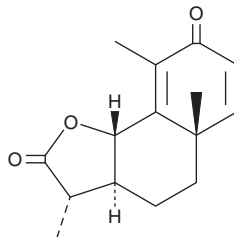
PRODUCT INFORMATION



Santonin

Item No. 27460

CAS Registry No.: 481-06-1
Formal Name: (3S,3aS,5aS,9bS)-3a,5,5a,9b-tetrahydro-3,5a,9-trimethyl-naphtho[1,2-b]furan-2,8(3H,4H)-dione
Synonyms: NSC 4900, (-)-Santonin, α -Santonin
MF: C₁₅H₁₈O₃
FW: 246.3
Purity: \geq 98%
UV/Vis.: λ_{max} : 242 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/Unidentified sp.



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

Laboratory Procedures

Santonin is supplied as a crystalline solid. A stock solution may be made by dissolving the santonin in the solvent of choice, which should be purged with an inert gas. Santonin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of santonin in these solvents is approximately 5 mg/ml. Santonin is also slightly soluble in ethanol.

Santonin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, santonin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Santonin has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Santonin is a sesquiterpenoid lactone that has been found in *Artemisia*.¹ It has been used as a building block in the synthesis of terpenoids. Santonin is an anthelmintic that induces paralysis of *A. lumbricoides*.²

References

1. Banerjee, A.K., Vera, W.J., and Gonzalez, N.C. Synthesis of terpenoid compounds from α -santonin. *Tetrahedron* **49**(122), 4761-4788 (1993).
2. Hermoso, R. and Monteoliva, M. Survival study of parasitic helminths. III. Effects of antihelminthic and other drugs on survival of *Ascaris lumbricoides*. *Rev. Iber. Parasitol.* **31**(4), 567-585 (1970).

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, 09/15/2023

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