

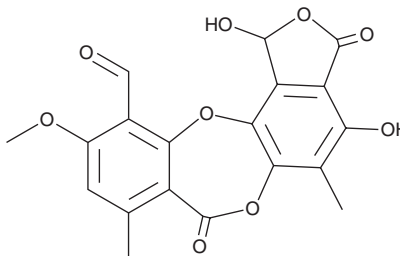
PRODUCT INFORMATION



Stictic Acid

Item No. 25017

CAS Registry No.: 549-06-4
Formal Name: 1,3-dihydro-1,4-dihydroxy-10-methoxy-5,8-dimethyl-3,7-dioxo-7H-isobenzofuro[4,5-b][1,4]benzodioxepin-11-carboxaldehyde
Synonym: NSC 87511
MF: C₁₉H₁₄O₉
FW: 386.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

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

Description

Stictic acid is a depsidone found in lichen.¹ It preferentially inhibits growth of HT-29 human colon adenocarcinoma cells over non-malignant cells (IC₅₀s = 29.29 and 2,478.4 µg/ml, respectively).²

References

1. Rao, P.S., Sarma, K.G., and Seshadri, T.R. The ultraviolet and infrared spectra of some lichen depsides and depsidones. *Proc. Ind. Acad. Sci.* **66**(1), 1-14 (1967).
2. Pejin, B., Iodice, C., Bogdanović, G., *et al.* Stictic acid inhibits cell growth of human colon adenocarcinoma HT-29 cells. *Arabian J. Chem.* **10**(Suppl. 1), S1240-S1242 (2017).

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, 12/19/2022

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