

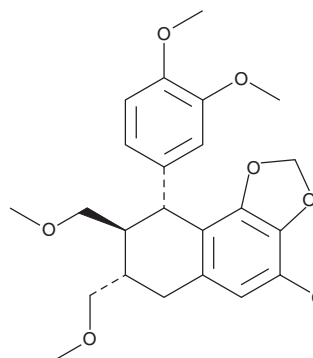
PRODUCT INFORMATION



Hypophyllanthin

Item No. 11715

CAS Registry No.: 33676-00-5
Formal Name: (7S,8S,9R)-9-(3,4-dimethoxyphenyl)-6,7,8,9-tetrahydro-4-methoxy-7,8-bis(methoxymethyl)-naphtho[1,2-d]-1,3-dioxole
Synonym: NSC 619044
MF: C₂₄H₃₀O₇
FW: 430.5
Purity: ≥98%
Supplied as: A crystalline 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

Hypophyllanthin is supplied as a crystalline solid. A stock solution may be made by dissolving the hypophyllanthin in the solvent of choice, which should be purged with an inert gas. Hypophyllanthin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of hypophyllanthin in these solvents is approximately 0.2, 10, and 30 mg/ml, respectively.

Hypophyllanthin 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

Hypophyllanthin is an active principle of *P. amarus*, a plant studied for its potential to treat hepatotoxicity.¹ In Caco-2 cells 100 μM hypophyllanthin has been shown to reversibly inhibit P-glycoprotein (also known as multidrug resistance protein 1 (MDR1)) function without effect on MDR2 activity.²

References

1. Murali, B., Amit, A., Anand, M.S., *et al.* An improved HPLC method for estimation of phyllanthin and hypophyllanthin in *Phyllanthus amarus*. *Journal of Natural Remedies* **1(1)**, 55-59 (2001).
2. Sukhaphirom, N., Vardhanabhuti, N., Chirdchupunsee, H., *et al.* Phyllanthin and hypophyllanthin inhibit function of P-gp but not MRP2 in Caco-2 cells. *J. Pharm. Pharmacol.* **65**, 292-299 (2012).

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, 10/24/2022

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