

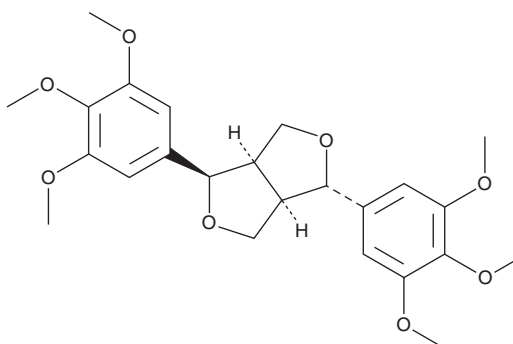
PRODUCT INFORMATION



Lirioresinol A dimethyl ether

Item No. 44940

CAS Registry No.: 24192-64-1
Formal Name: (1R,3aR,4S,6aR)-tetrahydro-1,4-bis(3,4,5-trimethoxyphenyl)-1H,3H-furo[3,4-c]furan
Synonyms: (+)-Epiyangambin, *epi*-Yangambin, NSC 172767
MF: C₂₄H₃₀O₈
FW: 446.5
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Magnolia officinalis*



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

Laboratory Procedures

Lirioresinol A dimethyl ether is supplied as a solid. A stock solution may be made by dissolving the lirioresinol A dimethyl ether in the solvent of choice, which should be purged with an inert gas. Lirioresinol A dimethyl ether is soluble in methanol and DMSO.

Description

Lirioresinol A dimethyl ether is a lignan that has been found in *H. nymphaeifolia* and has diverse biological activities.¹⁻³ It inhibits the aggregation of isolated rabbit platelets induced by platelet-activating factor (PAF), collagen, or arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) when used at a concentration of 100 μM.¹ Lirioresinol A dimethyl ether inhibits the proliferation of SK-MEL-147 melanoma cells (IC₅₀ = 13.09 μg/ml) and induces cell cycle arrest at the G₂/M phase in the same cells.² It is active against *L. amazonensis* and *L. braziliensis* intracellular amastigotes (IC₅₀s = 22.6 and 74.4 μM, respectively).³

References

1. Chen, J.-J., Chang, Y.-L., Teng, C.-M., *et al.* Anti-platelet aggregation alkaloids and lignans from *Hernandia nymphaeifolia*. *Planta Med.* **66**(3), 251-256 (1999).
2. Garcia, D.A., Pressete, C.G., Miranda, R., *et al.* Biological and metabolomics-guided isolation of tetrahydrofurofuran lignan from *Croton* spp. with antiproliferative activity against human melanoma cell line. *Fitoterapia* **177**, 106070 (2024).
3. Rebouças-Silva, J., Santos, G.F., Filho, J.M.B., *et al.* *In vitro* leishmanicidal effect of Yangambin and Epi-yangambin lignans isolated from *Ocotea fasciculata* (Nees) Mez. *Front Cell. Infect. Microbiol.* **12**, 1045732 (2022).

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, 01/27/2026

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