

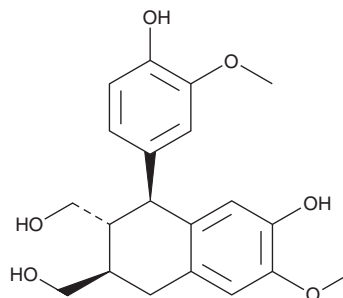
PRODUCT INFORMATION



Isolariciresinol

Item No. 36032

CAS Registry No.: 548-29-8
Formal Name: (1S,2R,3R)-1,2,3,4-tetrahydro-7-hydroxy-1-(4-hydroxy-3-methoxyphenyl)-6-methoxy-2,3-naphthalenedimethanol
Synonyms: (+)-Cyclolariciresinol, (+)-Isolariciresinol, (7'S,8'R,8R)-Isolariciresinol
MF: C₂₀H₂₄O₆
FW: 360.4
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Phyllanthus niruri*



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

Laboratory Procedures

Isolariciresinol is supplied as a solid. A stock solution may be made by dissolving the isolariciresinol in the solvent of choice, which should be purged with an inert gas. Isolariciresinol is soluble in acetone, chloroform, dichloromethane, DMSO, and ethyl acetate.

Description

Isolariciresinol is a lignan that has been found in *Cornus officinalis* and has diverse biological activities.^{1,2} It scavenges ABTS (Item No. 27317) radicals in a cell-free assay (IC₅₀ = 19.4 μM) and inhibits LPS-induced nitric oxide (NO) production in RAW 264.7 macrophages (IC₅₀ = 84.45 μM).¹ Isolariciresinol (20 μM) inhibits amyloid-β (1-42) (Aβ42; Item No. 20574) self-aggregation in a cell-free assay.²

References

1. Peng, Z., Wang, Y., He, J., et al. Chemical constituents and their antioxidant and anti-inflammatory activities from edible *Cornus officinalis* fruits. *Eur. Food Res. Technol.* **248**, 1003-1010 (2022).
2. Huang, X.-X., Xu, Y., Bai, M., et al. Lignans from the seeds of Chinese hawthorn (*Crataegus pinnatifida* var. major N.E.Br.) against β-amyloid aggregation. *Nat. Prod. Res.* **32(14)**, 1706-1713 (2018).

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, 12/18/2024

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