

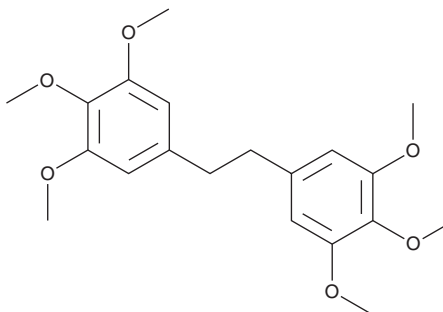
PRODUCT INFORMATION



Brittonin A

Item No. 44901

CAS Registry No.: 33284-74-1
Formal Name: 1,1'-(1,2-ethanediyl)bis[3,4,5-trimethoxy-benzene]
Synonyms: 1,2-Bis(3,4,5-trimethoxyphenyl)ethane, NSC 600172
MF: C₂₀H₂₆O₆
FW: 362.4
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

Brittonin A is supplied as a solid. A stock solution may be made by dissolving the brittonin A in the solvent of choice, which should be purged with an inert gas. Brittonin A is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Brittonin A is a methoxylated bibenzyl that has been found in the liverwort *F. inouei* and has anticancer activity.¹ It is cytotoxic to KB and vincristine-resistant KB/VCR cells (IC₅₀s = 42.1 and 33.7 μM, respectively), as well as K562 and adriamycin-resistant K562/A02 cells (IC₅₀s = 49.6 and 30.9 μM, respectively). Brittonin A (5 μM) increases vincristine-induced cell death in KB/VCR cells and adriamycin-induced cell death in K562/A02 cells.

Reference

1. Guo, D.X., Xiang, F., Wang, X.N., *et al.* Labdane diterpenoids and highly methoxylated bibenzyls from the liverwort *Frullania inouei*. *Phytochemistry* **71**(13), 1573-1578 (2010).

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, 11/20/2025

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