

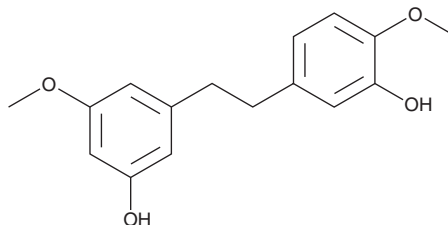
PRODUCT INFORMATION



Gigantol isomer-1

Item No. 46245

CAS Registry No.: 67884-30-4
Formal Name: 5-[2-(3-hydroxy-5-methoxyphenyl)ethyl]-2-methoxy-phenol
MF: $C_{16}H_{18}O_4$
FW: 274.3
Purity: $\geq 98\%$
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

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

Description

Gigantol isomer-1 is a bibenzyl derivative and gigantol isomer that has been found in *E. palustris*.¹ The levels of gigantol isomer-1 increase in *E. palustris* rhizomes after wounding.

Reference

1. Gehlert, R. and Kindl, H. Induced formation of dihydrophenanthrenes and bibenzyl synthase upon destruction of orchid mycorrhiza. *Phytochemistry* **30**(2), 457-460 (1991).

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, 07/29/2026

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