

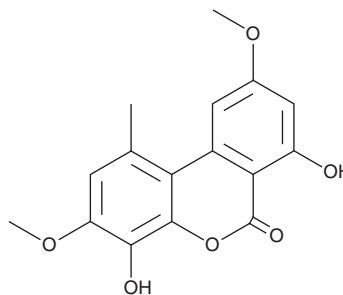
PRODUCT INFORMATION



Graphislactone A

Item No. 23643

CAS Registry No.: 52179-44-9
Formal Name: 4,7-dihydroxy-3,9-dimethoxy-1-methyl-6H-dibenzo[b,d]pyran-6-one
MF: $C_{16}H_{14}O_6$
FW: 302.3
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Fungus/*Cephalosporium* sp.



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

Laboratory Procedures

Graphislactone A is supplied as a solid. A stock solution may be made by dissolving the graphislactone A in the solvent of choice, which should be purged with an inert gas. Graphislactone A is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

Graphislactone A is an antioxidant produced by *Cephalosporium* IFB-E001 and *M. olivacea*.^{1,2} It exhibits dose-dependent scavenging activity against DPPH radicals and metal ion-dependent hydroxyl radicals *in vitro*.¹ Graphislactone A reduces peroxidation of human LDL and slows peroxidation of linoleic acid (Item Nos. 90150 | 21909) by 4.4-fold. It also inhibits acetylcholinesterase (AChE) with an IC_{50} value of $27 \mu\text{M}$.²

References

1. Song, Y.C., Huang, W.Y., Sun, C., *et al.* Characterization of graphislactone A as the antioxidant and free radical-scavenging substance from the culture of *Cephalosporium* sp. IFB-E001, an endophytic fungus in *Trachelospermum jasminoides*. *Biol. Pharm. Bull.* **28**(3), 506-509 (2005).
2. Hormazabal, E., Schmeda-Hirschmann, G., Astudillo, L., *et al.* Metabolites from *Microsphaeropsis olivacea*, an endophytic fungus of *Pilgerodendron uviferum*. *Z. Naturforsch. C.* **60**(1-2), 11-21 (2005).

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, 08/14/2024

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