

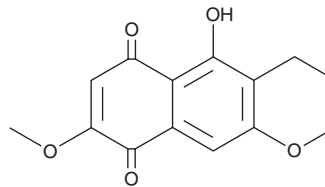
PRODUCT INFORMATION



6-Ethyl-2,7-dimethoxyjuglone

Item No. 28049

CAS Registry No.: 15254-86-1
Formal Name: 6-ethyl-5-hydroxy-2,7-dimethoxy-1,4-naphthalenedione
Synonym: 2,7-Dimethoxy-6-ethyljuglone
MF: C₁₄H₁₄O₅
FW: 262.3
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/Unidentified sp.



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

Laboratory Procedures

6-Ethyl-2,7-dimethoxyjuglone is supplied as a solid. A stock solution may be made by dissolving the 6-ethyl-2,7-dimethoxyjuglone in the solvent of choice, which should be purged with an inert gas. 6-Ethyl-2,7-dimethoxyjuglone is soluble in organic solvents such as ethanol, methanol, and DMSO.

Description

6-Ethyl-2,7-dimethoxyjuglone is a naphthoquinone fungal metabolite originally isolated from *H. toruloidea* and has antimicrobial activity.^{1,2} It is active against the bacteria *B. cereus* and *S. aureus* and the plant pathogenic fungi *Alternaria*, *Sclerotium*, and *Fusarium* in disk assays when used at a concentration of 50 µg per disk.²

References

1. van Eijk, G.W. and Roeymans, H.J. Naphthoquinone derivatives from the fungus *Hendersonula toruloidea*. *Experientia* **34(10)**, 1257-1258 (1978).
2. Sun, R., Gao, Y.-X., Shen, K.-Z., *et al.* Antimicrobial metabolites from the aquatic fungus *Delitschia corticola*. *Phytochem. Lett.* **4(2)**, 101-105 (2011).

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, 08/10/2023

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