

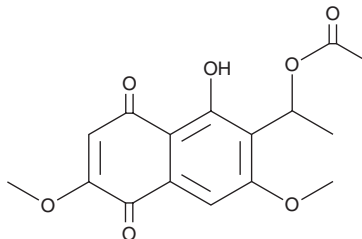
PRODUCT INFORMATION



2,7-Dimethoxy-6-(1-acetoxyethyl)juglone

Item No. 27267

CAS Registry No.: 68594-15-0
Formal Name: (-)-6-[1-(acetyloxy)ethyl]-5-hydroxy-2,7-dimethoxy-1,4-naphthalenedione
MF: C₁₆H₁₆O₇
FW: 320.3
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Zopfia* sp.



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

Laboratory Procedures

2,7-Dimethoxy-6-(1-acetoxyethyl)juglone is supplied as a solid. A stock solution may be made by dissolving the 2,7-dimethoxy-6-(1-acetoxyethyl)juglone in the solvent of choice, which should be purged with an inert gas. 2,7-Dimethoxy-6-(1-acetoxyethyl)juglone is soluble in organic solvents such as ethanol, methanol, DMSO, and dichloromethane.

Description

2,7-Dimethoxy-6-(1-acetoxyethyl)juglone is a naphthoquinone derivative originally isolated from the fungus *H. toruloidea*.¹

Reference

1. van Eijk, G.W. and Roeymans, H.J. Naphthoquinone derivatives from the fungus *Hendersonula toruloidea*. *Experientia* **34**(10), 1257-1258 (1978).

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, 08/23/2022

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