

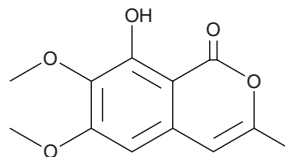
PRODUCT INFORMATION



6-O-Methylreticulol

Item No. 40221

CAS Registry No.: 24350-93-4
Formal Name: 8-hydroxy-6,7-dimethoxy-3-methyl-1H-2-benzopyran-1-one
Synonym: 8-hydroxy-6,7-dimethoxy-3-methyl Isocoumarin
MF: C₁₂H₁₂O₅
FW: 236.2
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥1 year
Item Origin: Bacterium/*Streptomyces mobaraensis*



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

Laboratory Procedures

6-O-Methylreticulol is supplied as a solid. A stock solution may be made by dissolving the 6-O-methylreticulol in the solvent of choice, which should be purged with an inert gas. 6-O-Methylreticulol is soluble in methanol, DMSO, and dimethyl formamide.

Description

6-O-Methylreticulol is an isocoumarin and a fungal metabolite that has been found in *S. mobaraensis*.¹

Reference

1. Saeed, A. and Ehsan, S. An efficient synthesis of 8-hydroxy-6,7-dimethoxy-3-methylisocoumarin (6-O-methylreticulol). *Chem. Heterocycl. Compd. (N Y)* **41(11)**, 1381-1385 (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, 11/08/2023

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