

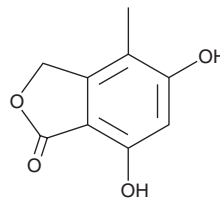
PRODUCT INFORMATION



5,7-Dihydroxy-4-methylphthalide

Item No. 43303

CAS Registry No.: 27979-57-3
Formal Name: 5,7-dihydroxy-4-methyl-1(3H)-isobenzofuranone
Synonym: DHMP
MF: $C_9H_8O_4$
FW: 180.2
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

5,7-Dihydroxy-4-methylphthalide is supplied as a solid. A stock solution may be made by dissolving the 5,7-dihydroxy-4-methylphthalide in the solvent of choice, which should be purged with an inert gas. 5,7-Dihydroxy-4-methylphthalide is slightly soluble (0.1-1 mg/ml) in ethanol and DMSO.

Description

5,7-Dihydroxy-4-methylphthalide is a precursor in the biosynthesis of mycophenolic acid (Item No. 21716) in fungi.¹ It is active against *C. neoformans* ($IC_{50} = 18.08 \mu\text{g/ml}$).² 5,7-Dihydroxy-4-methylphthalide inhibits LPS- or concanavalin A-induced proliferation of mouse splenic lymphocytes ($IC_{50}s = 3$ and $17 \mu\text{M}$, respectively).³

References

1. Hansen, B.G., Mnich, E., Nielsen, K.F., et al. Involvement of a natural fusion of a cytochrome P450 and a hydrolase in mycophenolic acid biosynthesis. *Appl. Environ. Microb.* **78**(14), 4908-4913 (2012).
2. Gao, J., Radwan, M.M., León, F., et al. Antimicrobial and antiprotozoal activities of secondary metabolites from the fungus *Eurotium repens*. *Med. Chem. Res.* **21**(10), 3080-3086 (2012).
3. Fujimoto, H., Fujimaki, T., Okuyama, E., et al. Immunomodulatory constituents from an ascomycete, *Microascus tardifaciens*. *Chem. Pharm. Bull. (Tokyo)* **47**(10), 1426-1432 (1999).

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, 04/28/2025

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