

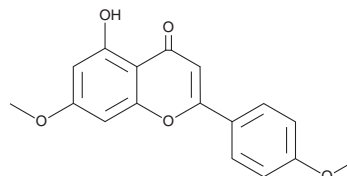
PRODUCT INFORMATION



7,4'-Di-O-methylapigenin

Item No. 35388

CAS Registry No.: 5128-44-9
Formal Name: 5-hydroxy-7-methoxy-2-(4-methoxyphenyl)-4H-1-benzopyran-4-one
Synonyms: Acacetin 7-methyl ether, Apigenin 7,4'-dimethyl ether, Dimethoxylapigenin, Genkwanin 4'-methyl ether, 5-Hydroxy-7,4'-dimethoxyflavone, 4'-Methoxytecto-chrysin
MF: C₁₇H₁₄O₅
FW: 298.3
Purity: ≥98%
UV/Vis.: λ_{max}: 269, 324 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Semisynthetic



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

Laboratory Procedures

7,4'-Di-O-methylapigenin is supplied as a solid. A stock solution may be made by dissolving the 7,4'-di-O-methylapigenin in the solvent of choice, which should be purged with an inert gas. 7,4'-Di-O-methylapigenin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 7,4'-di-O-methylapigenin in these solvents is approximately 1 and 3 mg/ml, respectively.

Description

7,4'-Di-O-methylapigenin is a flavonoid that has been found in *T. diffusa* and has diverse biological activities.¹⁻⁵ It inhibits monoamine oxidase B (MAO-B) but not MAO-A (IC₅₀s = 0.198 and >100 μM, respectively), as well as p38α MAPK but not JNK3 (IC₅₀s = 27.8 and >100 μM, respectively).^{1,2} 7,4'-Di-O-methylapigenin inhibits the production of nitric oxide (NO) in LPS-stimulated RAW 264.7 macrophages (IC₅₀ = 24.5 μM).³ It reduces the growth of *C. albicans* when used at a concentration of 45 μg/ml.⁴ 7,4'-Di-O-methylapigenin inhibits carrageenan-induced edema in rats (ED₂₅ = 75 mg/kg).⁵

References

1. Chaurasiya, N.D., Zhao, J., Pandey, P., et al. *Molecules* **24**(4), 810 (2019).
2. Goettert, M., Schattel, V., Koch, P., et al. *Chembiochem* **11**(18), (2010).
3. Tewtrakul, S. and Subhadhirasakul, S. *J. Ethnopharmacol.* **120**(1), 81-84 (2008).
4. Mangoyi, R., Midiwo, J., and Mukanganyama, S. *BMC Complement. Altern. Med.* **15**, 405 (2015).
5. Fourie, R.G. and Snyckers, F.O. *J. Nat. Prod.* **47**(6), 1057-1058 (1984).

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, 11/23/2022

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