

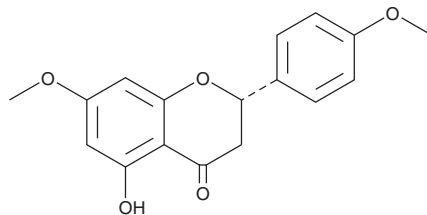
PRODUCT INFORMATION



4',7-Di-O-methylnaringenin

Item No. 36190

CAS Registry No.: 29424-96-2
Formal Name: 2S,3-dihydro-5-hydroxy-7-methoxy-2-(4-methoxyphenyl)-4H-1-benzopyran-4-one
MF: C₁₇H₁₆O₅
FW: 300.3
Purity: ≥98%
UV/Vis.: λ_{max}: 215, 227, 287 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Aglaia odorata*



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

Laboratory Procedures

4',7-Di-O-methylnaringenin is supplied as a solid. A stock solution may be made by dissolving the 4',7-di-O-methylnaringenin in the solvent of choice, which should be purged with an inert gas. 4',7-Di-O-methylnaringenin is soluble in DMSO.

Description

4',7-Di-O-methylnaringenin is a flavonoid aglycone that has been found in *Vitex quinata* and has potassium channel inhibitory and anticancer activities.^{1,2} It inhibits the voltage-gated potassium channel subtype K_v1.3 in a patch-clamp assay when used at a concentration of 30 μM.¹ 4',7-Di-O-methylnaringenin is cytotoxic to human LNCaP prostate, Lu1 lung, and MCF-7 breast cancer cells (EC₅₀s = 6.7, 4.7, and 1.1 μM, respectively).²

References

1. Teisseyre, A., Duarte, N., Ferreira, M.-J.U., *et al.* Influence of the multidrug transporter inhibitors on the activity of Kv1.3 voltage-gated potassium channels. *J. Physiol. Pharmacol.* **60**(1), 69-76 (2009).
2. Deng, Y., Chin, Y.-W., Chai, H.-B., *et al.* Phytochemical and bioactivity studies on constituents of the leaves of *Vitex quinata*. *Phytochem. Lett.* **4**(3), 213-217 (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, 11/07/2022

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