

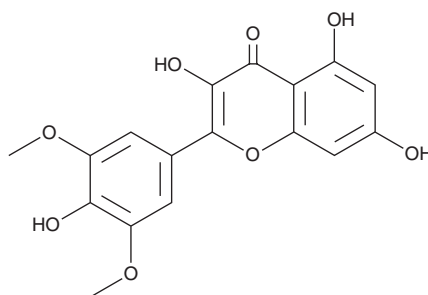
PRODUCT INFORMATION



Syringetin

Item No. 39249

CAS Registry No.: 4423-37-4
Formal Name: 3,5,7-trihydroxy-2-(4-hydroxy-3,5-dimethoxyphenyl)-4H-1-benzopyran-4-one
Synonym: Myricetin-3',5'-dimethyl ether
MF: C₁₇H₁₄O₈
FW: 346.3
Purity: ≥98%
UV/Vis.: λ_{max}: 253, 375 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Syringa oblata* Lindl



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

Laboratory Procedures

Syringetin is supplied as a solid. A stock solution may be made by dissolving the syringetin in the solvent of choice, which should be purged with an inert gas. Syringetin is soluble in DMSO.

Description

Syringetin is a methylated flavonol that has been found in *G. biloba* and has diverse biological activities.¹⁻³ It inhibits α-glucosidase and scavenges radicals in a Trolox equivalent antioxidant capacity (TEAC) assay (IC₅₀s = 36.8 and 12.3 μM, respectively).^{1,3} Syringetin (50 μM) reduces the proliferation of, induces cell cycle arrest at the G₂/M phase in, and decreases cyclin D1 and COX-2 levels in, Caco-2 colorectal adenocarcinoma cells.² It increases the lifespan of *C. elegans* when used at a concentration of 100 μM.³

References

1. Wu, B., Song, H.-P., Zhou, X., *et al.* Screening of minor bioactive compounds from herbal medicines by *in silico* docking and the trace peak exposure methods. *J. Chromatogr. A* **1436**, 91-99 (2016).
2. Gómez-Alonso, S., Collins, V.J., Vauzour, D., *et al.* Inhibition of colon adenocarcinoma cell proliferation by flavonols is linked to a G₂/M cell cycle block and reduction in cyclin D1 expression. *Food Chem.* **130(3)**, 493-500 (2012).
3. Büchter, C., Ackermann, D., Honnen, S., *et al.* Methylated derivatives of myricetin enhance life span in *Caenorhabditis elegans* dependent on the transcription factor DAF-16. *Food Funct.* **6(10)**, 3383-3392 (2015).

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, 10/09/2023

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