

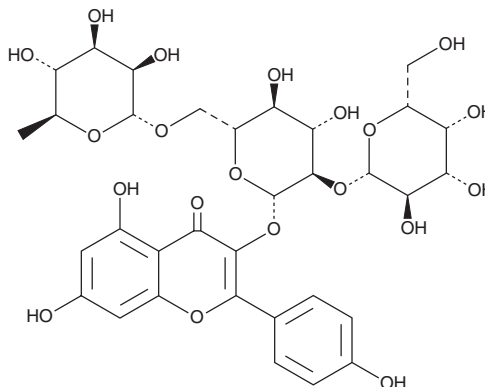
PRODUCT INFORMATION



Camelliaside A

Item No. 35628

CAS Registry No.: 135095-52-2
Formal Name: 3-[[O-6-deoxy- α -L-mannopyranosyl-(1 \rightarrow 6)-O-[β -D-galactopyranosyl-(1 \rightarrow 2)]- β -D-glucopyranosyl]oxy]-5,7-dihydroxy-2-(4-hydroxyphenyl)-4H-1-benzopyran-4-one
MF: C₃₃H₄₀O₂₀
FW: 756.7
Purity: \geq 98%
UV/Vis.: λ_{max} : 214, 263, 330 nm
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/*Camellia oleifera*



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

Laboratory Procedures

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

Description

Camelliaside A is a flavonoid that has been found in *C. sinensis* and has radical scavenging and enzyme inhibitory activities.¹⁻⁴ It scavenges superoxide radicals in a cell-free assay (IC₅₀ = 137.44 μ M).¹ Camelliaside A (1 μ M) inhibits recombinant human monoamine oxidase B (MAO-B) in a cell-free assay.⁴ It also inhibits 5-lipoxygenase (5-LO) in RBL-1 rat basophilic leukemia cells (IC₅₀ = 390 μ M).³

References

1. Kang, B., Chang, H., Na, Y.J., *et al.* Extract of enzyme-hydrolyzed green tea seed as potent melanin synthesis inhibitor. *Bull. Korean Chem. Soc.* **34**(7), 2199-2202 (2013).
2. Park, J.S., Rho, H.S., Kim, D.H., *et al.* Enzymatic preparation of kaempferol from green tea seed and its antioxidant activity. *J. Agric. Food Chem.* **54**(8), 2951-2956 (2006).
3. Sekine, T., Arai, Y., Ikegami, F., *et al.* Isolation of camelliaside C from "tea seed cake" and inhibitory effects of its derivatives on arachidonate 5-lipoxygenase. *Chem. Pharm. Bull. (Tokyo)* **41**(6), 1185-1187 (1993).
4. Shkondrov, A., Krasteva, I., Bucar, F., *et al.* A new tetracyclic saponin from *Astragalus glycyphyllos* L. and its neuroprotective and hMAO-B inhibiting activity. *Nat. Prod. Res.* **34**(4), 511-517 (2020).

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, 09/28/2022

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