

PRODUCT INFORMATION



Clitorin

Item No. 33211

CAS Registry No.: 55804-74-5

Formal Name: 3-[[O-6-deoxy- α -L-mannopyranosyl-(1 \rightarrow 2)-O-[6-deoxy- α -L-mannopyranosyl-(1 \rightarrow 6)]- β -D-glucopyranosyl]oxy]-5,7-dihydroxy-2-(4-hydroxyphenyl)-4H-1-benzopyran-4-one

Synonyms: Kaempferol 3-(2^G-rhamnosylrutinoside), Kaempferol 3-O-(2,6- α -L-dirhamnopyranosyl- β -D-glucopyranoside)

MF: C₃₃H₄₀O₁₉

FW: 740.7

Purity: \geq 98%

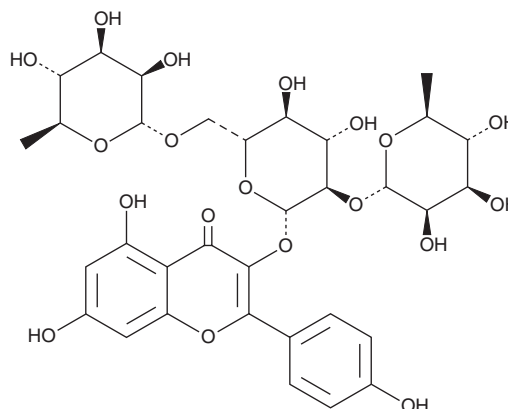
UV/Vis.: λ_{\max} : 265, 347 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: \geq 4 years

Item Origin: Plant/*Lysimachia christinae* Hance



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

Laboratory Procedures

Clitorin is supplied as a crystalline solid. A stock solution may be made by dissolving the clitorin in the solvent of choice, which should be purged with an inert gas. Clitorin is soluble in organic solvents such as chloroform and DMSO.

Description

Clitorin is a flavonol glycoside that has been found in *Clitoria*.¹ It has antioxidant activity in an online ABTS⁺-capillary electrophoresis-diode array detector (ABTS⁺-CE-DAD) antioxidant screening assay but is inactive in a DPPH (Item No. 14805) scavenging assay.^{2,3}

References

1. Kazuma, K., Noda, N., and Suzuki, M. Malonylated flavonol glycosides from the petals of *Clitoria ternatea*. *Phytochemistry* **62**(2), 229-237 (2003).
2. Ma, H., Li, J., An, M., et al. A powerful on line ABTS⁺-CE-DAD method to screen and quantify major antioxidants for quality control of Shuxuening Injection. *Sci. Rep.* **8**(1), 5441 (2018).
3. Chevalley, I., Marston, A., and Hostettmann, K. New phenolic radical scavengers from *Saxifraga cuneifolia*. *Pharm. Biol.* **38**(3), 222-228 (2000).

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, 12/13/2022

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