

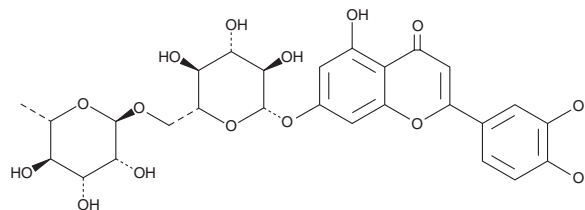
PRODUCT INFORMATION



Luteolin 7-O-Rutinoside

Item No. 34035

CAS Registry No.: 20633-84-5
Formal Name: 7-[[6-O-(6-deoxy- α -L-mannopyranosyl)- β -D-glucopyranosyl]oxy]-2-(3,4-dihydroxyphenyl)-5-hydroxy-4H-1-benzopyran-4-one
Synonyms: Luteolin 7-Rutinoside, Luteolin 7- β -Rutinoside, Scolymoside, Skolimoside
MF: C₂₇H₃₀O₁₅
FW: 594.5
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Synthetic



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

Laboratory Procedures

Luteolin 7-O-rutinoside is supplied as a solid. A stock solution may be made by dissolving the luteolin 7-O-rutinoside in the solvent of choice, which should be purged with an inert gas. Luteolin 7-O-rutinoside is soluble in the organic solvent DMSO.

Description

Luteolin 7-O-rutinoside is a polyketide-derived flavonoid glycoside that has been found in *V. bugulifolium* and has diverse biological activities.^{1-2,4} It scavenges DPPH (Item No. 14805) radicals in a cell-free assay when used at concentrations ranging from 1 to 50 μ M.¹ Luteolin 7-O-rutinoside is active against *B. subtilis*, *S. aureus*, *A. tumefaciens*, *M. luteus*, *E. coli*, and *P. aeruginosa* (MICs = 100-200 μ g/ml).² It is also active against *C. albicans*, *S. cerevisiae*, and *C. lusitaniae* (MICs = 100, 200, and 50 μ g/ml, respectively). Luteolin 7-O-rutinoside (57.6 μ g/animal) decreases hepatic and renal injury and increases survival in a mouse model of polyphosphate-induced lethal endotoxemia.³

References

1. Wang, W., Simon, J.E., Aviles, I.F., et al. Analysis of antioxidative phenolic compounds in artichoke (*Cynara scolymus* L.). *J. Agric. Food Chem.* **51**(3), 601-608 (2003).
2. Zhu, X., Zhang, H., and Lo, R. Phenolic compounds from the leaf extract of artichoke (*Cynara scolymus* L.) and their antimicrobial activities. *J. Agric. Food Chem.* **52**(25), 7272-7278 (2004).
3. Lee, I.-C. and Bae, J.-S. Anti-inflammatory effects of vicenin-2 and scolymoside on polyphosphate-mediated vascular inflammatory responses. *Inflamm. Res.* **65**(3), 203-212 (2016).
4. Fraga, C.G. *Plant Phenolics and Human Health: Biochemistry, Nutrition and Pharmacology*. 1st ed., John Wiley & Sons, Inc., Hoboken, New Jersey (2010).

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, 08/25/2023

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