

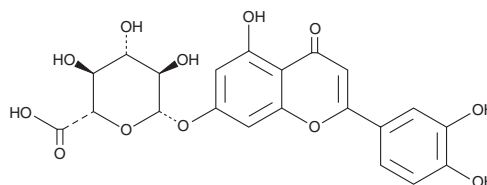
PRODUCT INFORMATION



Luteolin 7-O-Glucuronide

Item No. 33188

CAS Registry No.: 29741-10-4
Formal Name: 2-(3,4-dihydroxyphenyl)-5-hydroxy-4-oxo-4H-1-benzopyran-7-yl, β -D-glucopyranosiduronic acid
Synonym: Luteolin-7-O- β -D-glucuronide
MF: C₂₁H₁₈O₁₂
FW: 462.4
Purity: \geq 98%
UV/Vis.: λ_{max} : 256, 353 nm
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/*Lactuca sativa*



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

Laboratory Procedures

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

Description

Luteolin 7-O-glucuronide is a flavonoid that has been found in *S. chloroleuca* and *A. pilosa* and has diverse biological activities, including antioxidant, enzyme inhibitory, and anti-inflammatory properties.¹⁻³ It scavenges DPPH (Item No. 14805) radicals in a cell-free assay (IC₅₀ = 80.6 μ M).¹ Luteolin 7-O-glucuronide inhibits porcine α -amylase and yeast α -glucosidase *in vitro* (IC₅₀s = 61.5 and 14.7 μ M, respectively).³ It also inhibits rat aldose reductase *in vitro* (IC₅₀ = 0.7 μ M) and sorbitol accumulation in isolated rat lens by 91.8% when used at a concentration of 5 μ g/ml.¹ Luteolin 7-O-glucuronide inhibits LPS-induced cytokine release from RAW 264.7 macrophages in a dose-dependent manner.²

References

1. Kim, S.B., Hwang, S.H., Suh, H.-W., *et al.* Phytochemical analysis of *Agrimonia pilosa* Ledeb, its antioxidant activity and aldose reductase inhibitory potential. *Int. J. Mol. Sci.* **18**(2), 379 (2017).
2. Cho, Y.-C., Park, J., and Cho, S. Anti-inflammatory and anti-oxidative effects of luteolin-7-O-glucuronide in LPS-stimulated murine macrophages through TAK1 inhibition and Nrf2 activation. *Int. J. Mol. Sci.* **21**(6), 2007 (2020).
3. Asghari, B., Salehi, P., Sonboli, A., *et al.* Flavonoids from *Salvia chloroleuca* with α -amylase and α -glucosidase inhibitory effect. *Iran. J. Pharm. Res.* **14**(2), 609-615 (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, 11/29/2022

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