

PRODUCT INFORMATION

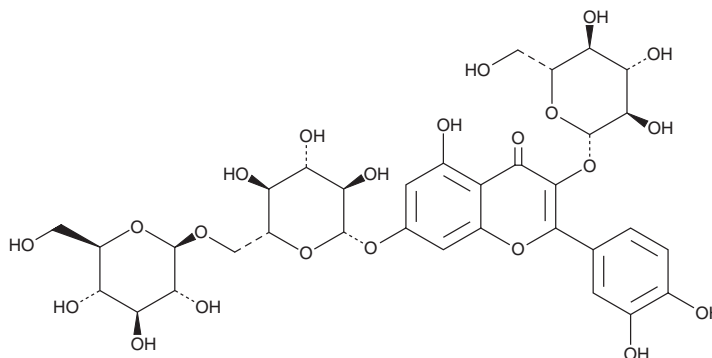


Quercetin 3-O-β-D-Glucose-7-O-β-D-Gentiobioside

Item No. 36130

CAS Registry No.: 60778-02-1
Formal Name: 2-(3,4-dihydroxyphenyl)-7-[(6-O-β-D-glucopyranosyl-β-D-glucopyranosyl)oxy]-3-(β-D-glucopyranosyloxy)-5-hydroxy-4H-1-benzopyran-4-one

MF: C₃₃H₄₀O₂₂
FW: 788.7
Purity: ≥98%
UV/Vis.: λ_{max}: 258, 363 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Desurainia sophia*



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

Laboratory Procedures

Quercetin 3-O-β-D-glucose-7-O-β-D-gentiobioside is supplied as a solid. A stock solution may be made by dissolving the quercetin 3-O-β-D-glucose-7-O-β-D-gentiobioside in the solvent of choice, which should be purged with an inert gas. Quercetin 3-O-β-D-glucose-7-O-β-D-gentiobioside is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of quercetin 3-O-β-D-glucose-7-O-β-D-gentiobioside in these solvents is approximately 3 and 2 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of quercetin 3-O-β-D-glucose-7-O-β-D-gentiobioside can be prepared by directly dissolving the solid in aqueous buffers. The solubility of quercetin 3-O-β-D-glucose-7-O-β-D-gentiobioside in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Quercetin 3-O-β-D-glucose-7-O-β-D-gentiobioside is a flavonoid glycoside that has been found in *L. apetalum* and has lipid-lowering activity.¹ It reduces sodium oleate-induced increases in triglyceride levels in HepG2 cells when used at a concentration of 30 μM.

Reference

1. Wang, S., Shi, P., Qu, L., *et al.* Bioactive constituents obtained from the seeds of *Lepidium apetalum* Willd. *Molecules* **22**(4), 540 (2017).

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, 12/19/2022

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