

PRODUCT INFORMATION



Apigenin-7-O-(2G-rhamnosyl)gentiobioside

Item No. 35893

CAS Registry No.: 174284-20-9
Formal Name: 7-[[O-6-deoxy- α -L-mannopyranosyl-(1 \rightarrow 2)-O-[β -D-glucopyranosyl-(1 \rightarrow 6)]]- β -D-glucopyranosyl]oxy]-5-hydroxy-2-(4-hydroxyphenyl)-4H-1-benzopyran-4-one

MF: C₃₃H₄₀O₁₉

FW: 740.7

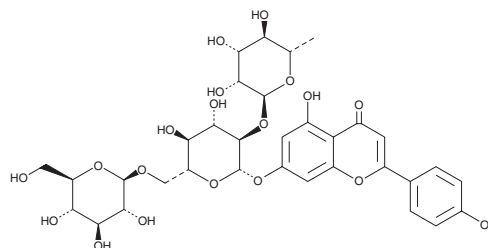
Purity: \geq 98%

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years

Item Origin: Plant/*Eleutherococcus sessiliflorus* (Rupr. & Maxim.) S.Y.Hu



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

Laboratory Procedures

Apigenin-7-O-(2G-rhamnosyl)gentiobioside is supplied as a solid. A stock solution may be made by dissolving the apigenin-7-O-(2G-rhamnosyl)gentiobioside in the solvent of choice, which should be purged with an inert gas. Apigenin-7-O-(2G-rhamnosyl)gentiobioside is soluble in chloroform. Apigenin-7-O-(2G-rhamnosyl)gentiobioside is slightly soluble in methanol and DMSO.

Description

Apigenin-7-O-(2G-rhamnosyl)gentiobioside is a flavonoid glycoside that has been found in *L. gracilipes*.¹

Reference

1. Kikuchi, M. and Matsuda, N. Flavone glycosides from *Lonicera gracilipes* var. *glandulosa*. *J. Nat. Prod.* **59**(3), 314-315 (1996).

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, 02/16/2024

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