

PRODUCT INFORMATION



(-)-Syringaresinol di-O-glucoside

Item No. 36672

CAS Registry No.: 66791-77-3
Formal Name: [(1R,3aS,4R,6aS)-tetrahydro-1H,3H-furo[3,4-c]furan-1,4-diyl]bis(2,6-dimethoxy-4,1-phenylene) bis-β-D-glucopyranoside
Synonym: (-)-Syringaresinol-4,4'-bis-O-β-D-glucopyranoside

MF: C₃₄H₄₆O₁₈
FW: 742.7

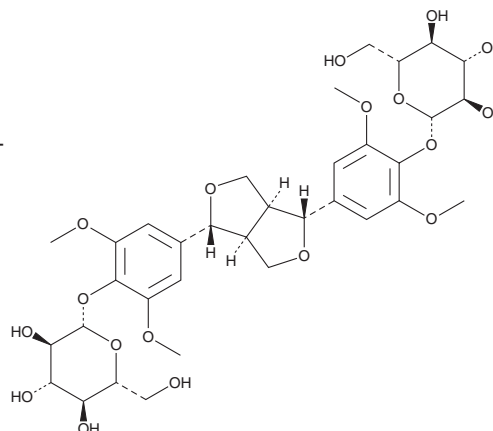
Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/*Eleutherococcus senticosus*



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

Laboratory Procedures

(-)-Syringaresinol di-O-glucoside is supplied as a solid. A stock solution may be made by dissolving the (-)-syringaresinol di-O-glucoside in the solvent of choice. (-)-Syringaresinol di-O-glucoside is soluble in organic solvents such as methanol, which should be purged with an inert gas. (-)-Syringaresinol di-O-glucoside is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

(-)-Syringaresinol di-O-glucoside is a phenolic glucoside that has been found in *D. aurantiacum* and has neuroprotective activity.¹ It reduces neurotoxicity induced by glutamate by 19.3% in PC12 cells when used at a concentration of 10 μM.

Reference

1. Xiong, L., Cao, Z.-X., Peng, C., *et al.* Phenolic glucosides from *Dendrobium aurantiacum* var. *denneanum* and their bioactivities. *Molecules* **18**(6), 6153-6160 (2013).

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, 12/21/2022

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