

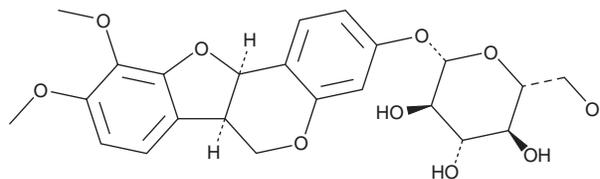
PRODUCT INFORMATION



9,10-Dimethoxypterocarpan-3-O-β-D-glucoside

Item No. 36192

CAS Registry No.: 94367-42-7
Formal Name: 6aR,11aR-dihydro-9,10-dimethoxy-6H-benzofuro[3,2-c][1]benzopyran-3-yl β-D-glucopyranoside
Synonyms: 7-β-D-glucopyranosyloxy-Astrapterocarpan, Methylnissolin 3-O-glucoside
MF: C₂₃H₂₆O₁₀
FW: 462.5
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Astragalus sinicus*



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

Laboratory Procedures

9,10-Dimethoxypterocarpan-3-O-β-D-glucoside is supplied as a solid. A stock solution may be made by dissolving the 9,10-dimethoxypterocarpan-3-O-β-D-glucoside in the solvent of choice, which should be purged with an inert gas. 9,10-Dimethoxypterocarpan-3-O-β-D-glucoside is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 9,10-dimethoxypterocarpan-3-O-β-D-glucoside in these solvents is approximately 2 mg/ml.

Description

9,10-Dimethoxypterocarpan-3-O-β-D-glucoside is a flavonoid glycoside that has been found in *A. membranaceus* and has anti-inflammatory activity.¹ It inhibits LPS-induced production of nitric oxide (NO) in RAW 264.7 macrophages (EC₅₀ = 14.7 μM).

Reference

1. Zhang, L.-J., Liu, H.-K., Hsiao, P.-C., *et al.* New isoflavonoid glycosides and related constituents from astragali radix (*Astragalus membranaceus*) and their inhibitory activity on nitric oxide production. *J. Agric. Food Chem.* 59(4), 1131-1137 (2011).

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