

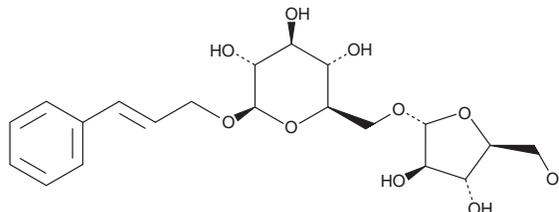
PRODUCT INFORMATION



Rosarin

Item No. 35214

CAS Registry No.: 84954-93-8
Formal Name: (2E)-3-phenyl-2-propen-1-yl
6-O- α -L-arabinofuranosyl- β -D-
glucopyranoside
MF: C₂₀H₂₈O₁₀
FW: 428.4
Purity: \geq 98%
UV/Vis.: λ_{max} : 251 nm
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/*Rhodiola rosea*



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

Laboratory Procedures

Rosarin is supplied as a solid. A stock solution may be made by dissolving the rosarin in the solvent of choice, which should be purged with an inert gas. Rosarin is soluble in DMSO.

Description

Rosarin is a glycoside that has been found in *R. rosea* and has anti-inflammatory and immune regulatory activities.¹ It prevents LPS-induced increases in induced nitric oxide synthase (iNOS), TNF- α , IL-1 β , and IL-6 levels in BV-2 microglia when used at concentrations of 10 and 50 μ g/ml. Rosarin reduces LPS-induced nitric oxide (NO) production in BV-2 microglia. It inhibits the growth of Jurkat T cells induced by concanavalin A (Item No. 14951) or anti-CD3 and anti-CD28 antibodies (IC₅₀s = 60 and 68 μ M, respectively) and increases TRAIL levels in isolated mouse splenic T cells when used at concentrations of 0.5, 5, and 50 μ M.²

References

1. Lee, Y., Jung, J.-C., Jang, S., *et al.* Anti-inflammatory and neuroprotective effects of constituents isolated from *Rhodiola rosea*. *Evid. Based Complement. Alternat. Med.* 514049 (2013).
2. Marchev, A.S., Dimitrova, P., Koycheva, I.K., *et al.* Altered expression of TRAIL on mouse T cells via ERK phosphorylation by *Rhodiola rosea* L. and its marker compounds. *Food Chem. Toxicol.* **108(Pt B)**, 419-428 (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, 02/07/2024

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