

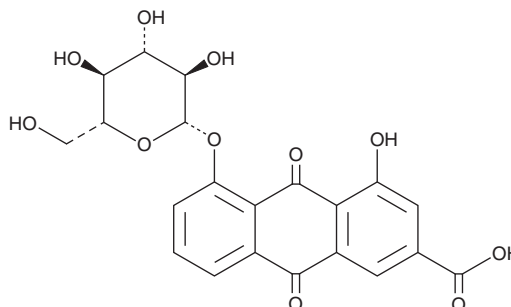
PRODUCT INFORMATION



Rhein-8-glucoside

Item No. 28844

CAS Registry No.: 34298-86-7
Formal Name: 5-(β-D-glucopyranosyloxy)-9,10-dihydro-4-hydroxy-9,10-dioxo-2-anthracenecarboxylic acid
Synonym: Rhein 8-O-β-D-glucopyranoside
MF: C₂₁H₁₈O₁₁
FW: 446.4
Purity: ≥98%
UV/Vis.: λ_{max}: 227, 284, 428 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Rheum palmatum* L.



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

Laboratory Procedures

Rhein-8-glucoside is supplied as a crystalline solid. A stock solution may be made by dissolving the rhein-8-glucoside in the solvent of choice, which should be purged with an inert gas. Rhein-8-glucoside is slightly soluble in DMSO.

Description

Rhein-8-glucoside is an anthraquinone glycoside that has been found in rhubarb.¹ It increases the rate of sennoside A (Item No. 24969) metabolism by mouse intestinal bacteria *in vitro* when used at concentrations ranging from 0.03 to 1 mM. Rhein-8-glucoside increases the purgative activity of sennoside A in mice in a dose-dependent manner. It also inhibits human protein tyrosine phosphatase (PTP1B; IC₅₀ = 11.5 μM).² Rhein-8-glucoside is metabolized to rhein (Item No. 17345) by intestinal bacteria.¹

References

1. Takayama, K., Tsutsumi, H., Ishizu, T., *et al.* The influence of rhein 8-O-β-D-glucopyranoside on the purgative action of sennoside A from rhubarb in mice. *Biol. Pharm. Bull.* **35(12)**, 2204-2208 (2012).
2. Li, S., An, T.-Y., Li, J., *et al.* PTP1B inhibitors from *Saussurea lappa*. *J. Asian Nat. Prod. Res.* **8(3)**, 281-286 (2006).

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, 12/08/2022

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