

PRODUCT INFORMATION

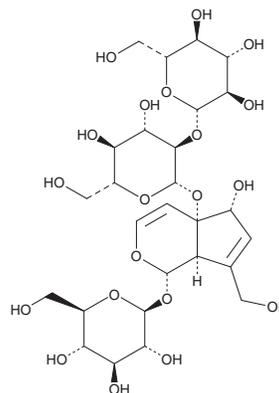


Rehmannioside D

Item No. 29034

CAS Registry No.: 81720-08-3
Formal Name: (1S,4aS,5R,7aR)-1-(β-D-glucopyranosyloxy)-5,7a-dihydro-5-hydroxy-7-(hydroxymethyl)cyclopenta[c]pyran-4a(1H)-yl 2-O-β-D-glucopyranosyl-β-D-glucopyranoside

MF: C₂₇H₄₂O₂₀
FW: 686.6
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Rehmanniae Radix*



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

Laboratory Procedures

Rehmannioside D is supplied as a crystalline solid. A stock solution may be made by dissolving the rehmannioside D in the solvent of choice, which should be purged with an inert gas. Rehmannioside D is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of rehmannioside D in these solvents is approximately 20 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of rehmannioside D can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of rehmannioside D in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Rehmannioside D is an iridoid glycoside that has been found in *R. glutinosa*.¹

Reference

1. Matsumoto, M., Shoyama, Y., Nishioka, I., *et al.* Identification of viruses infected in *Rehmannia glutinosa* Libosch. var *purpurea* Makino and effect of virus infection on root yield and iridoid glycoside contents. *Plant Cell Rep.* **7(8)**, 636-638 (1989).

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, 11/23/2022

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