

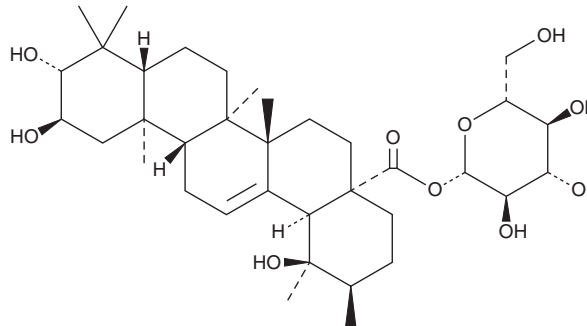
PRODUCT INFORMATION



Rosamultin

Item No. 39939

CAS Registry No.: 88515-58-6
Formal Name: 2 α ,3 β ,19-trihydroxy-urs-12-en-28-oic acid, β -D-glucopyranosyl ester
Synonyms: Tormentic Acid 28-O-glucoside ester, Ursolazuroside 1
MF: C₃₆H₅₈O₁₀
FW: 650.8
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Argentina anserina* (L.) Rydb.



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

Laboratory Procedures

Rosamultin is supplied as a solid. A stock solution may be made by dissolving the rosamultin in the solvent of choice, which should be purged with an inert gas. Rosamultin is sparingly soluble (1-10 mg/ml) in ethanol and slightly soluble (0.1-1 mg/ml) in DMSO.

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 rosamultin can be prepared by directly dissolving the solid in aqueous buffers. Rosamultin is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Rosamultin is a triterpenoid glycoside that has been found in *R. rugosa* and has diverse biological activities.¹⁻⁴ It inhibits HIV-1 protease by 52.9% in a cell-free assay when used at a concentration of 100 μ M.¹ Rosamultin (0.01 nM) reduces hydrogen peroxide-induced apoptosis, increases in malondialdehyde (MDA) levels, and decreases in superoxide dismutase (Sod), catalase (Cat), and glutathione peroxidase (Gpx) activities in H9c2 rat cardiomyocytes.² *In vivo*, rosamultin (12 mg/kg) increases survival, prevents decreases in body weight, and protects against splenic and intestinal damage in irradiated mice.³ Rosamultin (30 mg/kg, p.o.) inhibits acetic acid-induced writhing in mice.⁴

References

1. Park, J.C., Kim, S.C., Choi, M.R., *et al.* Anti-HIV protease activity from rosa family plant extracts and rosamultin from *Rosa rugosa*. *J. Med. Food* **8**(1), 107-109 (2005).
2. Zhang, L., Liu, Y., Li, J.Y., *et al.* Protective effect of rosamultin against H₂O₂-induced oxidative stress and apoptosis in H9c2 cardiomyocytes. *Oxid. Med. Cell. Longev.* 8415610 (2018).
3. Liu, N., Niu, M., Luo, S., *et al.* Rosamultin ameliorates radiation injury via promoting DNA injury repair and suppressing oxidative stress *in vitro* and *in vivo*. *Chem. Biol. Interact.* **393**, 110938 (2024).
4. Jung, H.-J., Nam, J.-H., Choi, J., *et al.* 19 α -hydroxyursane-type triterpenoids: Antinociceptive anti-inflammatory principles of the roots of *Rosa rugosa*. *Biol. Pharm. Bull.* **28**(1), 101-104 (2005).

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, 08/20/2024

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