

PRODUCT INFORMATION



Raddeanin A

Item No. 40235

CAS Registry No.: 89412-79-3
Formal Name: (3 β)-3-[(O-6-deoxy- α -L-mannopyranosyl-(1 $\rightarrow$ 2)-O- β -D-glucopyranosyl-(1 $\rightarrow$ 2)- α -L-arabinopyranosyl)oxy]-olean-12-en-28-oic acid

Synonym: NSC 382873

MF: C₄₇H₇₆O₁₆

FW: 897.1

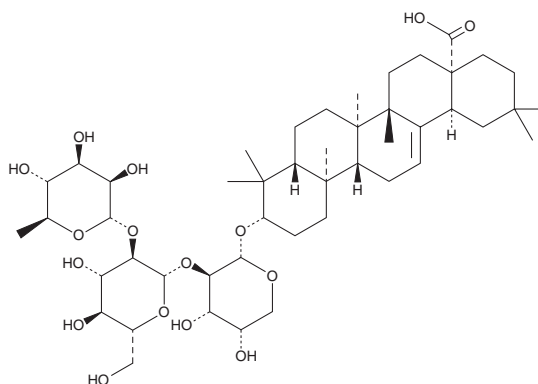
Purity: $\geq 98\%$

Supplied as: solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Plant/*Anemone raddeana*



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

Laboratory Procedures

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

Description

Raddeanin A is a triterpenoid saponin that has been found in *A. raddeana* and has diverse biological activities.¹⁻³ It binds to amyloid- β (1-42) (A β 42; K_d = 27.63 μ M).² Raddeanin A (4 μ M) inhibits the invasion and migration of, and induces apoptosis and cell cycle arrest at the G₀/G₁ phase in, HeLa cervical cancer cells.¹ It improves learning and memory in the Morris water maze in mice in the 3xTg model of Alzheimer's disease when administered at a dose of 10 mg/kg.² Raddeanin A (50 mg/kg) reduces neurological deficits, lesion volume, edema, and microglia lesion infiltration in a mouse model of acute cerebral hemorrhage induced by intracerebral injection of bacterial collagenase VII.³

References

- Shen, X., Li, L., He, Y., *et al.* Raddeanin A inhibits proliferation, invasion, migration and promotes apoptosis of cervical cancer cells via regulating miR-224-3p/Slit2/Robo1 signaling pathway. *Aging (Albany NY)* 13(5), 7166-7179 (2021).
- Liu, M.H., Tang, Y., Qu, L.Q., *et al.* Raddeanin A isolated from *Anemone raddeana* Regel improves pathological and cognitive deficits of the mice model of Alzheimer's disease by targeting β -amyloidosis. *Phytomedicine* 122, 155121 (2024).
- Hei, B., Ouyang, J., Zhou, J., *et al.* Raddeanin A (RA) reduced acute inflammatory injury in mouse experimental cerebral hemorrhage by suppression of TLR4. *Int. J. Med. Sci.* 19(8), 1235-1240 (2022).

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, 02/29/2024

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