

PRODUCT INFORMATION



Hederasaponin B

Item No. 30641

CAS Registry No.: 36284-77-2

Formal Name: 3β-[[2-O-(6-deoxy-α-L-mannopyranosyl)-α-L-arabinopyranosyl]oxy]-olean-12-en-28-oic acid, O-6-deoxy-α-L-mannopyranosyl-(1→4)-O-β-D-glucopyranosyl-(1→6)-β-D-glucopyranosyl ester

Synonyms: Hederacoside B, Hederagenin B

MF: C₅₉H₉₆O₂₅

FW: 1,205.4

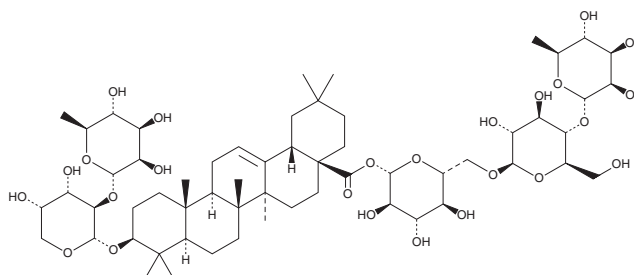
Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/Rhizoma *Anemones Raddeanae*



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

Laboratory Procedures

Hederasaponin B is supplied as a solid. A stock solution may be made by dissolving the hederasaponin B in the solvent of choice, which should be purged with an inert gas. Hederasaponin B is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of hederasaponin B in these solvents is approximately 30 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 hederasaponin B can be prepared by directly dissolving the solid in aqueous buffers. The solubility of hederasaponin B in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Hederasaponin B is a triterpene saponin that has been found in *H. helix* and has antiviral and antioxidant activities.^{1,2} It inhibits the cytopathic effect induced by the enterovirus 71 (EV71) genotypes C3 and C4a (EC₅₀s = 24.77 and 41.77 μg/ml, respectively), as well as reduces expression of the viral structural capsid protein VP2, in Vero cells.¹ Hederasaponin B (100, 150, and 200 μg/ml) inhibits superoxide generation in human neutrophils.²

References

1. Song, J., Yeo, S.-G., Hong, E.-H., et al. Antiviral activity of hederasaponin B from *Hedera helix* against enterovirus 71 subgenotypes C3 and C4a. *Biomol. Ther. (Seoul)* **22**(1), 41-46 (2014).
2. Chen, X., Lu, J., He, W., et al. Antiperoxidation activity of triterpenoids from rhizome of *Anemone raddeana*. *Fitoterapia* **80**(2), 105-111 (2009).

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, 12/12/2022

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