

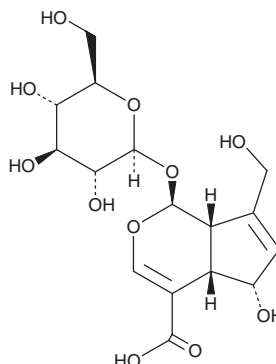
PRODUCT INFORMATION



Deacetylasperulosidic Acid

Item No. 25100

CAS Registry No.: 14259-55-3
Formal Name: 1-(β-D-glucopyranosyloxy)-1S,4aS,5S,7aS-tetrahydro-5-hydroxy-7-(hydroxymethyl)-cyclopenta[c]pyran-4-carboxylic acid
Synonym: 10-Deacetylasperulosidic Acid
MF: C₁₆H₂₂O₁₁
FW: 390.3
Purity: ≥98%
UV/Vis.: λ_{max}: 235 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Deacetylasperulosidic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the deacetylasperulosidic acid in the solvent of choice, which should be purged with an inert gas. Deacetylasperulosidic acid is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of deacetylasperulosidic acid in these solvents is approximately 10 and 5 mg/ml, respectively.

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 deacetylasperulosidic acid can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of deacetylasperulosidic acid in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Deacetylasperulosidic acid is a monoterpene glycoside originally isolated from *D. macropodum* that has antioxidant activity.^{1,2} It decreases serum malondialdehyde levels and increases superoxide dismutase activity in rats without affecting serum glutathione peroxidase activity when administered at doses of 30 and 50 mg/kg per day for seven days.²

References

1. Inoue, H., Ueda, S., Hirabayashi, M., *et al.* Studies on the monoterpene glucosides. IV. Monoterpene glucosides of *Daphniphyllum macropodum* MIQ. *Yakugaku Zasshi* **86(10)**, 943-947 (1966).
2. Ma, D.L., Chen, M., Su, C.X., *et al.* *In vivo* antioxidant activity of deacetylasperulosidic acid in noni. *J. Anal. Methods Chem.* **2013**, 804504 (2013).

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, 10/05/2022

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