

PRODUCT INFORMATION

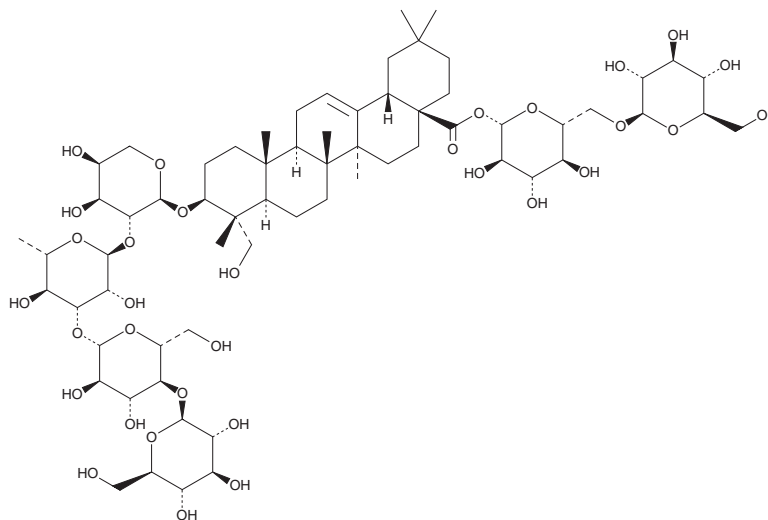


Macranthoidin B

Item No. 31533

CAS Registry No.: 136849-88-2
Formal Name: (3 β ,4 α)-3-[(O- β -D-glucopyranosyl-(1 \rightarrow 4)-O- β -D-glucopyranosyl-(1 \rightarrow 3)-O-6-deoxy- α -L-mannopyranosyl-(1 \rightarrow 2)- α -L-arabinopyranosyl)oxy]-23-hydroxy-olean-12-en-28-oic acid, 6-O- β -D-glucopyranosyl- β -D-glucopyranosyl ester

Synonym: Macranthoiside I
MF: C₆₅H₁₀₆O₃₂
FW: 1,399.5
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/*Lonicera japonica*



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

Laboratory Procedures

Macranthoidin B is supplied as a crystalline solid. A stock solution may be made by dissolving the macranthoidin B in the solvent of choice, which should be purged with an inert gas. Macranthoidin B is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of macranthoidin 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 macranthoidin B can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of macranthoidin 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

Macranthoidin B is a triterpenoid saponin that has been found in *Lonicerae* and has anticancer activity.¹ It increases the production of reactive oxygen species (ROS) and induces apoptosis in HCT116 cells when used at concentrations ranging from 20 to 400 μ M.

Reference

1. Fan, X., Rao, J., Zhang, Z., *et al.* Macranthoidin B modulates key metabolic pathways to enhance ROS generation and induce cytotoxicity and apoptosis in colorectal cancer. *Cell Physiol. Biochem.* **46(4)**, 1317-1330 (2018).

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, 12/08/2022

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