

PRODUCT INFORMATION



Amorfrutin B

Item No. 19618

CAS Registry No.: 78916-42-4
Formal Name: 3-[(2E)-3,7-dimethyl-2,6-octadien-1-yl]-2-hydroxy-4-methoxy-6-(2-phenylethyl)-benzoic acid

MF: C₂₆H₃₂O₄

FW: 408.5

Purity: ≥97%

Supplied as: A solid

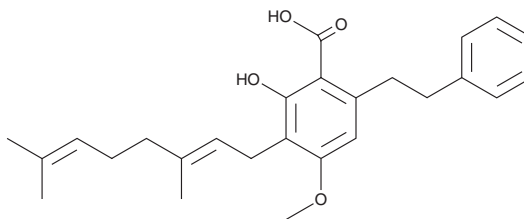
Storage: -20°C

Stability: ≥4 years

Special Conditions: Keep cool and dry. Protect from light.

Item Origin: Plant/*Amorpha fruticosa*

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



Laboratory Procedures

Amorfrutin B is supplied as a solid. A stock solution may be made by dissolving the amorfrutin B in the solvent of choice. Amorfrutin B is soluble in organic solvents such as ethanol and DMSO, which should be purged with an inert gas.

Description

Amorfrutin B is a partial agonist of the peroxisome proliferator-activated receptor γ (PPAR γ ; $K_i = 19$ nM and $EC_{50} = 73$ nM) that was first isolated from *A. fruticosa*.¹ It binds to PPAR α and PPAR β/δ with K_i values of 2.6 and 1.7 μ M, respectively.¹ In insulin-resistant mice, 100 mg/kg/day amorfrutin B is reported to improve insulin sensitivity, glucose tolerance, and plasma lipid concentrations after two weeks of treatment.¹

Reference

1. Weidner, C., Wowro, S.J., Freiwald, A., *et al.* Amorfrutin B is an efficient natural peroxisome proliferator-activated receptor gamma (PPAR γ) agonist with potent glucose-lowering properties. *Diabetologia* **56**(8), 1802-1812 (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, 11/08/2022

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