

Genipin

UCP2 inhibitor

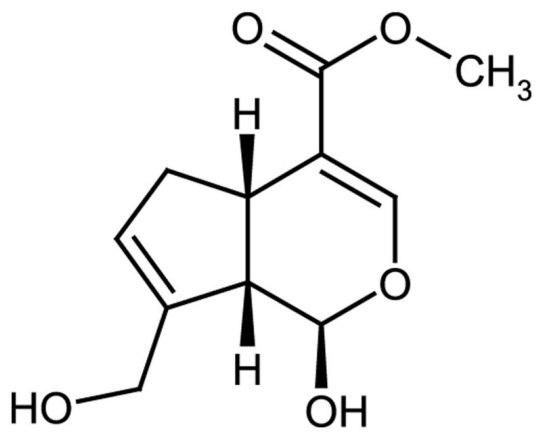
Cell permeable inhibitor of uncoupling protein 2 (UCP2). Increases glucose-stimulated insulin secretion, mitochondrial membrane potential and ATP levels in pancreatic island cells. Induces apoptosis. Protein cross-linking agent. Anti-inflammatory and anti-angiogenic.

Citations: 5

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (+)-Genipin, Methyl (1S,2R,6S)-2-hydroxy-9-(hydroxymethyl)-3-oxabicyclo[4.3.0]nona-4,8-diene-5-carboxylate

Appearance White to off-white crystalline solid.

CAS 6902-77-8

Couple Target UCP

Couple Type Inhibitor

Formula $C_{11}H_{14}O_5$

MW 226.2

Purity ≥98% (HPLC)

RTECS GY5828000

Solubility Soluble in DMSO (50mg/ml), dimethyl formamide (25mg/ml), 100% ethanol (5mg/ml), methanol or ethyl acetate; insoluble in water.

Source Isolated from *Gardenia jasminoides* Ellis.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com