

Puerarin

5-HT_{2c} inhibitor

Biologically active isoflavone. Affects serotonin levels and platelet aggregation in blood cells. 5-HT_{2c} antagonist. Antibacterial. Displays cardioprotective effects. Induces apoptosis.

Citations: 8

[View Online »](#)

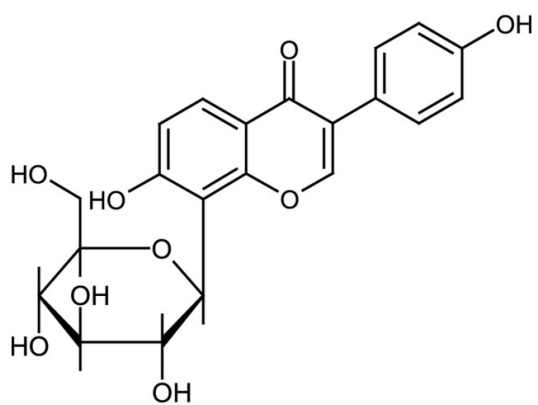
Ordering Information

[Order Online »](#)

ALX-350-249-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 8-(β-D-Glucopyranosyl-7-hydroxy-3- (4-hydroxyphenyl)-4H-1-benzopyran-4-one, NPI-031G

Appearance White to off-white powder.

CAS 3681-99-0

Couple Target Serotonin receptor

Couple Type Inhibitor, Ligand

Formula $C_{21}H_{20}O_9$

MW 416.4

Purity ≥98% (HPLC)

RTECS UO5216000

Solubility Soluble in methanol.

Source Isolated from *Kudzu* root.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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