

(-)-Arctigenin

Antitumor agent

Lignan derivative, which shows antitumor, anti-inflammatory, immunomodulatory and neuroprotective activities. Down-regulates anti-apoptotic protein Bcl-X_L. Potent inhibitor of HIV type-I integrase and DNA topoisomerase II. Shows phytoestrogenic and cytotoxic properties against Hep62 cells. Antagonist for platelet activating factor and Ca²⁺. Blocks the activation of Akt (protein kinase B; PKB) induced by glucose starvation.

Citations: 9

[View Online »](#)

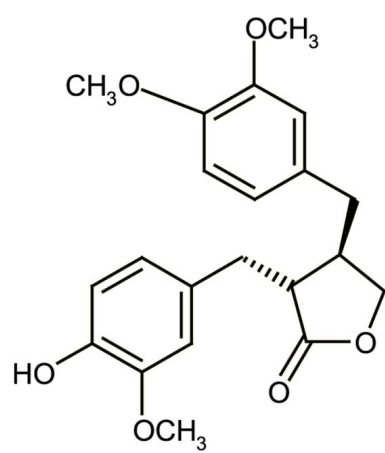
Ordering Information

[Order Online »](#)

ALX-350-312-M025	25mg
------------------	------

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 (3R,4R)-4-[(3,4-Dimethoxyphenyl)methyl]dihydro-3-[(4-hydroxy-3-methoxyphenyl)methyl]-2-(3H)-furanone

Appearance Colorless crystals.

CAS 7770-78-7

Couple Target HIV, Topoisomerase

Couple Type Inhibitor

Formula $C_{21}H_{24}O_6$

Identity Determined by NMR and MS.

MW 372.4

Purity ≥97% (HPLC)

Solubility Soluble in DMSO (34mg/ml, heating, sonication); sparingly soluble in water.

Source Isolated from *Arctium lappa*.

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-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

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

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