

(-)-Arctiin

Antitumor agent

Lignan derivative with antitumor activities. Down-regulates anti-apoptotic protein Bcl-X_L. Potent inhibitor of HIV type-I integrase and DNA topoisomerase II. Shows phytoestrogenic activity and is an antagonist for platelet activating factor and Ca²⁺.

Citations: 8

[View Online »](#)

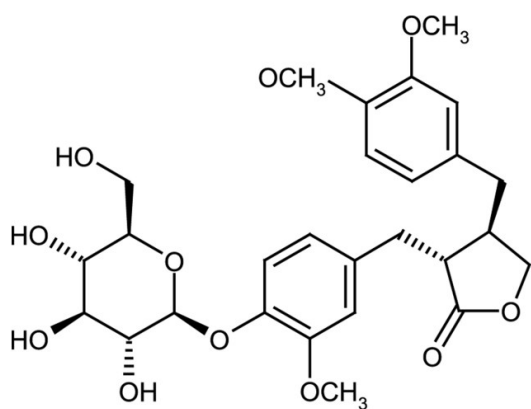
Ordering Information

[Order Online »](#)

ALX-350-318-M025	25mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
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]-3-[[4-(β-D-glucopyranosyloxy)-3-methoxyphenyl]methyl]dihydro-2(3H)-furanone
Appearance	White to off-white solid.
CAS	20362-31-6
Couple Target	HIV, Topoisomerase
Couple Type	Inhibitor
Formula	$C_{27}H_{34}O_{11}$
MW	534.6
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (5mg/ml); sparingly soluble in water.
Source	Isolated from <i>Arctium lappa</i> .

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-eu@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-
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)