

Magnolol

Antitumor and anti-angiogenic agent

Potent and highly tolerable antitumor and antiangiogenic compound. Anxiolytic, anti-thrombotic and antibacterial. Shows central depressant effects *in vivo*. Inhibitor of nitric oxide (NO) and TNF- α production in LPS-activated macrophages by the suppression of inducible nitric oxide synthase (iNOS; NOS II) expression ($IC_{50}=16.8\mu M$). Isomeric with honokiol (Prod. No. ALX-350-350).

Citations: 6

[View Online »](#)

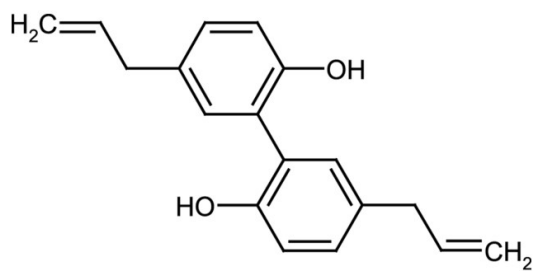
Ordering Information

[Order Online »](#)

ALX-350-352-M005	5mg
ALX-350-352-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	5,5'-Diallyl-2,2'-biphenyldiol
Appearance	White to light yellow powder.
CAS	528-43-8
Couple Target	NOS
Couple Type	Inhibitor
Formula	$C_{18}H_{18}O_2$
MI	14: 5697
MW	266.3
Purity	≥98% (HPLC)
RTECS	DV5105500
Solubility	Soluble in DMSO, 100% ethanol or methanol.
Source	Isolated from <i>Magnolia officinalis</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-
be@enzolifesciences.com](mailto:info-be@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)