

# Honokiol

## iNOS inhibitor

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- $\alpha$  production in LPS-activated macrophages by the suppression of inducible nitric oxide synthase (iNOS; NOS II) expression ( $IC_{50}$ =6.4 $\mu$ M). Isomeric with magnolol (Prod. No. ALX-350-352). Inhibits NF- $\kappa$ B activation. Potent scavenger of superoxide and peroxy radicals.

Citations: 9

[View Online »](#)

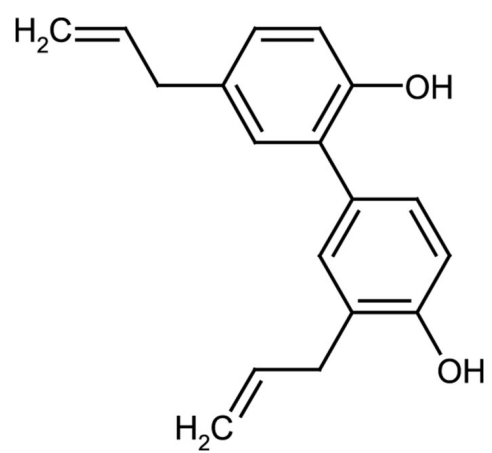
## Ordering Information

[Order Online »](#)

ALX-350-350-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	5,3'-Diallyl-2,4'-biphenyldiol, 5,3'-Diallyl-2,4'-dihydroxydiphenyl
Appearance	White to off-white powder.
CAS	35354-74-6
Couple Target	NOS
Couple Type	Inhibitor
Formula	$C_{18}H_{18}O_2$
MI	14: 4742
MW	266.3
Purity	≥98% (HPLC)
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-  
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)