

# Radicicol

## HSP90 inhibitor

Antifungal macrocyclic lactone antibiotic with antimalarial activity. Potent inhibitor of HSP90. Binds more strongly to HSP90 (nanomolar affinity) than to Grp94. Also binds to yeast HSP90, *E. coli* HtpG and TRAP-1. Non-competitive inhibitor of ATP citrate lyase. Anti-angiogenic. Specifically inhibits the interaction between HIF-1 $\alpha$ /Arnt heterodimer and the hypoxia-responsive element (HRE), reducing VEGF expression. Protein tyrosine kinase inhibitor, e.g. p60<sup>V-src</sup> (IC<sub>50</sub>=0.27nM). Inhibitor of cyclooxygenase-2 (COX-2) expression (IC<sub>50</sub>=27nM) without affecting COX-1 expression in LPS-stimulated macrophages. Induces the differentiation of HL-60 cells into macrophages, blocking cell cycle at G1 and G2. Suppressor of NIH 3T3 cell transformation by diverse oncogenes such as *mos*, *ras* and *src* in part by blocking the key signal transduction intermediates such as MAP kinase and GAP-associated p62. Inhibitor of AP-1-, NF- $\kappa$ B- and serum response factor (SRF)-mediated transcription (e.g. expression of iNOS). Represses the transcriptional function of the estrogen receptor. Inhibits archeal growth and DNA topoisomerase VI (a Topo IIB family topoisomerase). Blocks replication of negative-strand RNA viruses. Causes morphological reversion of ras-transformed cells and leads to selective depletion of Raf kinase. Displays anti-angiogenic activity. Binds to and inhibits Hsp 90 molecular chaperone.

Citations: 28

[View Online »](#)

## Ordering Information

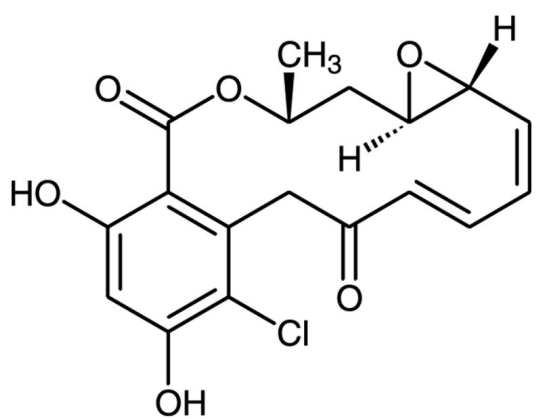
[Order Online »](#)

BML-EI285-0001

1mg

## 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	Monorden, 1aS-(1aR*,2Z,4E,14*,15aR*)]-8-Chloro-1a,14,15,15a-tetrahydro-9,11-dihydroxy-14-methyl-6H-oxireno[e][2]benzoxacyclotetradecin-6,12(7H)-dione
Appearance	White to off-white or yellowish solid. May appear adhered to vial.
CAS	12772-57-5
Couple Target	HSP90
Couple Type	Inhibitor
Formula	$C_{18}H_{17}ClO_6$
MI	14: 6253
MW	364.8
Purity	≥97% (HPLC)
RTECS	RR1105000
Solubility	Soluble in DMSO, methanol or acetone.
Source	Isolated from <i>Humicola fuscoatra</i> .



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)