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. Noncompetitive inhibitor of ATP citrate lyase. Anti-angiogenic. Specifically inhibits the interaction between HIF-1α/Arnt heterodimer and the hypoxiaresponsive element (HRE), reducing VEGF expression. Protein tyrosine kinase inhibitor, e.g. $p60^{V-}$ src (IC₅₀=0.27nM). Inhibitor of cyclooxygenase-2 (COX-2) expression (IC $_{50}$ =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-κB- 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 »

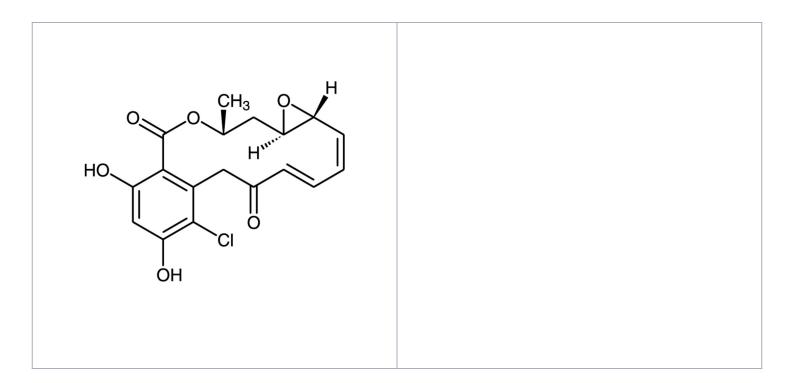
Ordering Information

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

Monorden, 1aS-(1aR*,2Z,4E,14*,15aR*)]-8-Chloro-**Alternative Name**

> 1a,14,15,15a-tetrahydro-9,11-dihydroxy-14-methyl-6Hoxireno[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 C₁₈H₁₇CIO₆

ΜI 14: 6253

MW 364.8

Purity ≥97% (HPLC)

RTECS RR1105000

Solubility Soluble in DMSO, methanol or acetone.

Source



Last modified: May 29, 2024

European Sales Office ENZO LIFE SCIENCES (ELS) AG

Phone: +41 61 926 8989 eu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 fr@enzolifesciences.com

Isolated from Humicola fuscoatra.

Phone: +49 7621 5500 526 de@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 uk@enzolifesciences.com