

# Picropodophyllin

## Insulin receptor inhibitor

A potent and selective inhibitor of insulin-like growth factor 1 receptor (IGF-1R,  $IC_{50}=6$  nM). It efficiently blocks IGF-1R activity, reduces pAKT and pERK1/2, induces apoptosis in IGF-1R-positive tumor cells and causes complete tumor regression in xenografted and allografted mice.

Picropodophyllin downregulates IGF-1R by interfering with the action of  $\beta$ -arrestin 1/MDM2. It causes tumor regression and attenuates invasiveness of uveal melanoma cells. It displays beneficial effects on tumor growth, angiogenesis, bone disease and survival in a mouse multiple myeloma model. It attenuates intimal hyperplasia after vascular injury. Inhibition of IGF-1R is non-competitive with ATP.

Citations: 8

[View Online »](#)

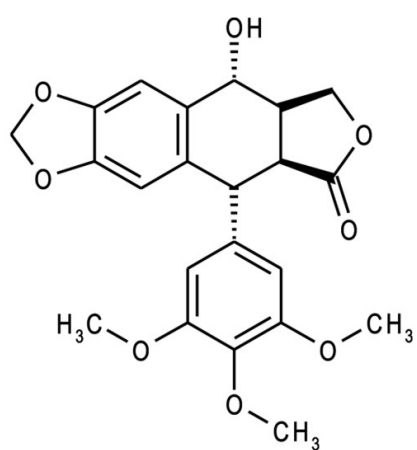
## Ordering Information

[Order Online »](#)

|                |     |
|----------------|-----|
| BML-EI372-0001 | 1mg |
|----------------|-----|

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Appearance** White solid

**CAS** 477-47-4

**Couple Target** IGF receptor

**Couple Type** Inhibitor

**Formula**  $C_{22}H_{22}O_8$

**MW** 414.4

**Purity** ≥98% (HPLC)

**Solubility** Soluble in DMSO (10mg/ml, warm) or dimethyl formamide (25mg/ml)

**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)