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 β -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

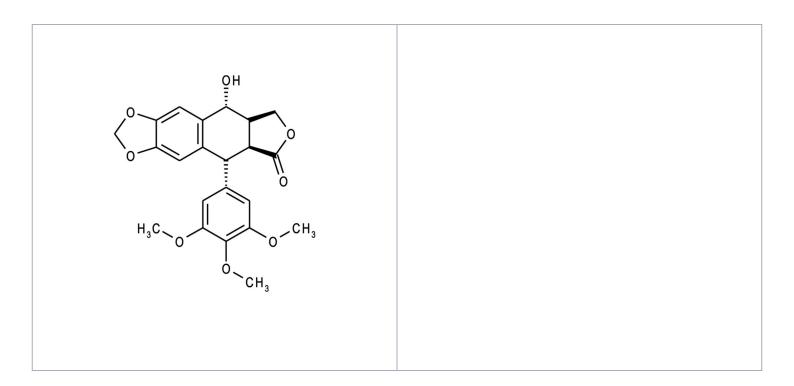
Order Online »

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

