# Betulinic acid (high purity)

**Antitumor and anti-HIV agent** 

Antitumor and anti-HIV agent. Induces apoptosis by activating mitochondrial permeability transition (MPT). Inhibits NF- $\kappa$ B activation and NF- $\kappa$ B-regulated gene expression induced by carcinogens and inflammatory stimuli. Decreases expression of Bcl-2 and cyclin D1. Potent proteasome activator. Triggers CD95- and p53-independent apoptosis via activation of caspases. It induces mitochondrial permeability pore opening and the release of cytochrome c and apoptosis-inducing factor. Displays anti-HIV activity. EC<sub>50</sub> for inhibion of HIV-1 replicaton in H9 lymphocytes: 1.4 $\mu$ M.

Citations: 22

View Online »

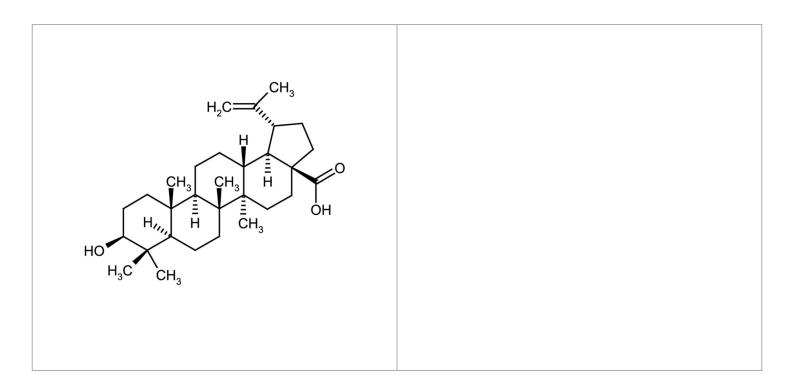
## **Ordering Information**

**Order Online** »

ALX-350-277-M00	5mg	
ALX-350-277-M02	25mg	

Manuals, SDS & CofA

**View Online** »



### **Handling & Storage**

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

**Handling** Keep cool and dry.

Long Term Storage +4°C

**Shipping** Ambient Temperature

#### Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name3β-Hydroxy-20(29)-lupaene-28-oic acid

**Appearance** White to off-white powder.

**CAS** 472-15-1

Formula  $C_{30}H_{48}O_3$ 

**MW** 456.7

Purity ≥97% (HPLC)

**Solubility** Soluble in DMSO.

Source Isolated from *Platanus acerifolia* (plane) tree bark.

Last modified: May 29, 2024