# **Anacardic acid**

Cell permeable salicylic acid analog that acts as a potent, non-competitive inhibitor of p300 and PCAF (p300/CBP-associated factor) histone acetyltransferase (HAT) activities (IC $_{50}$ ~8.5µM and ~5µM, respectively). Displays a variety of other biological activities, such as antibacterial, antimicrobial, prostaglandin synthase inhibition, tyrosinase, and lipoxygenase inhibition.Inhibits SUMOylation of RanGAP1-C2 with an IC $_{50}$  = 2.2µM

Citations: 14

View Online »

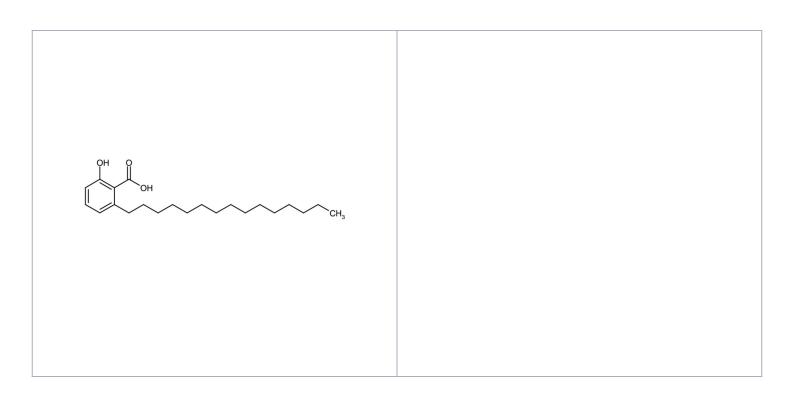
## **Ordering Information**

**Order Online** »

ALX-270-381-M005	5mg
ALX-270-381-M025	25mg

Manuals, SDS & CofA

**View Online** »



### **Handling & Storage**

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

stable for up to 3 months when stored at -20°C.

**Handling** Protect from light. Packaged under inert gas. After reconstitution, prepare aliquots and

store at -20°C.

Long Term Storage -20°C

**Shipping** Ambient Temperature

#### Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name 2-Hydroxy-6-pentadecylbenzoic acid, 6-Pentadecylsalicylic

acid, AA

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

CAS 16611-84-0

Couple Target HAT, Lipoxygenase, Tyrosinase

Couple Type Inhibitor

Formula  $C_{22}H_{36}O_3$ 

MI 14: 621

MW 348.5

Purity ≥95% (HPLC)

**Soluble** in 100% ethanol, methanol, DMSO,

dichloromethane or ethylacetate.

**Source** Synthetic.



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 info-

eu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

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