

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} ~8.5 μ M and ~5 μ M, 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} = 2.2 μ M

Citations: 14

[View Online »](#)

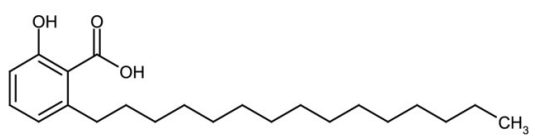
Ordering Information

[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)
Solubility	Soluble in 100% ethanol, methanol, DMSO, dichloromethane or ethylacetate.
Source	Synthetic.



Last modified: May 29, 2024

ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@enzolifesciences.com

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

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

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

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