
Anacardic Acid

(Catalog No. M37000)

Alternative Names

AA, 2-Hydroxy-6-pentadecylbenzoic Acid, 6-Pentadecylsalicylic Acid

Description

Anacardic Acid is a pentadecane aliphatic chain containing hydroxylcarboxylic acid. The acid is an effective inhibitor of the activities of p300 and p300/CBP associated factor histone acetyltransferase. Additionally, it has been shown to potently inhibit HAT-dependent transcription and SUMOylation, as well as decrease expression of NF- κ B regulated gene products including, invasion, proliferation, inflammation, and cell survival. Research suggests that anacardic acid may be a potent activator of ARK-1 (Aurora kinase A) mediated phosphorylation of histone H3, and a novel activator of Aurora kinase. Anacardic Acid has demonstrated bactericidal activity against methicillin-resistant *Staphylococcus aureus* (MRSA).

CAS Number

16611-84-0

Chemical Formula

$C_{22}H_{36}O_3$

Molecular Weight

348.52

Size

5 mg

Purity

$\geq 95\%$

Solubility

Soluble in 100% ethanol, methanol, DMSO, dichloromethane or ethylacetate

Storage Temp

-20°C

Ordering Information:

| Products | Size | Cat. No. |
|----------------|------|-----------|
| Anacardic Acid | 5 mg | M37000 -1 |

This product is for research purposes only. Not intended for use in diagnostic procedures.