

Garcinol

(Catalog No. M35001)

Alternative Names

Camboginol

Description

Garcinol is a polyisoprenylated benzophenone derivative from the rind of *Garcinia indica* fruit and is a potent inhibitor of histone acetyltransferases p300 (IC₅₀= ~7 μm) and PCAF (IC₅₀= ~5 μm) both in vitro and in vivo. p300, also known as E1A binding protein p300, regulates cell growth and division through interacting with numerous transcription factors, and the p53 associated transcriptional coactivator PCAF (P300/CBP-associated factor) plays a direct role in transcriptional regulation. Reports show that Garcinol represses the p300-mediated acetylation of p53 in vivo, promotes apoptosis through caspase-3 activation and predominantly down-regulates global gene expression in HeLa cells. Researchers found that Garcinol exhibits dose-dependent cancer cell-specific growth inhibition in ER-positive MCF-7 and ER-negative MDA-MB-231 breast cancer cells. Additional research shows that Garcinol exhibits anti-carcinogenic and anti-inflammatory properties through selectively suppressing Prostaglandin E2 synthesis and 5-lipoxygenase product formation.

CAS Number

78824-30-3

Chemical Formula

C₃₈H₅₀O₆

Molecular Weight

602.8

Size

10 mg

Purity

95%

Solubility

Soluble in DMSO (25 mg/ml), Ethanol (25 mg/ml), DMF (25 mg/ml), Water (nil)

Storage Temp

-20°C

Ordering Information:

Products	Size	Cat. No.
Garcinol	10 mg	M35001 -1

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