

Garcinol

Epigenetic Enzyme Inhibitor

Catalog # K311-903C

Lot # P3320-18

CAS # 78824-30-3

Product Description

Molecular Formula: C₃₈H₅₀O₆

Appearance: Yellow powder

Molecular Weight: 602.8

Purity: >97% (TLC); NMR (Conforms)

Solubilization: May be dissolved in DMSO (25 mg/ml); or ethanol (25 mg/ml)

Alias

Camboginol, 3-(3,4-dihydroxybenzoyl)-4-hydroxy-8,8-dimethyl-1,7-bis(3-methyl-2-buten-1-yl)-5-[(2S)-5-methyl-2-(1-methylethenyl)-4-hexen-1-yl]-bicyclo[3.3.1]non-3-ene-2,9-dione

Specific Activity

HAT Inhibitor

Storage and Stability

Store desiccated as supplied at -20°C for up to 2 years.

Store solutions at -20°C for up to 2 months.

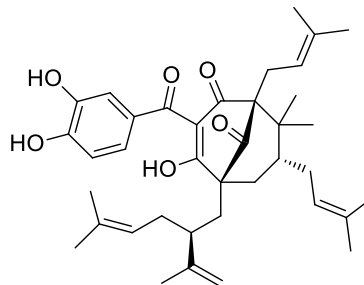
Scientific Background

Potent inhibitor of histone acetyl transferases p300 (IC₅₀=7 μM) and pCAF (IC₅₀=5 μM)¹. Induces apoptosis in various cell lines². Promotes expansion of human hematopoietic stem cells³. Cell permeable.

References

1. K Balasubramanyam et al. J. Biol. Chem. 2004 279:33716
2. MA Parasramka et al. Nutr. Cancer 2011 63:456
3. T Nishino et al. Nutr. PLoS One 2011 6:e24298

Molecular Structure



Garcinol

Epigenetic Enzyme Inhibitor

Catalog # K311-903C

CAS # 78824-30-3

Lot # P3320-18

Purity >97%

Format Yellow powder

Solubilization May be dissolved in DMSO (25 mg/ml); or ethanol (25 mg/ml)

Storage & Shipping Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 2 months.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.