Garcinol

Catalog No: 14076, 14077 Format: 50 mg, 10 mg

Chemical Properties:

$$\label{eq:main_state} \begin{split} &\mathsf{MW} = 602.8\\ &\mathsf{C}_{38}\mathsf{H}_{50}\mathsf{O}_6\\ &\mathsf{CAS}\ 78824\text{-}30\text{-}3\\ &\mathsf{Physical}\ \mathsf{Properties:}\ \ \mathsf{Yellow}\ \mathsf{solid}\\ &\mathsf{Names:}\ \mathsf{Garcinol} \end{split}$$

Pharmacology: Potent inhibitor of histone acetyltransferases p300 (IC_{50} = 7 M) and pCAF (IC_{50} = 5 M)(ref 1). Induces apoptosis in various cell lines (ref 2). Promotes expansion of human hematopoietic stem cells (ref 3). Cell permeable.

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

Fluorescent Properties: N/A

Quality Control: >97% (TLC); NMR (Conforms)

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. PLoS One 2011, 6:e24298

Storage and Guarantee: Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 2 months. This product is guaranteed for 6 months from date of arrival.



