Sirtinol



Catalog No: 14074, 14075 Format: 25 mg, 5 mg

Chemical Properties:

$$\begin{split} &\mathsf{MW} = 394.5\\ &\mathsf{C}_{26}\mathsf{H}_{22}\mathsf{N}_2\mathsf{O}_2\\ &\mathsf{CAS} \ 410536\text{-}97\text{-}9\\ &\mathsf{Physical Properties:} \ \ \mathsf{Yellow \ powder}\\ &\mathbf{Names:} \ \mathsf{Sirtinol}, \ 2\text{-}[(2\text{-}Hydroxynaphthalen\text{-}1\text{-}ylmethylene)amino]\text{-}N\text{-}(1\text{-}phenethyl)benzamide} \end{split}$$

Pharmacology: Selective sirtuin inhibitor which does not effect HDAC1 (IC_{50} = 131 M, 38 M and 68 M for SIRT1, SIRT2 and Sir2P respectively) (ref 1). Induces apoptotic and autophagic cell death in MCF-7 human breast cancer cell lines (ref 2). Suppresses survivin expression which correlates with decreased viability in HT29 colon cancer cells (ref 3). Displays pro-inflammatory effects in adipocytes (ref 4).

Solubilization: May be dissolved in DMSO (35 mg/ml); Ethanol (10 mg/ml)

Fluorescent Properties: N/A

Quality Control: >98% (TLC); NMR (Conforms)

References:

- 1. CM Grosinger et al. J. Biol. Chem. 2001, 276:38837
- 2. J Wang et al. Int. J. Oncol. 2012, 41:1101
- 3. YF Hsu et al. Biochim. Biophys. Acta 2012, 1820:104
- 4. QQ Lin et al. Cytokine 2012, 60:447

Storage and Guarantee: Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 1 month. This product is guaranteed for 6 months from date of arrival.

