

# Cambinol

(Catalog No. M47000)

## Alternative Names

5-[(2-hydroxy-5,8-dihydronaphthalen-1-yl)methyl]-6-phenyl-2-thioxo-2,3-dihy.

## Description

Cambinol is a cell-permeable  $\beta$ -naphthol compound that inhibits the NAD-dependent deacetylase activity of hSIRT1 and hSIRT2 (IC<sub>50</sub> = 56  $\mu$ M and 59  $\mu$ M, respectively) in a substrate-, but not NAD-, competitive manner. It inhibits SIRT5 deacetylase activity only at much higher concentrations (IC<sub>50</sub> >300  $\mu$ M) and exhibits no inhibition against class I and II HDACs. Cambinol-induced apoptosis in BCL6-expressing Burkitt lymphoma cell lines has been attributed to the hyperacetylation of BCL6 and p53. Shown to effectively inhibit xenografted Daudi Burkitt lymphoma growth in mice in vivo.

## CAS Number

14513-15-6

## Chemical Formula

C<sub>21</sub>H<sub>16</sub>N<sub>2</sub>O<sub>2</sub>S

## Molecular Weight

360.4

## Size

5 mg

## Purity

>95%

## Solubility

Soluble in DMSO or EtOH

## Storage Temp

-20°C

## Ordering Information:

### Products

Cambinol

### Size

5 mg

### Cat. No.

M47000 -1

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