

## Sirtinol

Epigenetic Enzyme Inhibitor

### Catalog # S35-904B

Lot # P3320-35

CAS # 410536-97-9

### Product Description

Molecular Formula: C<sub>26</sub>H<sub>22</sub>N<sub>2</sub>O<sub>2</sub>

Appearance: Yellow powder

Molecular Weight: 394.5

Purity: >98% (TLC); NMR (Conforms)

Solubilization: May be dissolved in DMSO (35 mg/ml) or ethanol (10 mg /ml)

Storage: Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 1 month.

### Alias

2-[[[2-Hydroxy-1-naphthalenyl)methylene]amino]-N-(1-phenylethyl)benzamide

### Specific Activity

SIRT Inhibitor

### Storage and Stability

Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 1 month.

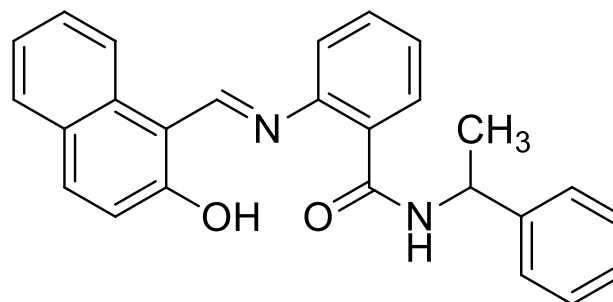
### Scientific Background

Selective sirtuin inhibitor which does not effect HDAC1, IC<sub>50</sub> = 131, 38 and 68 μM for SIRT1, SIRT2 and Sir2P respectively<sup>1</sup>. Induces apoptotic and autophagic cell death in MCF-7 human breast cancer cells<sup>2</sup>. Suppresses survivin expression which correlates with decreased viability in HT29 colon cancer cells<sup>3</sup>. Displays pro-inflammatory effects in adipocytes<sup>4</sup>.

### 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

### Molecular Structure



## Sirtinol

Epigenetic Enzyme Inhibitor

Catalog #	S35-904B
CAS #	410536-97-9
Lot #	P3320-35
Purity	>98% (TLC)
Format	Yellow powder
Solubilization	May be dissolved in DMSO (35 mg/ml) or ethanol (10 mg /ml)
Storage & Shipping	Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 1 month.

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

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