

Splitomicin

Epigenetic Enzyme Inhibitor

Catalog # S35-904G

Lot # P3320-36

CAS # 5690-03-09

Product Description

Molecular Formula: C₁₃H₁₀O₂

Appearance: White powder

Molecular Weight: 198.2

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

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

Alias

1,2-Dihydro-3H-naphtho[2,1-b]pyran-3-one

Specific Activity

Sirt1 Inhibitor

Storage and Stability

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

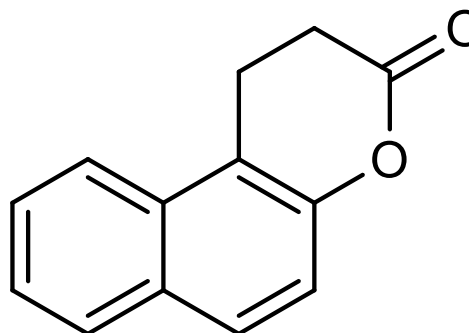
Scientific Background

Initially shown to inhibit Sir2p (the yeast Sirt1 homolog) and create a conditional phenocopy of a sir2 deletion mutant in *S. cerevisiae*¹. Cells grown in the presence of the drug have silencing defects at telomeres, silent mating-type loci and the ribosomal DNA¹. Inhibits all class III HDACs (sirtuins) and alleviates gene silencing in Fragile X mental retardation syndrome². Reverses ischemic preconditioning-induced cardioprotection in a mouse model³.

References

1. A Bedalov et al. Proc. Natl. Acad. Sci. USA 2001 98:15113
2. R Biacsi et al. PLoS Genet. 2008 4(3):e1000017
3. SM Nadtochiy et al. Cardiovasc. Res. 2011 89:643

Molecular Structure



Splitomicin

Epigenetic Enzyme Inhibitor

Catalog #	S35-904G
CAS #	5690-03-09
Lot #	P3320-36
Purity	>98% (TLC)
Format	White powder
Solubilization	May be dissolved in DMSO (>25 mg/ml) or ethanol (>25 mg/ml)
Storage & Shipping	Store desiccated as supplied at ambient temperature 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
www.signalchem.com

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