

---

## Splitomicin

(Catalog No. M45001)

---

### Alternative Names

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

### Description

Splitomicin is a cell-permeable lactone that acts as a selective inhibitor of NAD<sup>+</sup> dependent histone deacetylase activity of Sir2 protein. It creates a conditional phenocopy of a sir2 deletion mutant in *S. cerevisiae* and sensitizes mammalian cells to a variety of DNA-damaging agents by abrogating Sir2p activity on p53. Acts by either altering or blocking access to the acetylated histone binding pocket.

### CAS Number

5690-03-9

### Chemical Formula

C<sub>13</sub>H<sub>10</sub>O<sub>2</sub>

### Molecular Weight

198.2

### Size

5 mg

### Purity

≥98%

### Solubility

DMSO

### Storage Temp

-20°C

### Ordering Information:

Products	Size	Cat. No.
Splitomicin	5 mg	M45001 -1

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