Antimycin A

Antibiotic

Antimycin A is a bacterial metabolite complex produced by *Streptomyces*. It is composed majorly of Antimycin A1, A2, A3, and A4. Antimycin A is an antibiotic. It is capable of inducing apoptosis in mammalian cells, which is not prevented by the presence of Bcl-2. Antimycin A acts as an inhibitor of cellular respiration and more specifically, the electron transport chain of oxidative phosphorylation via its binding to cytochrome c reductase.

Citations: 29

View Online »

Ordering Information

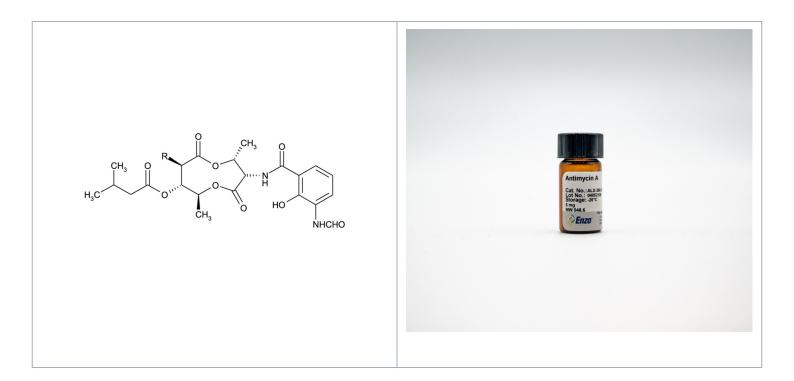
Order Online »

ALX-380-075-M005	5mg
ALX-380-075-M010	10mg

Manuals, SDS & CofA

View Online »

- Antimycin A complex composed mostly of Antimycin A1, A2, A3, and A4
- Used for the study of the mitochondrial function



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Protect from light.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Antipiricullin, Virosin

Appearance White to faint yellow powder.

CAS 1397-94-0

Identity Identity determined by NMR.

MI 14: 714 (A1) & 715 (A3)

RTECS CD0350000

Solubility Soluble in 100% ethanol (50mg/ml).

Source Isolated from *Streptomyces sp.* Mixture of antimycin A's.

Last modified: May 29, 2024

