

# 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

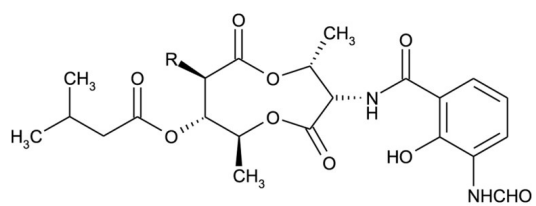
[Order Online »](#)

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	<b>14:</b> 714 (A1) & 715 (A3)
RTECS	CD0350000
Solubility	Soluble in 100% ethanol (50mg/ml).
Source	Isolated from <i>Streptomyces</i> sp. Mixture of antimycin A's.



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)