

# Oligomycin

F-ATPase inhibitor

Macrolide antibiotic that inhibits membrane bound mitochondrial ATPase (F<sub>1</sub>), preventing phosphoryl group transfer. Induces apoptosis.

Citations: 16

[View Online »](#)

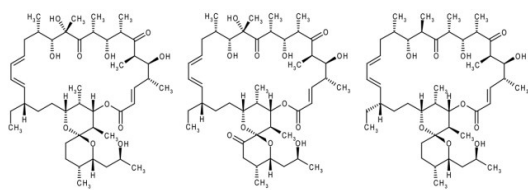
## Ordering Information

[Order Online »](#)

|                  |      |
|------------------|------|
| ALX-380-037-M005 | 5mg  |
| ALX-380-037-M010 | 10mg |

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Long Term Storage** +4°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Appearance** White to faint yellow powder.

**CAS** 1404-19-9

**Couple Target** ATPase

**Couple Type** Inhibitor

**MI** **14:** 6833

**MW** 791.1 (Oligomycin A)

**Purity** ≥90% (HPLC, TLC)

**Purity Detail** Mixture of oligomycin A (major component), B, and C.

**RTECS** RK3325000

**Solubility** Soluble in 100% ethanol, methanol (10mg/ml) or acetone (50mg/ml).

**Source** Isolated from *Streptomyces diastatochromogens*.

**Last modified:** May 29, 2024



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)