

# Rapamycin

An mTOR inhibitor that modulates cell growth, proliferation, and autophagy - frequently used in signaling pathway research and immunosuppressive studies.

Rapamycin is a macrocyclic-triene antibiotic with potent immunosuppressant activity. Rapamycin forms a complex with FKBP12 that binds to an effector, thus inhibiting IL-2 and other growth promoting lymphokines. The effectors were identified as FRAP (FKBP12 rapamycin-associated protein) and RAFT1 (rapamycin and FKBP12 target). Rapamycin/FKBP complex does not inhibit FRAP PI 4-kinase activity, but does inhibit FRAP autophosphorylation. Rapamycin induces inhibition of p70<sup>S6k</sup>, p33<sup>cdk2</sup> and p34<sup>cdc2</sup>. Rapamycin selectively blocks signaling leading to the activation of p70/85 S6 kinase. Rapamycin complexed with FKBP12 binds to and inhibits the activity of the mammalian target of rapamycin (mTOR) complex 1 (mTORC1). Rapamycin can inhibit cell growth, enhance apoptosis, and activates autophagy in a variety of cell types.

Citations: 127

[View Online »](#)

## Ordering Information

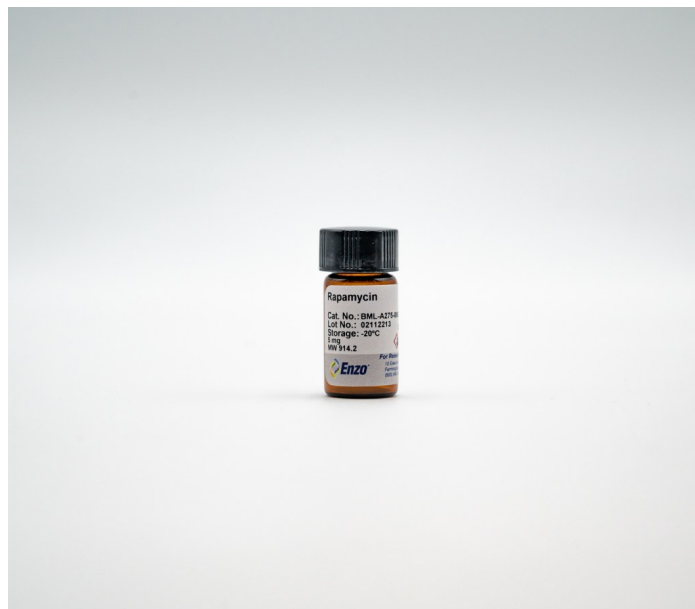
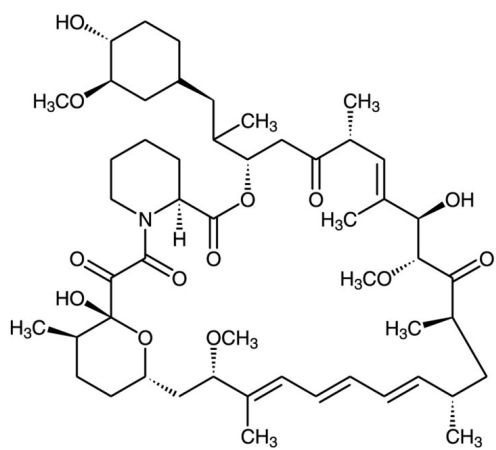
[Order Online »](#)

BML-A275-0005	5mg
BML-A275-0025	25mg

## Manuals, SDS & CofA

[View Online »](#)

- Antibiotic with immunosuppressant and anti-cancer properties
- Commonly used to study autophagy in a variety of cells
- Highly cited



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Sirolimus, AY-22,989, RAPA, Rapamune
Appearance	Off-white or yellow solid.
CAS	53123-88-9
Couple Target	Immunophilin, mTOR
Couple Type	Inhibitor, Ligand
Formula	$C_{51}H_{79}NO_{13}$
MI	14: 8114
MW	914.2
Purity	≥98% (HPLC)
RTECS	VE6250000
Solubility	Soluble in chloroform, DMSO (25mg/ml), or 100% methanol.
Source	Isolated from <i>Streptomyces hygroscopus</i>

## Technical Info / Product Notes

Replacement for **ADI-HPK-109**.



Last modified: May 29, 2024

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

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