

Rapamycin

Immunosuppressant

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^{S6k}, p33^{cdk2} and p34^{cdc2}. 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: 126

[View Online »](#)

Ordering Information

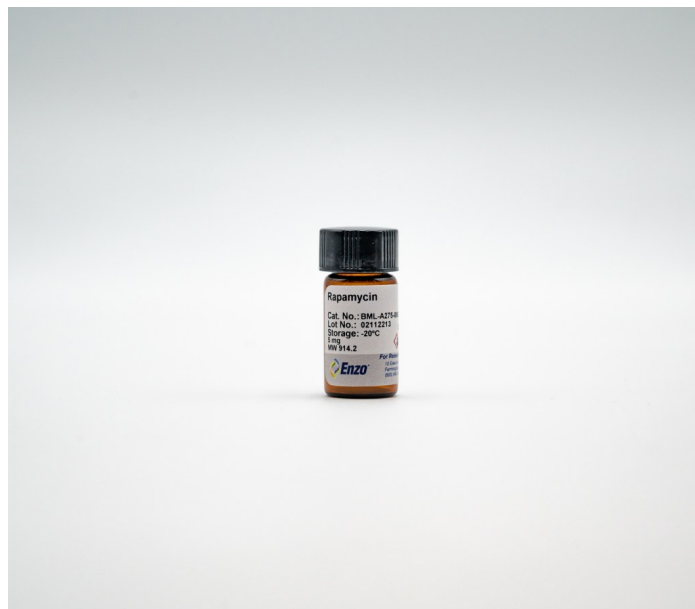
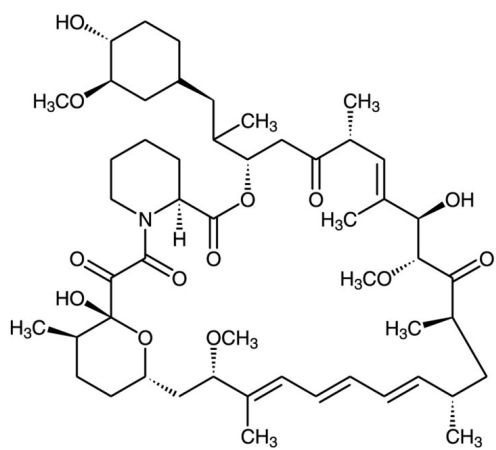
[Order Online »](#)

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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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