

# Rapamycin

Macrolide originally produced by *Streptomyces*; mTOR inhibitor.

An antifungal antibiotic that has immunosuppressive activity. It blocks cytokine-mediated signal transduction pathways.

## Ordering Information

[Order Online »](#)

LKT-R0161-M001	1mg
LKT-R0161-M010	10mg
LKT-R0161-M025	25mg
LKT-R0161-M100	100mg

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Protect from light. Hygroscopic.

**Long Term Storage** +4°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Sirolimus, AY-22989

**CAS** 53123-88-9

**Couple Type** Inhibitor

**Formula**  $C_{51}H_{79}NO_{13}$

**MI** 13: 8202

**MW** 914.17

**Purity** ≥99%

**Solubility** Soluble in ether, chloroform, acetone, methanol, DMF, DMSO (20 mg/mL), and ethanol (up to 2 mM). Sparingly soluble in hexane and petr ether. Insoluble in water.



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-be@enzolifesciences.com](mailto:info-be@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)