

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-M010	10mg
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

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@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