

# FK506

## Immunosuppressant

Immunosuppressant that blocks T cell proliferation *in vitro* by inhibiting the generation of several lymphokines, especially IL-2. Shown to inhibit the activity of FK506 binding protein and to inhibit the  $\text{Ca}^{2+}$ -dependent phosphatase and protein phosphatase 2B (PP2B; calcineurin). Activates NF- $\kappa$ B through phosphorylation and degradation of I $\kappa$ B $\alpha$ .

Citations: 27

[View Online »](#)

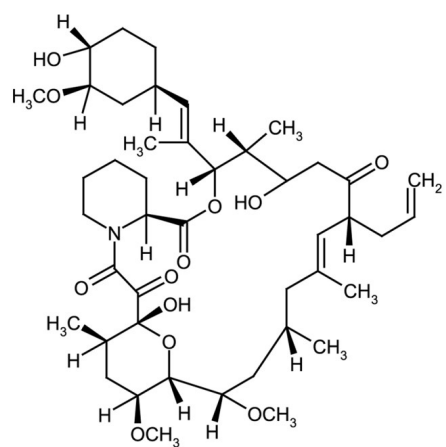
## Ordering Information

[Order Online »](#)

ALX-380-008-M001	1mg
ALX-380-008-M005	5mg
ALX-380-008-M025	25mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Handling** Protect from light.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Tacrolimus

**Appearance** White to off-white crystalline powder.

**CAS** 104987-11-3

**Couple Target** FKBP, Serine/threonine-protein phosphatase

**Couple Type** Inhibitor, Ligand

**Formula**  $C_{44}H_{69}NO_{12}$

**Identity** FT-IR

**MI** 14: 9025

**MW** 804.0

**Purity** ≥98% (HPLC)

**RTECS** KD4200000

**Solubility** Soluble in DMSO, 100% ethanol, methanol, ethyl acetate or acetone; insoluble in water.

**Source** Isolated from *Streptomyces tsukubaensi*.



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-be@enzolifesciences.com)

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