

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 Ca²⁺-dependent phosphatase and protein phosphatase 2B (PP2B; calcineurin). Activates NF-κB through phosphorylation and degradation of IκBα.

Citations: 27

[View Online »](#)

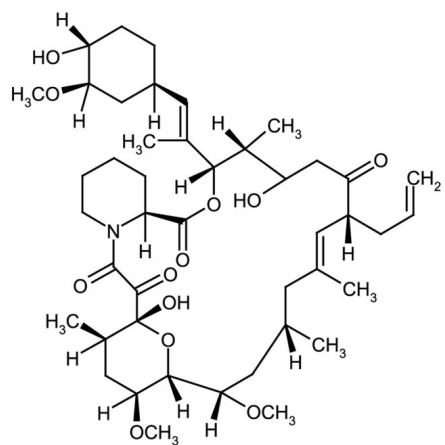
Ordering Information

[Order Online »](#)

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



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)

Isolated from *Streptomyces tsukubaensis*.

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)

Last modified: May 29, 2024