

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 Ca^{2+} -dependent phosphatase and protein phosphatase 2B (PP2B; calcineurin). Activates NF- κ B through phosphorylation and degradation of I κ B α .

Citations: 27

[View Online »](#)

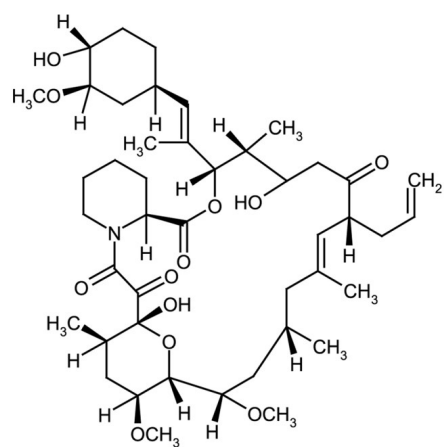
Ordering Information

[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



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-
fr@enzolifesciences.com](mailto:info-fr@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

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.